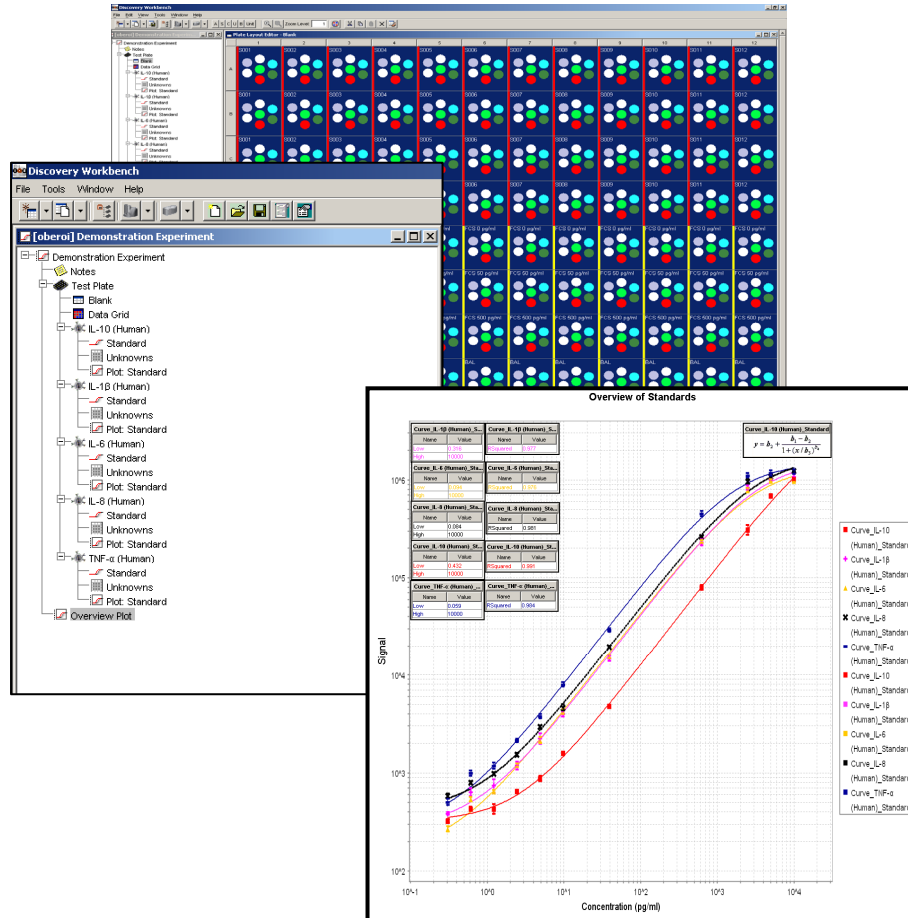


Data Analysis Toolbox

User's Guide



For Software Version 3.0



Meso Scale Discovery™
A division of Meso Scale Diagnostics, LLC.

Disclaimer of Warranty and Liability

MESO SCALE DIAGNOSTICS, LLC. (MSD) MAKES NO REPRESENTATION OR WARRANTY OF ANY KIND, AND HEREBY EXPRESSLY EXCLUDES AND DISCLAIMS ANY AND ALL REPRESENTATIONS AND WARRANTIES WITH REGARD TO THIS USER GUIDE (GUIDE), INCLUDING, BUT NOT LIMITED TO, ANY IMPLIED WARRANTY OF MERCHANTABILITY, SUITABILITY OR FITNESS FOR A PARTICULAR PURPOSE, DATA ACCURACY, SYSTEM INTEGRATION, QUIET TITLE OR NONINFRINGEMENT. NO STATEMENT IN THIS GUIDE SHALL BE INTERPRETED TO GRANT OR EXTEND ANY WARRANTY ON THE PRODUCTS DESCRIBED HEREIN. MSD SHALL NOT BE LIABLE FOR ERRORS OR OMISSIONS CONTAINED HEREIN OR FOR ANY DAMAGES OF ANY KIND, INCLUDING BUT NOT LIMITED TO DIRECT, INDIRECT, SPECIAL, INCIDENTAL, CONSEQUENTIAL OR PUNITIVE, WHETHER ARISING IN CONTRACT, TORT, STRICT LIABILITY OR OTHERWISE, THAT MAY BE INCURRED IN CONNECTION WITH THE FURNISHING, PERFORMANCE OR USE OF THIS GUIDE.

Changes in Publication

The information contained in this guide is subject to change without notice. We welcome comments on this publication and suggestions for future editions.

No Reproduction

No part of this guide may be duplicated, reproduced, stored in a retrieval system, translated, transcribed, or transmitted in any form or by any means without the express prior written permission of MSD. This guide shall be returned to MSD within two (2) business days following any request by MSD.

Meso Scale Discovery, Meso Scale Diagnostics, Discovery Workbench, Multi-Array, Multi-Spot, Sector, Sulfo-TAG, TAG, and MSD-Design are trademarks of Meso Scale Diagnostics, LLC. Unauthorized use of any Meso Scale Diagnostics, LLC. trademark, service mark or logo may be a violation of federal and state trademark laws.

All other trademarks are the property of their respective owners.

© 2006. Meso Scale Diagnostics, LLC. Gaithersburg, Maryland.

Web Site: <http://www.meso-scale.com/>

E-mail: sales@meso-scale.com

Table of Contents

1.	Introduction	1
1.1.	Multi-Array® Technology	1
1.2.	Intended Audience.....	1
1.3.	How to Use This Guide	1
2.	Data Analysis Toolbox Overview	2
3.	Definitions and Descriptions of Common Views	3
3.1.	Plate Data History.....	3
3.2.	Experiment View Window	5
3.3.	Loading, Saving, and Exporting	8
3.3.1.	Loading Plate Data into the Plate Data History.....	8
3.3.2.	Saving and Exporting Experiment.....	8
3.3.3.	Saving and Exporting Experiment Templates.....	9
3.3.4.	Saving and Exporting Plate Layouts	10
3.4.	Plate Layouts.....	11
3.4.1.	Step 1. Create New Layout.....	11
3.4.2.	Step 2. Assign Assays (Analytes).....	12
3.4.3.	Step 3. Format Plate Layout	13
3.4.4.	Step 4. Saving Plate Layouts	15
3.5.	Analyze Data and Modify Experiment	17
3.5.1.	Creating an Experiment	17
3.5.2.	Editing an Experiment.....	18
3.6.	Reports and Presenting Data	20
3.7.	Quick Reference Guide	22
4.	Practice Examples for Discovery Workbench Training	23
4.1.	Example 1: Quick Start Guide.....	23
4.1.1.	Step 1: Import the data file into the Plate Data History.	23
4.1.2.	Step 2: Import and Edit an Existing Plate Layout	24
4.1.3.	Step 3. Create an Experiment Using the Plate Layout	27
4.1.4.	Step 4. Generate a PDF report.	27
4.2.	Example 2: One Spot IL-17 Cytokine Data.....	28
4.2.1.	Step 1: Import the data file into the Plate Data History.	28
4.2.2.	Step 2: Create Plate Layout.....	29
4.2.3.	Step 3. Create an Experiment Using the Plate Layout	40
4.2.4.	Step 4. Manipulating and Configuring the Experiment.....	41

4.2.5.	Step 5. Saving, reporting, and copying plots and tables.....	48
4.2.6.	Advanced Functions: Excluding Data.....	52
4.2.7.	Advanced Functions: Detection Limits	52
4.3.	Example 3: 10-spot Mouse Cytokine Panel	55
4.3.1.	Step 1: Import the data file into the Plate Data History.	55
4.3.2.	Step 2: Create Plate Layout.....	56
4.3.3.	Step 3. Create an Experiment Using the Plate Layout	63
4.3.4.	Step 4. Manipulating and Configuring the Experiment.....	66
5.	Frequently Asked Questions	71
5.1.	Layouts.....	71
5.1.1.	How do I assign assays?	71
5.1.2.	How do I use an assay that is not in the list?	71
5.1.3.	How do I assign replicates?	71
5.1.4.	How do I enter unusual dilution patterns?.....	71
5.1.5.	How do I name my samples?	71
5.2.	Experiments	72
5.2.1.	How do I get my data into a new experiment?.....	72
5.2.2.	I have an experiment that I like, how do I reuse it?	72
5.2.3.	How do I share the templates/experiments with other users?	72
5.2.4.	How do I create an experiment template?	72
5.3.	Miscellaneous	73
5.3.1.	Where did the menu item go?	73
5.4.	Data Analysis.....	73
5.4.1.	How do I exclude data?.....	73
5.4.2.	How do I control detection limits?.....	74
5.4.3.	How do I control the fitting?	74
5.5.	Reporting and Working with Other Applications.....	74
5.5.1.	How do I print out the experiment?	74
5.5.2.	How do I remove items from the printed output?	74
5.5.3.	Can I get a text file containing the data?	74
5.5.4.	How do I get a plot/table into MS Word or MS Powerpoint? ..	75
6.	Appendix - Fitting functions.....	77
6.1.	Four Parameter Logistic Fit.....	77
7.	Technical Support.....	80

1. Introduction

1.1. Multi-Array® Technology

Meso Scale Discovery (MSD) develops, manufactures and markets solutions for complex biological assays, providing cost effective and valuable information to scientists in drug discovery, therapeutic screening, and basic life science research. MSD's solution-based product portfolio is based on Multi-Array technology, a proprietary combination of patterned arrays and electrochemiluminescence detection, enabling large numbers of measurements with exceptional sensitivity, wide dynamic range and convenience.

MSD markets detection instrument systems, a line of custom microplates for use with these instruments, and a proprietary line of reagents. MSD also provides services to customers. MSD® Multi-Array® plates are available in 96-, 384-, and 1536-well format, and in regular or high binding capacity formats. They may be purchased uncoated or coated with avidin, and custom surface treatments and coatings also are available.

MSD also markets its proprietary Multi-Spot® plates that have patterned arrays within each well of a microplate. Multiple analytes are measured simultaneously within a single well, increasing throughput and enabling novel assay formats. Custom surface treatments, array preparations, and coatings are also available for these plates.

For more details and information about applications, please visit the MSD web site at www.meso-scale.com.

1.2. Intended Audience

This guide is for all users of the Sector Instruments and describes how to analyze data from Multi-Array plates run on these instruments. Users should understand general computer/ Microsoft® Windows® terminology, and be familiar with standard laboratory practices. The intended users of the Sector instruments are those conducting research in the Life Sciences.

The Sector Imager 2400, Sector Imager 6000 and Sector PR share many features. Throughout this manual, the reader should assume that a feature is common to the analysis of data derived from Multi-Array plates run on any of these instruments unless otherwise noted. The user should refer to the appropriate manual for guidance on running Sector instruments and acquiring data from them.

1.3. How to Use This Guide

This guide is divided into chapters containing main topics and subsections. Use the Table of Contents and the Index to find topics of interest quickly. The Figures list and Tables list indicate the locations of all images and tables that enhance understanding of written information within this guide. The Appendix contains supplemental information regarding shortcut keys, instrument specifications, safety symbols and labels, and robotics integration. A Glossary is also included.

2. Data Analysis Toolbox Overview

The MSD Discovery Workbench 3.0 Data Analysis Toolbox includes simple straightforward method for analyzing data generated from MULTI-ARRAY plates run on the MSD SECTOR Instruments. Data can be analyzed as it is collected or after it has been stored in the Plate Data History. The data analysis functions allow simple analysis of complex data sets. Experiments are created and stored in the Plate Data History database along with the plate data. The software uses 4-parameter logistic curve fit to fit standards and calculates concentrations for unknowns and controls. Tables, graphs, notes, and reports can be quickly generated and published and shared through PDFs or presented in Powerpoint, Word, or Excel documents. User tracking and change control allows experiments to be published in an unalterable format and assigned to an owner for traceability. The Data Analysis Toolbox can be installed at the user's desktop in addition to the instrument for convenient data analysis. Plate layouts, experiment templates, experiments, and plate data can be shared between users.

The DISCOVERY WORKBENCH 3.0 has a simple workflow for data analysis:

1. Run Plate
2. Create Plate Layout
3. Analyze Plate Data
4. Modify tables and plots
5. Print report

Alternatively, once plate layouts and experiment templates have been created, the template can be selected when running the plate to automatically create an experiment.

This manual describes the features of the software and has two examples that walk the user through analysis of a single spot assay and a multiplex plate.

For more information on the MULTI-ARRAY technology visit our website at www.meso-scale.com.

3. Definitions and Descriptions of Common Views

3.1. Plate Data History

Plate Data History: All of the data (plates and experiments) are stored in the Workbench's Plate Data History. Experiments and plates can be retrieved through the Plate Data History. New to the Discovery Workbench is the Experiment View tab in the Plate Data History, which tracks all of the experiments.

Library – The library stores plate layouts and experiment templates that can be used to analyze data.

Plate Data File – A specific text file generated by Discovery Workbench. The file can be generated by opening a plate in the Plate Data History and exporting to a text file. A text file is automatically created when a plate is read.

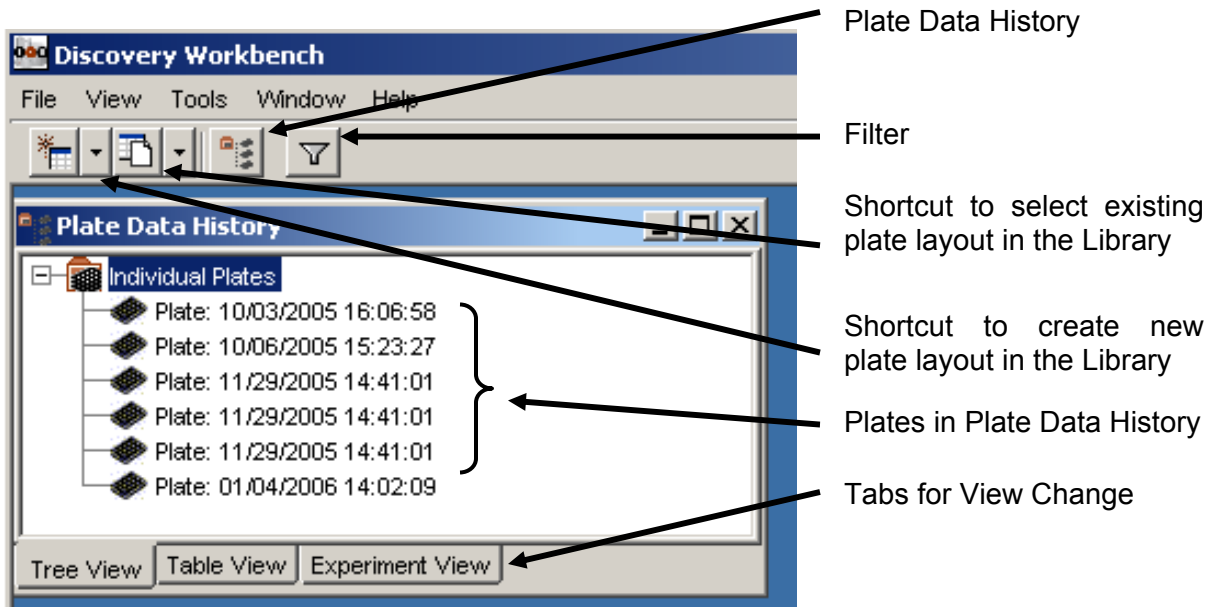


Figure 1. Plate Data History: Tree View.

Filter – Opens a new filter dialog to filter what is viewed in the Plate Data History. Filtering can be done by owner, plate type, date created, plate name, or barcode.

Individual Plates – Plates that are run as single plates are stored under the *Individual Plates* section. If plates are run as stacks, the plates will be stored under the *Runs* section.

Experiment - The analysis containing a layout, ECL results from one or more plates, tables and plots generated from a plate data file. The analyzed data and reports are stored in an experiment.

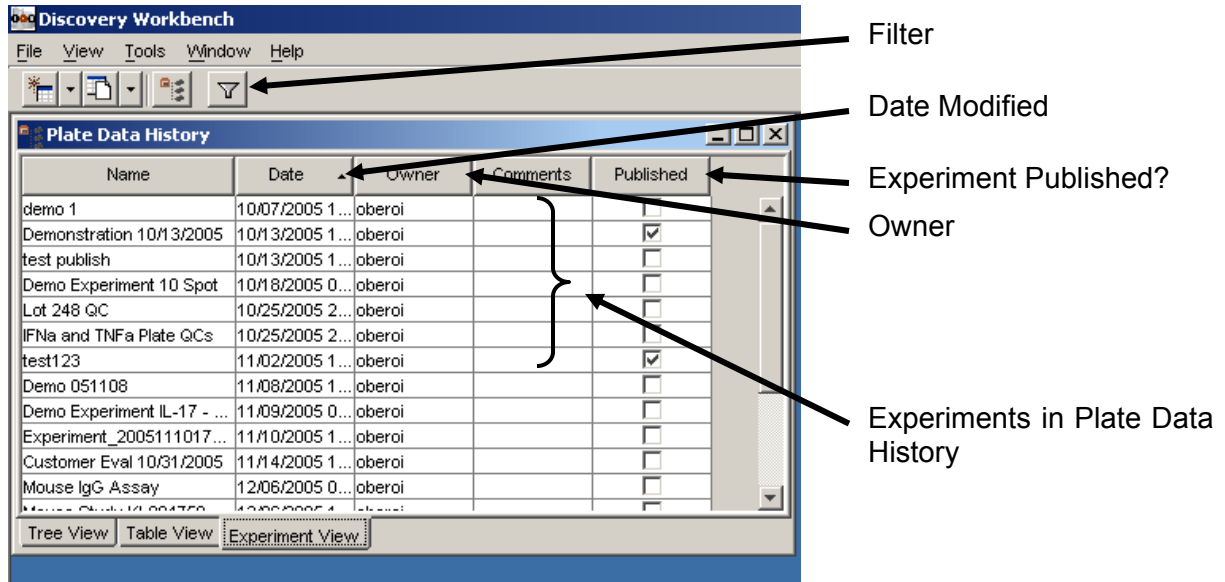


Figure 2. Plate Data History: Experiment View.

Experiment Name – The experiment name is defined by the user when an experiment is created.

Experiment Date – The date the experiment was last modified.

Experiment Owner – The operator who made the last modification to the experiment.

Experiment Comments – Area for comments.

Published – Once an experiment has been published, it can no longer be modified. The owner and the experiment date are locked in the database.

3.2. Experiment View Window

The experiment window is automatically named by the owner and the experiment name. The experiment is a collection of tables and plots that are viewed in a tree view. The tree can be expanded or collapsed at different branch points.

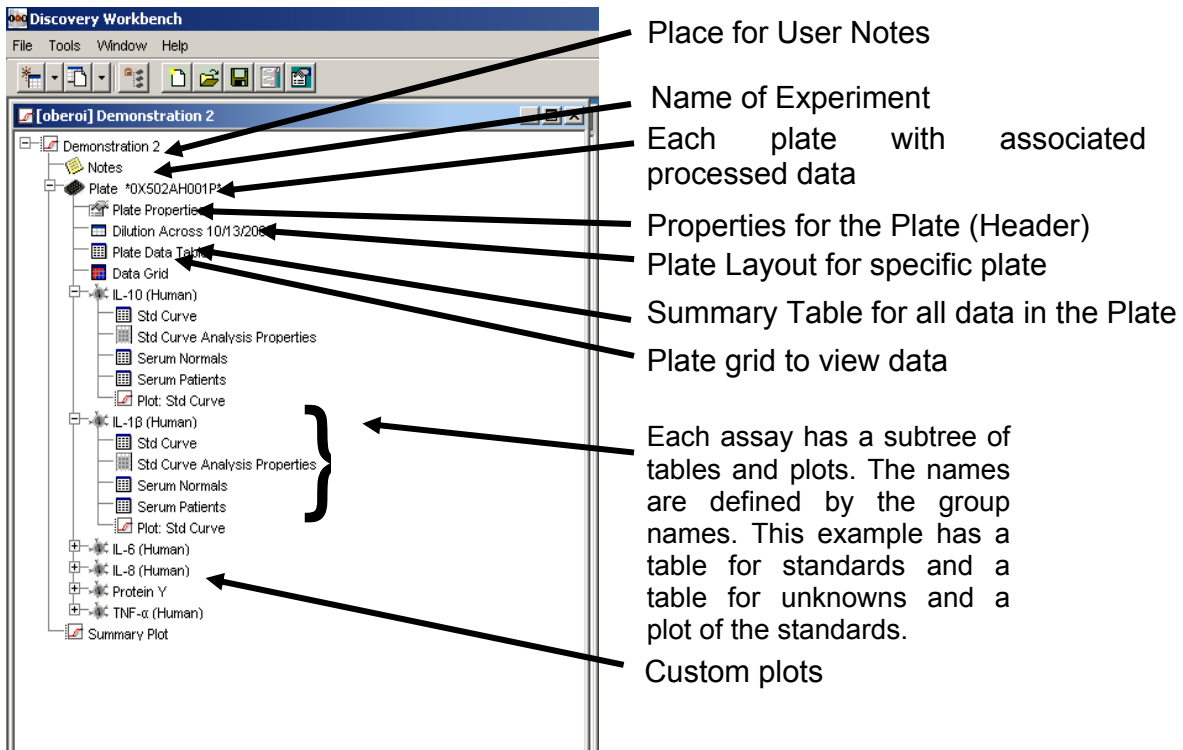


Figure 3. Overview of Experiment Tree View

Plate Data – Each plate forms a branch of a tree that can be collapsed. Under each plate there are plate properties, the plate layout, Plate Data Table, Data Grid, and analyzed data for each assay within the plate.

Plate Properties – Contains the information from the plate header.

Plate Layout – In this example the plate layout is called *Dilution Across 10/13/2005*. The plate layout is stored locally in the experiment and can be modified without modification of the plate layout in the Library.

Plate Data Table – A table of all of the results from the plate in one location. This table can be exported to Excel and used with Pivot tables to quickly make summary tables.

Data Grid – Plate layout view of the data, sample, and well properties.

Custom Plots – Custom plots can be created and associated with the experiment.

Plate Layout (Editor) – The mapping of the types of assays and types of samples on the plate. The assay part of the layout describes which assays are on which spots and in which well. The sample part of the layout describes the standards, controls and unknowns with corresponding concentrations and dilutions.

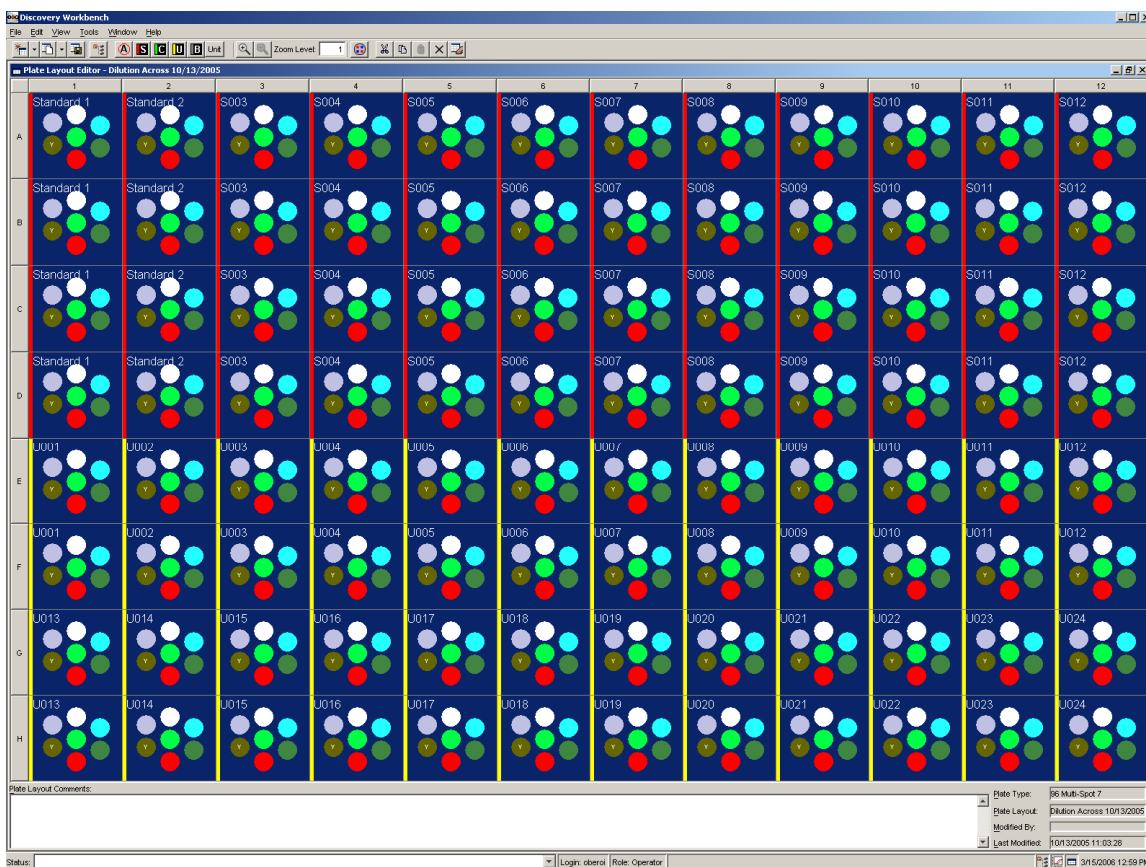


Figure 4. Plate Layout Editor View and the Plate Layout Toolbar

A: Assign Assay – Assignment for the assays in the well. For a MULTI-SPOT plate, there will be multiple assay assignments per well.

S: Assign Standards – Assignment for wells having a known concentration of analyte/target that is used to generate a calibration (standard) curve.

C: Assign Controls – Assignment for wells having a known concentration of analyte/target that is used to assess accuracy of the standard curve.

U: Assign Unknowns – Assignment for wells containing a sample that is being measured.

B: Assign Blanks – Assignment for wells containing a background control that can be subtracted or divided from other wells.

Unit – Assignment of the unit for the samples and standards. The default unit is pg/ml.

Groups – For each sample type, groups of wells can be defined. The groups allow for multiple standard curves and multiple control or unknown populations. A separate analysis will be done for each group, including a separate table and plot (if it is a standard).

Assay Subtree – For each assay there is an auto-generated set of tables and plots. For each assay, a table will be created for each group of samples (standards, controls, unknowns, or blanks). For each set of standards, a table of the curve fitting analysis properties is generated and a plot for each standard curve.

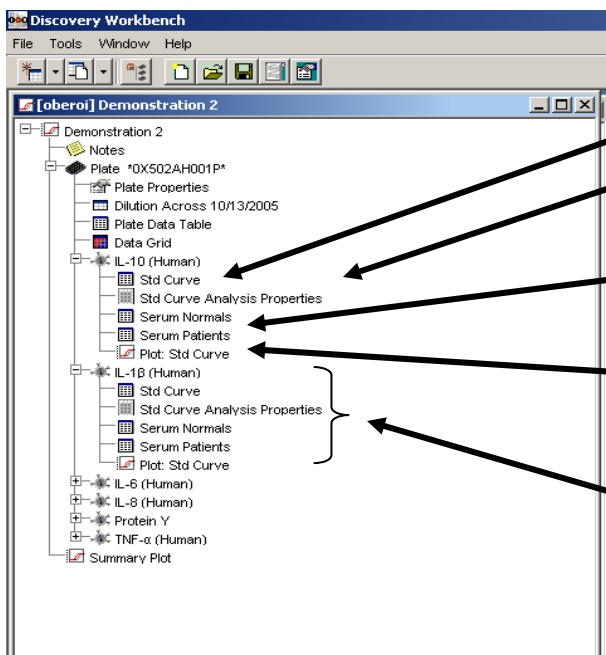


Table for standards
(called Std Curve)

Analysis properties for the curve fit

Two table for unknowns (two
groups Serum Normals and
Serum Patients)

Plot for standards

Each assay has a subtree of
tables and plots. The names
are defined by the group
names. This example has a
table for standards and a table
for unknowns and a plot of the
standards.

Plate Data Table – Table formatted as a spreadsheet containing all of the data for a plate in an Experiment. Each table can be exported to a text file or to a clipboard to be imported into other applications such as Microsoft Excel.

Plot – The plot is a line graph of the Standards and the curve fit generated for that Standard group.

Experiment Template – A blank experiment that is stored without data. The template contains all the formatting of the tables, graphs, and plate layouts. When running a plate, an experiment template can be used to generate an experiment automatically.

Fit Profiles – The algorithm and initial parameters used to fit the standard curve.

3.3. Loading, Saving, and Exporting

3.3.1. Loading Plate Data into the Plate Data History

The **Plate Data History** cannot be shared between computers, so when using the Discovery Workbench on a computer not connected to an instrument, the plate data file must be loaded into the **Plate Data History**. Alternatively, experiments, templates, plate layouts, and profiles can be shared between computers using the import and export feature for experiment.

To import plate data, the plate data must be exported in the correct format from the instrument. To create the correct format, from the instrument's **Plate Data History**, export the plate data using the default Export Format. This file contains all of the information needed to transfer the plate data and plate properties to another Discovery Workbench. The file will be named with the MSD Barcode and a time stamp of the time the plate was run (for example: oW502AI274._2006-02-06-133756.txt)

To load the plate data into the **Plate Data History**, open the **Plate Data History** window. When the window is selected, select the File menu from the toolbar and select *Import->Plate Data File*. If there is an error in the import, try to re-export the data from the instrument. Usually a custom export format is used to save data generated during a plate run but in this case the standard format should be chosen to insure all of the data is transferred completely.

Once the plate has been loaded, the plate data view will open up and the plate should be inserted into the Plate Data History.

3.3.2. Saving and Exporting Experiment

Experiments are saved in the **Plate Data History**. When an experiment is closed, the experiment can be saved or a copy can be saved using the *Save As* function from the File Menu.

If an experiment needs to be shared between Discovery Workbench installations, an experiment can be exported from *File->Export Experiment*. This option is available only when an experiment window is open and selected. When an experiment is exported, it is saved as a .zip file. This zip file stores all of the experimental data including the graphs, tables, plate layout and experiment. The exported experiment can be loaded into the **Plate Data History** by selecting *Import->Experiment* from the File menu.

One can export just the data from an experiment can be exported to a text file using the **Export Data to Text File** function in an experiment. Right-click on the Experiment name or the plate tree and select *Export Data to Text File*. If the function is selected from the Experiment, all the plates in the experiment will be exported. If it is selected from a specific plate, only that plate will be selected to be exported.

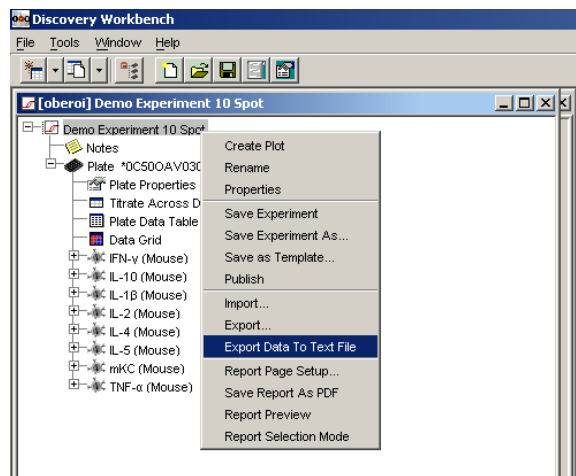


Figure 5. View of the Export Data to Text File drop-down menu from the Experiment window.

The **Export to Text File** dialog box will open up. This is similar to the **Export Data** function from the instrument or Plate Data History. The data can be exported in the default format, or a custom format to the Output Path. Different than the Export Data function, the data type that is exported can be selected. If the text file is to contain the calculated concentrations for each of the wells, select Calc Concentration. If the text file is to contain the ECL signal, select ECL Signals from the Data Type dropdown menu.

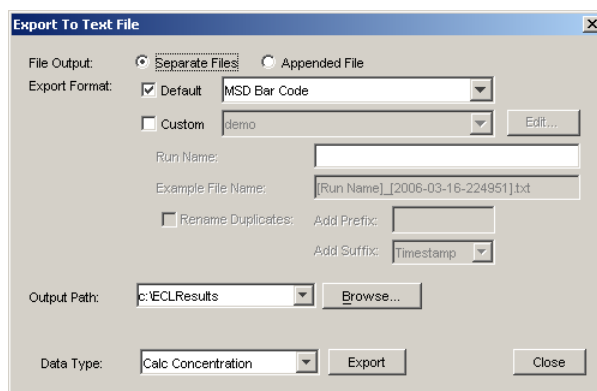


Figure 6. Export to Text File Dialog Box.

If the experiment has been published, the data cannot be exported to a text file.

3.3.3. Saving and Exporting Experiment Templates

Experiment templates can be saved by right clicking on the experiment in the experiment window (see Figure 5). The experiment template is stored in the Library in the database.

To export a template close the experiment once it is saved as a template, you can open up the template from *File->Open Experiment Template*. Once the template is open, the template can be exported from the File menu.

Templates can be imported into the Library using the *File->Import->Template* function when the Plate Data History window is selected.

3.3.4. Saving and Exporting Plate Layouts

Plate layouts can be saved in multiple locations. A library of Plate Layouts is stored in the Discovery Workbench database. The plate layouts are stored by plate type. Once a plate layout is used in an experiment, a copy of that plate layout is saved with the experiment. Modifications to the plate layout in the experiment do not change the plate layout in the Library unless the **Save Plate Layout in Library As** function is selected from the *File* menu. Plate Layouts can be shared between Discovery Workbench installation through the export and import functions.

3.4. Plate Layouts

3.4.1. Step 1. Create New Layout

Plate layouts are specific to an instrument type and a plate type. Plate layouts are stored in the Library. New plate layouts can be created by selecting *File->New Plate Layout in Library* or by clicking the New Plate Layout icon on the Discovery Workbench toolbar.



Figure 7. Discovery Workbench Toolbar: New Plate Layout in Library Icon

Once a new plate layout has been selected, the New Plate Layout dialog box will open. Select the plate type and provide a name for the plate layout. After clicking New, the Plate Layout Editor will open up with a Plate View. An example of a 7-spot plate view is shown in Figure 8.

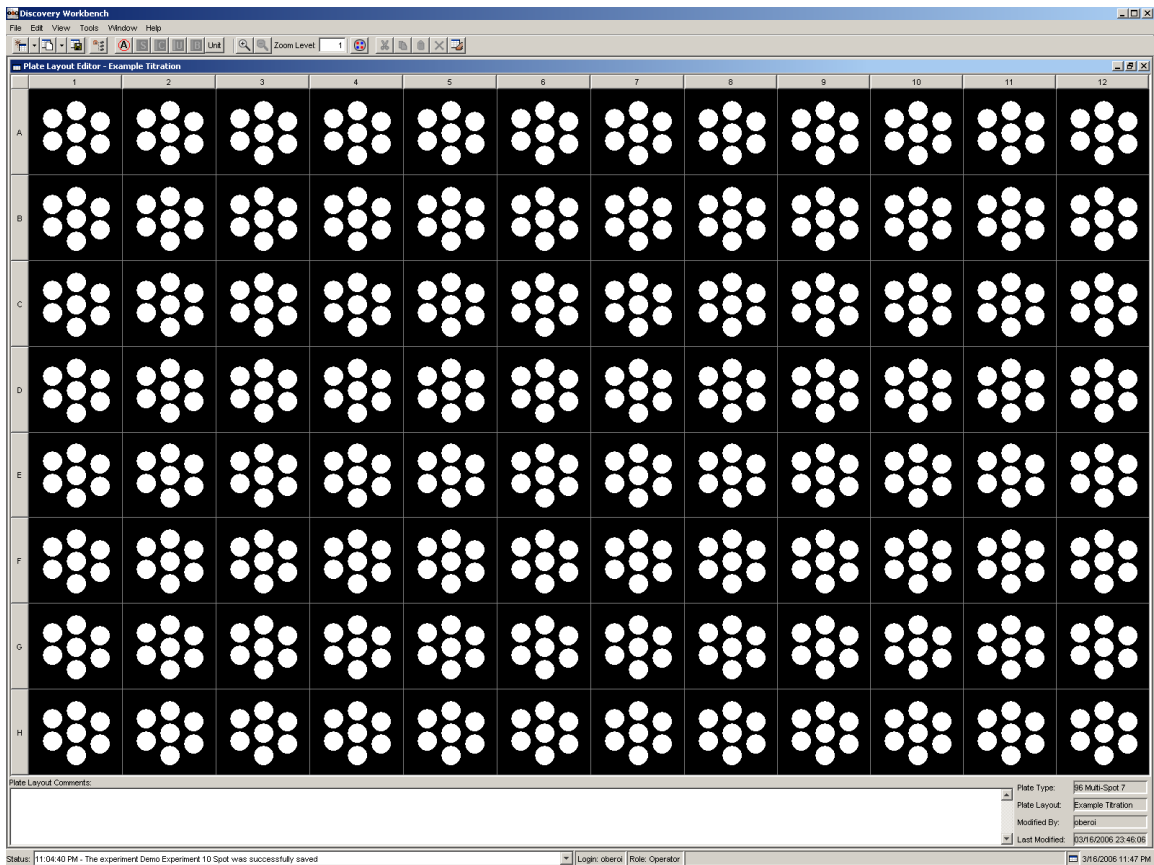


Figure 8. Plate Layout Editor View

3.4.2. Step 2. Assign Assays (Analytes)

An assay must be assigned to each spot that is to be analyzed (assays are not assigned to BSA blank spots). Selecting the **A** icon in the toolbar, opens up the Assign Assay dialog from which assays can be assigned to each spot. Alternatively, the assays can be assigned by selecting *Edit->Assign Assays*.

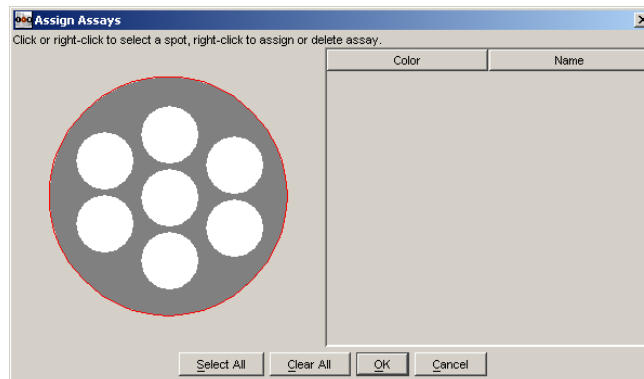


Figure 9. Assign Assay Dialog Box

Right clicking on a spot opens up a menu for selecting an assay. Discovery Workbench is preloaded with many of the assays available through MSD. New assay names can be selected through the **New** function in the pull down menu (Figure 10).

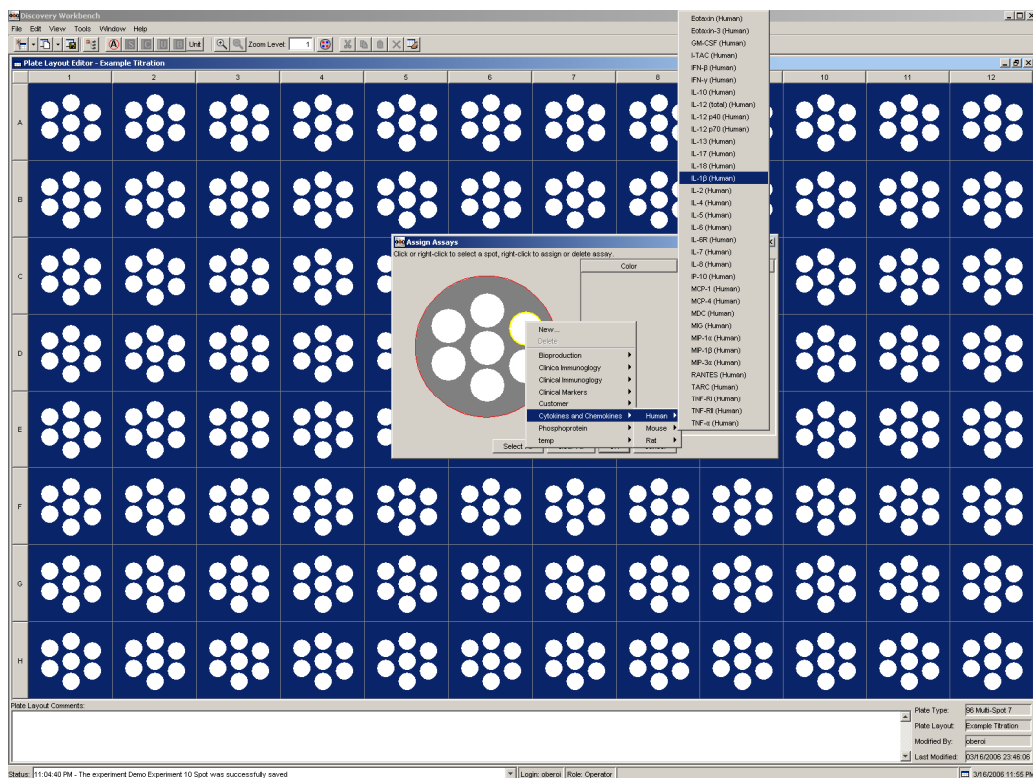


Figure 10. Assign Assay pull down menu.

Once the assays have been assigned, the spots will be color-coded for the assays.

3.4.3. Step 3. Format Plate Layout

Assigning Standards, Controls, Unknowns, and Blanks (sample types) can be done through either the toolbar or the *Edit* menu as shown below.

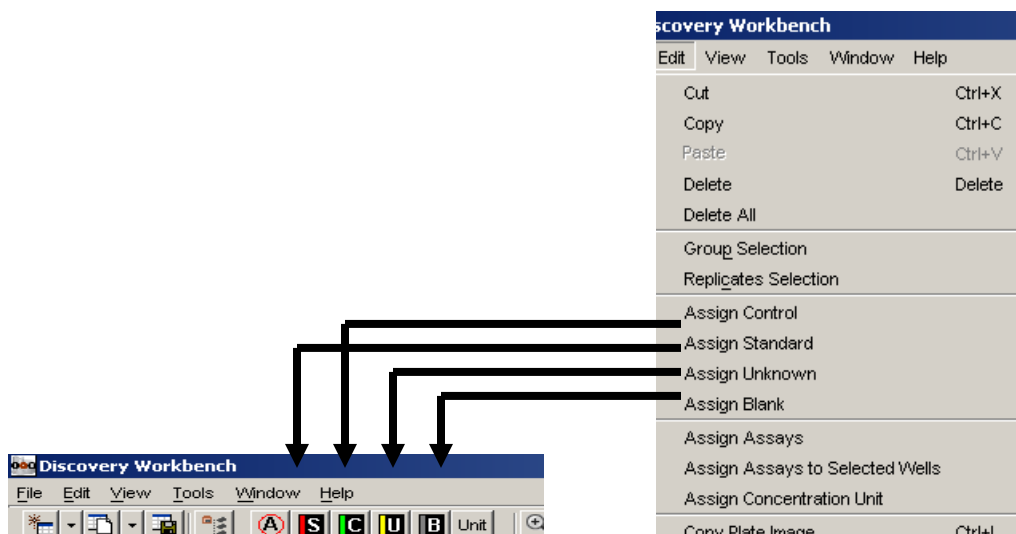


Figure 11. Assign Standards, controls, unknowns, blanks, units, and assays toolbar.

To assign wells to a particular sample type, highlight the wells in the plate view then assign the appropriate sample type. It is best to start assigning standards, follow with controls and finish with unknowns. That way you can take advantage of the autofill options to set which standard applies to which wells.

- 1) Select a set of wells. The set does not have to be all together. You can press and hold the *control button* and *click on wells*, or *press and hold the shift* and *click on wells* to generate any pattern of selected wells.
- 2) Right click on the layout and pick assign standard, assign control, or assign unknown as desired. Alternatively, the icon in the toolbar can be selected or the function from *Edit->Assign*.
- 3) A dialog will appear which is used to define the samples. The upper part of the dialog is used to make assignments and the lower table shows the results and allows direct editing of values.
- 4) Pick a group name. For many users this can simply be any name you wish to use. The name appears in the output and can also be used to identify samples that have the same name but were processed differently, etc. Items with the same sample name and the same assay but in different groups are analyzed separately.
- 5) Use the replicate choices as needed. NOTE: replicates can also be created by using copy and paste on the layout. This may be more efficient than using the replicates buttons in some cases.
- 6) If you have concentration/dilutions that follow simple factors, the autofill options can be used to load much of the concentration/dilution information.
- 7) The table can be directly edited as desired. NOTE: changes in replication options or use of the auto fill options will overwrite this table, even if it contains hand entered sample names or concentrations.

8) Click OK and repeat as needed for the other sample types or groups

An example of setting up a plate layout is shown in the practice examples.

A key component of assigning standards, controls, or unknowns is the group they are associated with. Default group names will be assigned, but one can form multiple groups and have multiple standard curves.

When assigning standards and controls, each assay can have a different concentration. This can be accomplished by selecting/deselecting the **Apply To All Assays** checkbox and selecting which assay to apply the concentrations to. The default setting is that all the assays have the same concentrations.

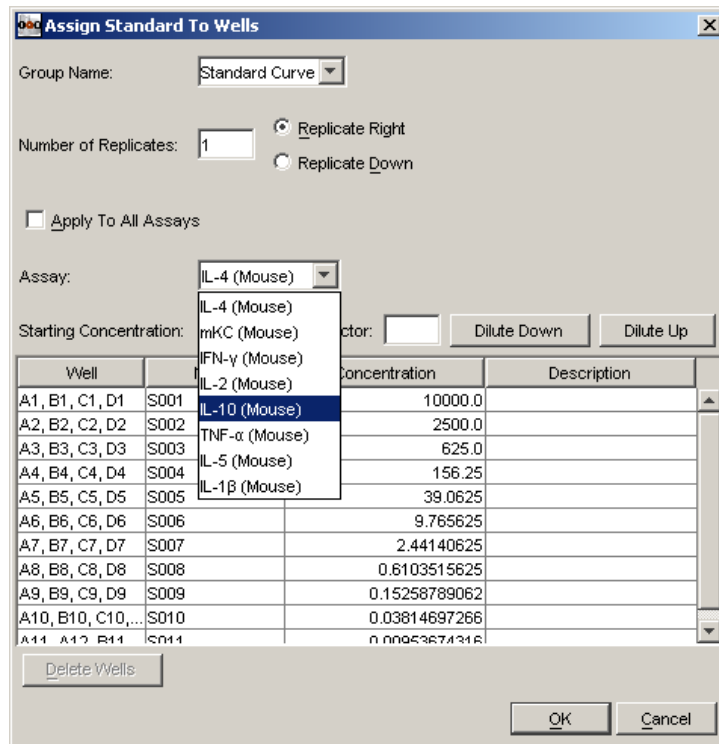


Figure 12. Selecting a concentrations for a specific assay when assigning concentrations to standards.

3.4.4. Step 4. Saving Plate Layouts

New or edited Plate Layouts should be saved in the library before proceeding to analyze data. To save a Plate Layout either close the Plate Layout window and you will be prompted to save the Plate Layout or select the *File->Save Plate Layout in Library* or use the shortcut button on the toolbar.

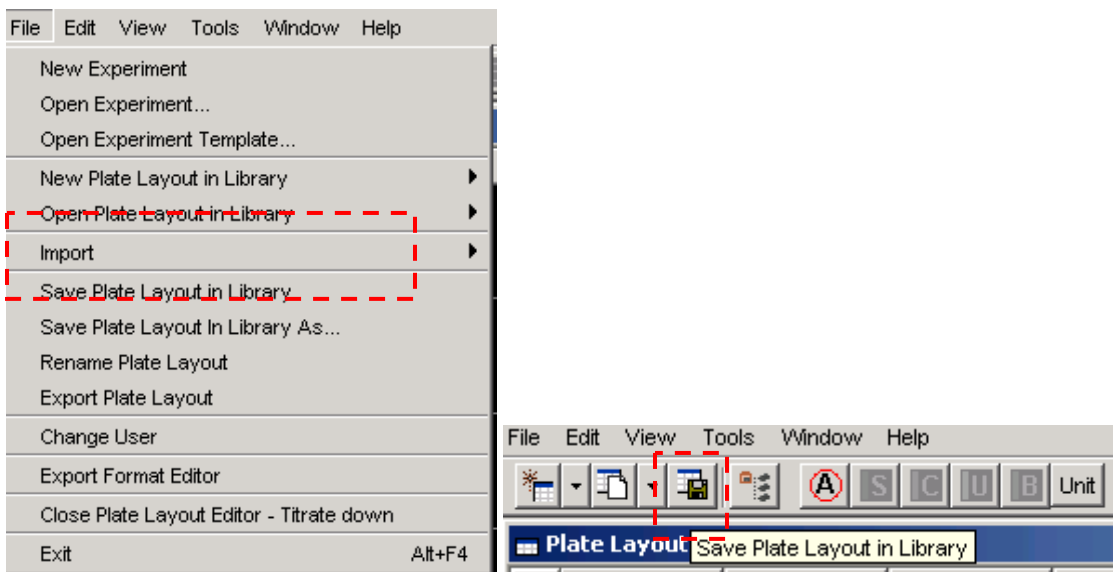


Figure 13. Saving a Plate Layout using either the pulldown from the File menu or the toolbar at the top of the window.

3.5. Analyze Data and Modify Experiment

A plate can be quickly analyzed once a Plate Layout has been created or edited to reflect the distribution of samples and assays on the plate.

3.5.1. Creating an Experiment

To create an Experiment, highlight the plate to be analyzed in the Plate Data History. Select *File->Analyze Plates*. You may also right click on a highlighted plate, and then select Analyze Plates. A pull down menu with the available Plates Layouts for that plate type will be shown. Select the appropriate Plate Layout and click OK.

An experiment window will open automatically.

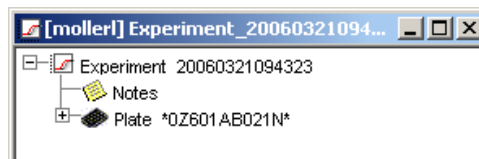


Figure 14. The Experiment window.

The tree is expanded by clicking on the +.

Experiments, plates and groups can be renamed by right clicking on the item and selecting **Rename** from the menu.

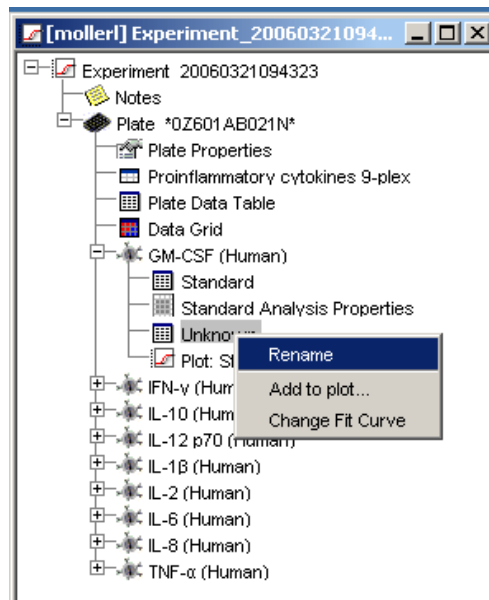


Figure 15. Renaming an Experiment, Plate or Group.

3.5.2. Editing an Experiment

Results are presented in Tables and Plots within an Experiment. If required, Tables and Plots may be customized further.

Editing a Table

To edit a Table, first open the Table by double-clicking it in the Tree view.

Sorting

Clicking on any of the column headers will sort the data in the Table from low to high or high to low according to that header.

Sample	Well	Concentration (pg/ml)	Signal	Mean	CV	Calc. Concentration (pg/ml)	Calc. Conc. Mean (pg/ml)	Calc. Conc. CV
S001	A02	2,500	969,359	973,806.5	0.646	2481	2493	0.682
	A01		978,254			2505		
S002	B02	625	257,707	257,932	0.123	630	631	0.126
	B01		258,157			631		
S003	C01	156.25	65,796	65,864.5	0.147	157	157	0.149
	C02		65,933			158		
S004	D02	39.062	17,049	16,991.5	0.479	39.8	39.7	0.49
	D01		16,934			39.6		
S005	E01	9.766	4,152	4,152.5	0.017	9.15	9.15	0.018
	E02		4,153			9.15		
S006	F02	2.441	1,307	1,338.5	3.33	2.46	2.53	4.11
	F01		1,370			2.61		
S007	G02	0.61	584	525	15.9	0.774	0.638	30.3
	G01		466			0.501		
S008	H01	0	242	244.5	1.45	0	0	N/A
	H02		247			0		

Sample	Well	Concentration (pg/ml)	Signal	Mean	CV	Calc. Concentration (pg/ml)	Calc. Conc. Mean (pg/ml)	Calc. Conc. CV
S008	H01	0	242	244.5	1.45	0	0	N/A
	H02		247			0		
S007	G02	0.61	584	525	15.9	0.774	0.638	30.3
	G01		466			0.501		
S006	F02	2.441	1,307	1,338.5	3.33	2.46	2.53	4.11
	F01		1,370			2.61		
S005	E01	9.766	4,152	4,152.5	0.017	9.15	9.15	0.018
	E02		4,153			9.15		
S004	D02	39.062	17,049	16,991.5	0.479	39.8	39.7	0.49
	D01		16,934			39.6		
S003	C01	156.25	65,796	65,864.5	0.147	157	157	0.149
	C02		65,933			158		
S002	B02	625	257,707	257,932	0.123	630	631	0.126
	B01		258,157			631		
S001	A02	2,500	969,359	973,806.5	0.646	2481	2493	0.682
	A01		978,254			2505		

Figure 16. Sorting a Table. In this example, the data in the Standards Table are sorted from high to low (top) and low to high (bottom) concentration

Adding and Removing Columns

To add or remove columns, right click anywhere on the column titles then select or deselect columns from the menu. Columns can be moved around by dragging them with the left mouse button.

Excluding Data

To exclude data from a Table, highlight the data to be excluded, right click and choose **Exclude Selected**. To reverse this process, highlight the data to be re-included, right click and choose **Include Selected**.

Editing a Plot

To edit a Plot, either right click the Plot in the Tree View and select **Edit Plot Properties** from the menu or open the Plot by double-clicking it in the Tree view, then go to **Edit > Edit Plot**. In this window, the Plot name can be modified, error bars and detection ranges can be turned on or off and the shape and color of the data points for the curve can be edited.

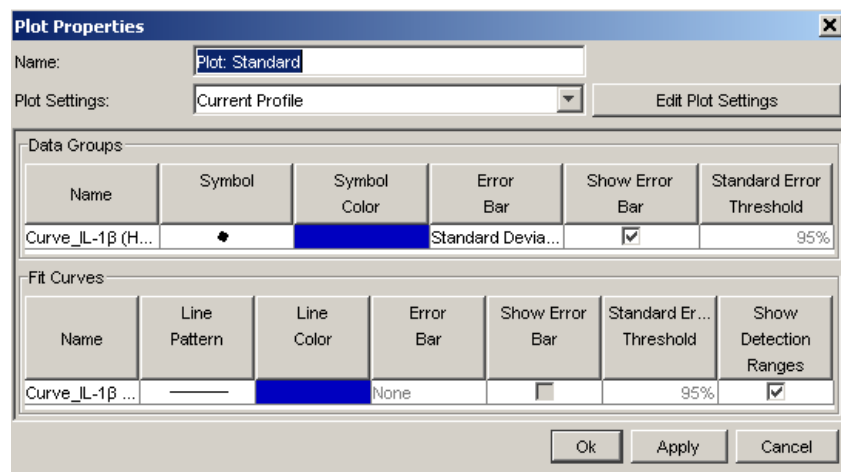


Figure 17. Editing Plot Properties: Plot name, line and symbol color, symbol shape, error bars, displaying detection range.

To edit axes and Plot Legend, click on the Edit Plot Settings button. This window may also be accessed when a plot is open from the **Edit > Plot Settings** menu or by right clicking the Plot in the Tree View and selecting **Edit Plot Settings** from the menu.

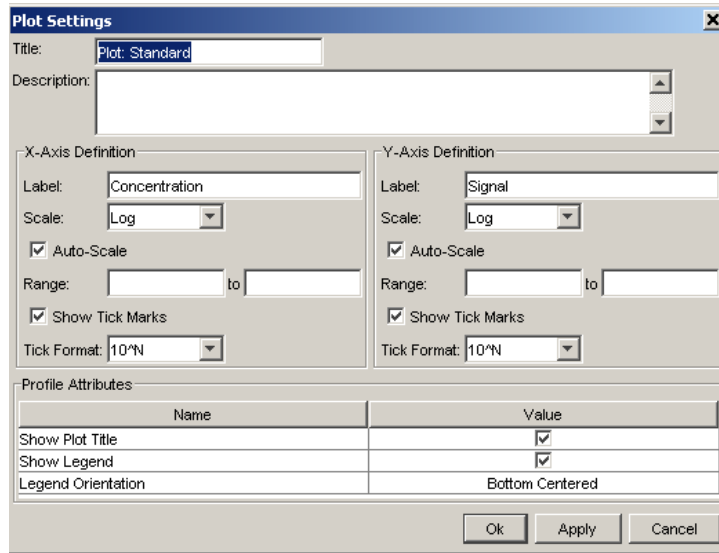


Figure 18. Editing Plot Settings: Plot title, Axes, Legend orientation

3.6. Reports and Presenting Data

3.6.1. Creating a Report

To compile all data from an experiment in a PDF report, highlight the experiment name, right click and select **Save Report As PDF** from the menu. The content of a PDF reports can be customized by first selecting sections to be included using the **Report Selection Mode** in the experiment right click menu.

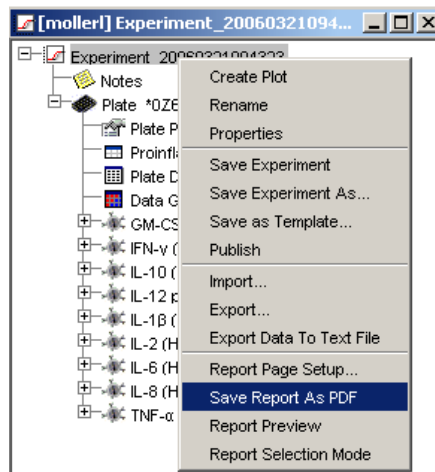


Figure 19. Creating a PDF report.

3.6.2. Presenting Data

The data from an experiment can be transferred to other software programs.

To copy a Table, open the Table by double clicking it in the Tree menu. In the **Edit** menu, select **Copy Table to Clipboard** to copy the data for further manipulation or **Copy Image to Clipboard** to copy an image of the Table.

To copy a Plot, open the Plot by double clicking it in the Tree menu. In the **Edit** menu, select **Copy Plot** to copy an image of the Plot. Plots or tables copied to the clipboard can then be pasted into another application.

Shortcut buttons are also available to copy the Table to the clipboard (Figure 19).

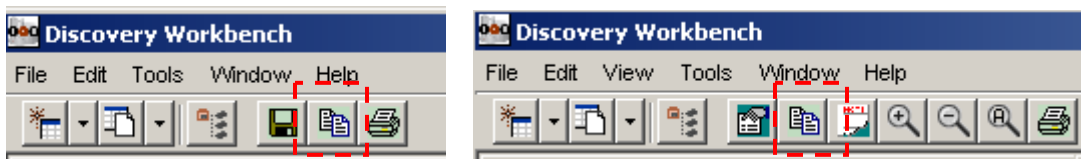


Figure 20. Shortcuts for copying a Table (left) or Plot (right) to the clipboard.

3.7. Quick Reference Guide



Create new plate layout. Use the drop down menu to select the appropriate instrument.



Edit existing plate layout. Use the drop down menu to select a plate layout from the libraries.



Plate data history. Contains data of previously run plates and experiments.



Assign assay



Assign standard



Assign control



Assign unknown



Assign Blank



Assign concentration units



Save plate layout



Spot Map

4. Practice Examples for Discovery Workbench Training

4.1. Example 1: Quick Start Guide

This example highlights the basic steps for analyzing MSD data using an existing Plate Layout.

4.1.1. Step 1: Import the data file into the Plate Data History.

Open MSD Discovery Workbench. Open the Plate Data History by selecting the icon, or selecting it from the **Tools** menu at the top. Look for a plate that was created on **11/08/2005 11:15:50**.



If the data is *not* in the **Plate Data History** of the Workbench DAT, then you must import the plate data file in the default export format. The default file name consists of the plate barcode and a time stamp for the plate run.

With the **Plate Data History** open, select **File->Import->Plate Data File** then select the file *Example1.txt*. The text file was exported from the run using the default MSD barcode export format. This plate will now be imported into the Plate Data History.

A new plate should open up and the data window should open and a plate created on *11/08/2005 11:15:50* should be created in the Plate Data History. The plate type is 96 and it was run on a Sector Imager 6000.

4.1.2. Step 2: Import and Edit an Existing Plate Layout

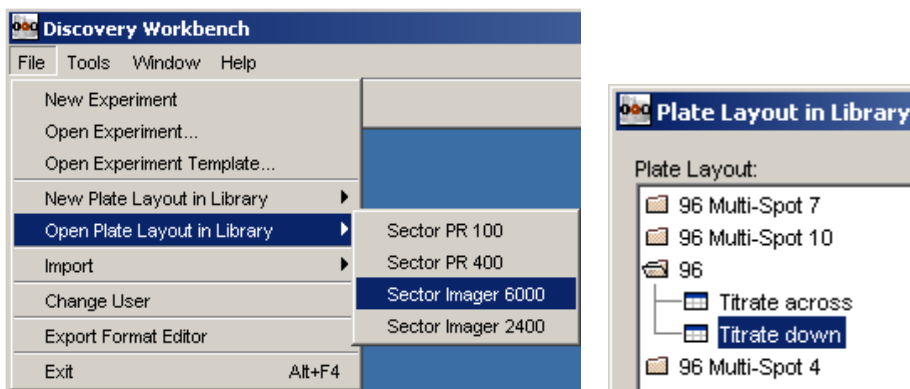
The next step is to import and edit an existing plate layout to reflect the contents of the plate. Plate layouts and experiment templates are stored in a library in Discovery Workbench.

Step 2.1 Import Layout.

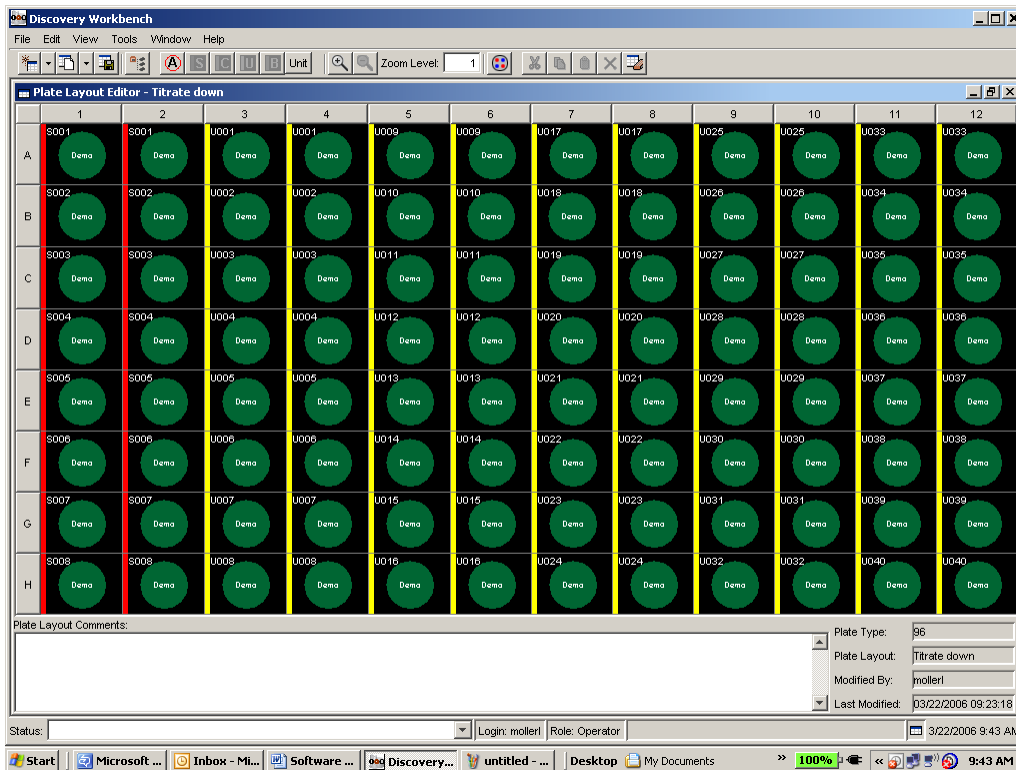
- From the **File** menu, select **File->Import ->Plate Layout**. Select *Titrate Down Layout.zip*

Step 2.2: Select Layout

- From the **File** menu, select **File->Open Plate Layout in Library->Sector Imager 6000**. Select *96* and the plate layout named *Titrate Down* from the **Plate Layout** library menu and click **Open**.



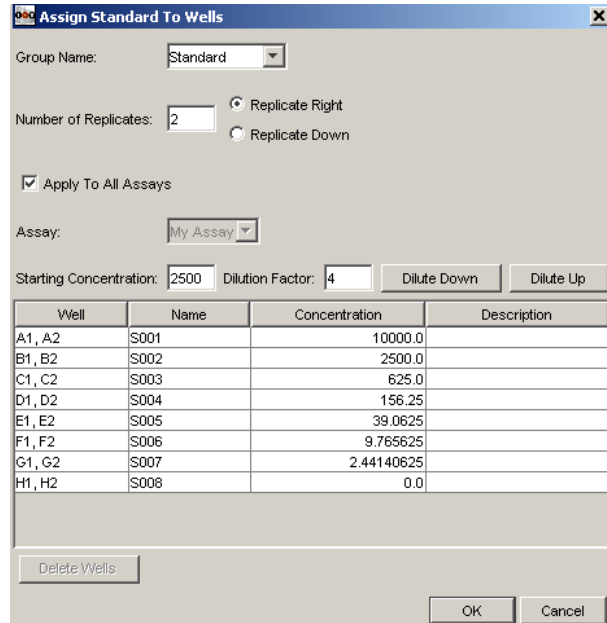
A plate layout should appear as shown below:



Step 2.3: Edit Plate Layout

In this plate layout, the standards for the calibration curve are in duplicate in columns 1 and 2 and samples (unknown) are in duplicate in columns 3-12. The top calibrator is set at 10,000 pg/ml and the other calibrators are serial 1:4 dilutions. The calibrators in the plate data file are different from those in this plate layout – they are 4 fold dilutions starting at 2500 pg/ml with the lowest calibrator set to 0 pg/ml.

- To edit the standards, highlight columns 1-2 and click on the **S** button on the toolbar. This will open the Standard window shown below.



- To change the concentrations of the standards, enter **2** in the Number of Replicates field, **2500** in the Starting Concentration field and **4** in the Dilution Factor field and click dilute down. In the concentration column, double click on the H1, H2 row and change the value to 0 then click OK.

Step 2.4 Save the Plate Layout

- Save the Plate Layout by going to **File->Save Plate Layout in Library** and then close the window by selecting the **x** in the upper right part of the plate view window.

4.1.3. Step 3. Create an Experiment Using the Plate Layout

Now that you have modified the plate layout in the Library, the plate can be analyzed.

- Highlight the plate to be analyzed in the Plate Data History (For this exercise the plate is: *11/08/2005 11:15:50*). Select **File->Analyze Plates**. You may also right click on the highlighted plate, then select Analyze Plate.
- A pull down menu with the available Plates Layouts for that plate type will be shown.

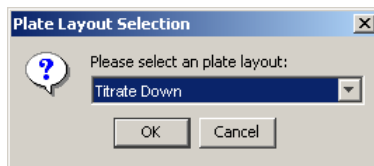


Figure 21. Pulldown window for selecting the Plate Layout to be used for analyzing an experiment.

- Select *Titrare Down* as the plate layout. Click OK, and an Experiment is automatically created as shown below. Expand the tree by clicking on the +.
- To open a table or plot within the tree, double click on the item and a new window displaying that item will open.

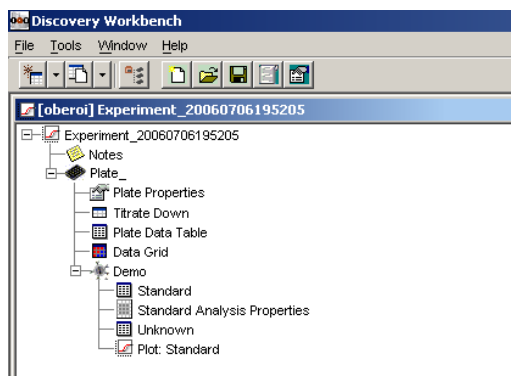


Figure 22. Experiment tree created in Example 1.

4.1.4. Step 4. Generate a PDF report.

A PDF report of the Experiment can be saved for printing or data sharing purposes.

For this experiment, select **File -> Save Report as PDF** and save the PDF as ***Experiment1report.PDF***. The report should be 17 pages long.

Close the experiment. You may choose to save or not save the experiment for this exercise.

4.2. Example 2: One Spot IL-17 Cytokine Data

This example demonstrates procedures involved in the use of MSD Workbench Data Analysis Tools using example data from a single spot plate.

4.2.1. Step 1: Import the data file into the Plate Data History.

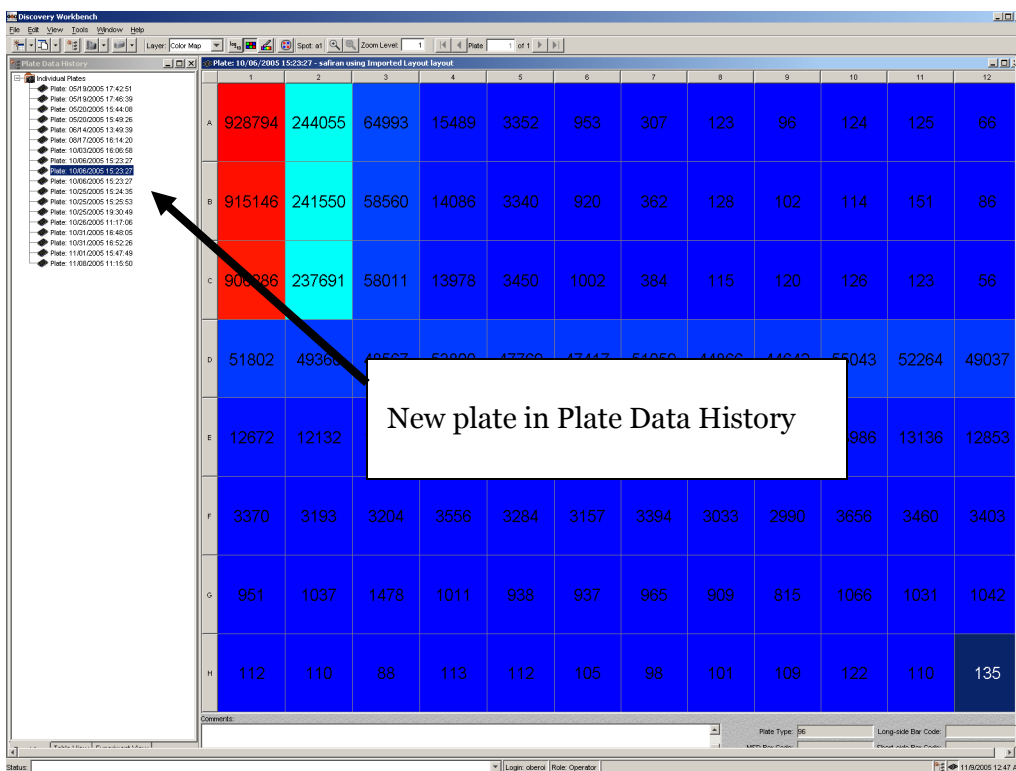
If the data is not in the **Plate Data History** of the Workbench DAT, then you must import the plate data file in the default export format. The default file name consists of the plate barcode and a time stamp for the plate run.

- Open MSD Discovery Workbench. Open the Plate Data History by selecting the icon, or selecting it from the **Tools** menu at the top.



Once the Plate Data History is open, select **File->Import->Plate Data File** then select the file *Example2.txt*. The text file was exported from the run using the default MSD barcode export format. This plate will now be imported into the Plate Data History.

A new plate should open up and the data window should open and a plate created on **10/06/2005 15:23:27** should be created in the Plate Data History. The plate type is 96 and it was run on a Sector Imager 6000.



The data is a titration of human cytokine IL-17 standards in rows A-C with unknown samples below. The standard curve consists of a 1:4 dilution series of calibrator from 10,000 pg/ml to 0.153 pg/ml across the top 3 rows in triplicate. The bottom 5 rows contain controls. For the first part of the example, we will treat these as unknowns assayed in triplicate across.

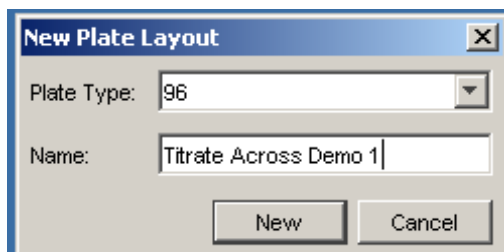
4.2.2. Step 2: Create Plate Layout

As a good practice, close the plate window before continuing. Having too many (greater than 20) windows open in Discovery Workbench could reduce processing speed.

The next step is to generate the plate layout to analyze the plate. Plate layouts and experiment templates are stored in a library in Discovery Workbench.

Step 2.1: Select Plate Type

- From the **File** menu, select **File->New Plate Layout in Library->Sector Imager 6000**. Select 96 from the **Plate Type** dropdown menu. For the name, type in: *Titrate Across Demo 1*. Click on the New Button.



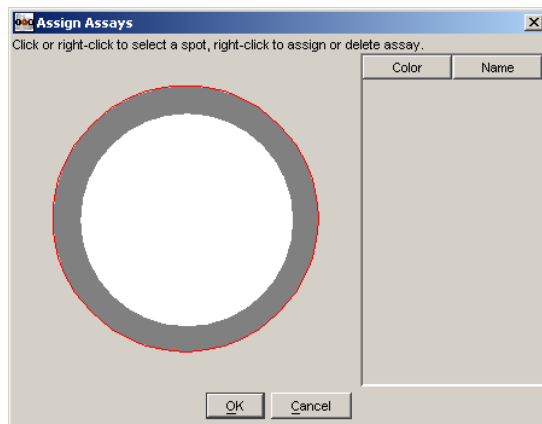
This should open a 1-spot plate view with the spot being white. The Toolbar should appear on the top of the window as shown below:



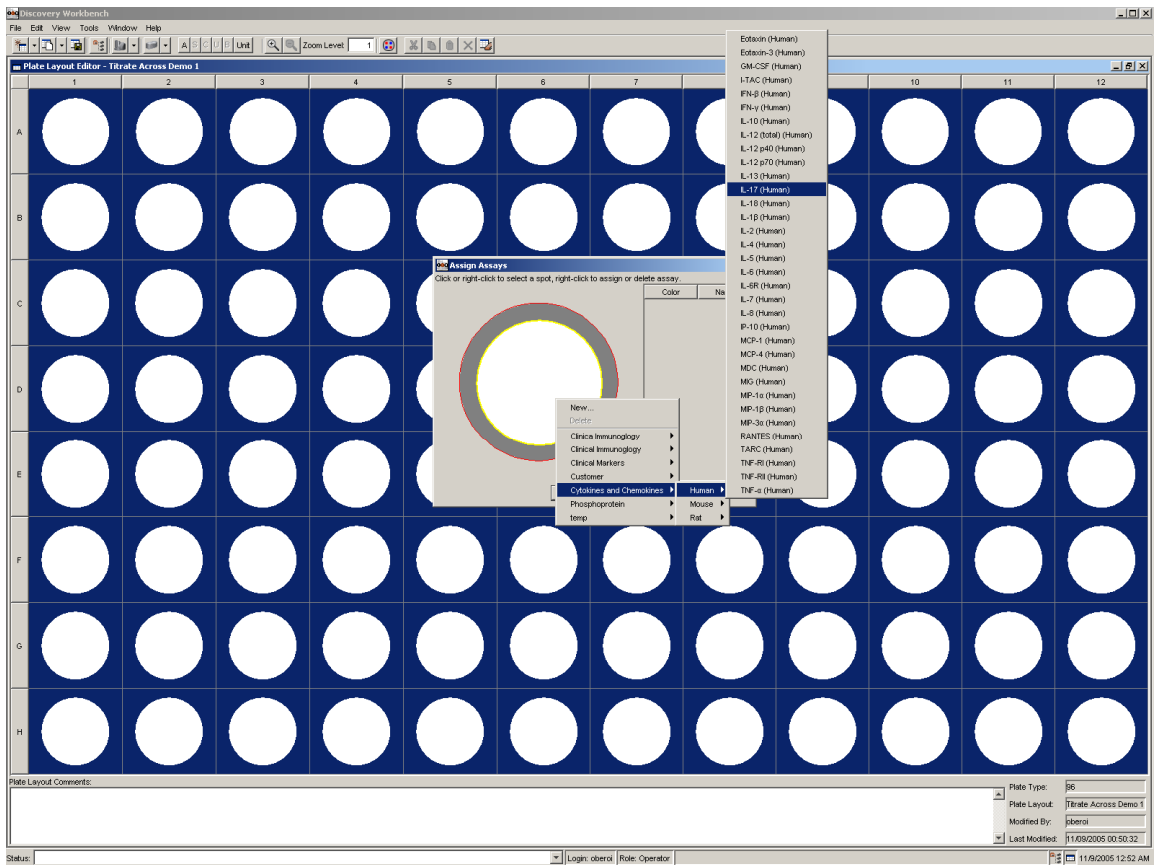
Step 2.2 Assign Assays (A)

The **A** button in the toolbar is used to assign the assays for the spots.

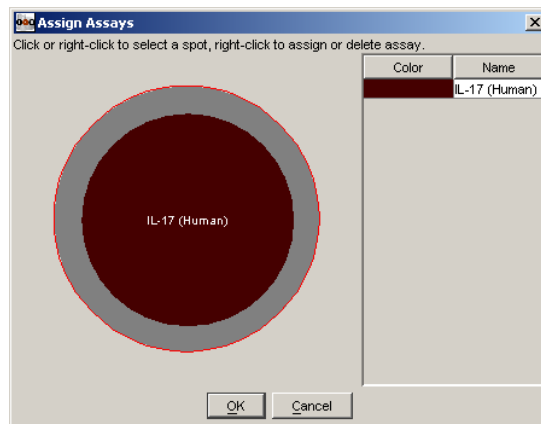
- Left click on the **A**, and an **Assign Assays** window will open as shown below:



- To assign an assay to the spot, right click on the spot and a drop down menu will open up. From the drop down menu, select **Cytokines and Chemokines->Human->IL-17 (Human)**.



After you have assigned the assay, the **Assign Assays** window should look like the one below.



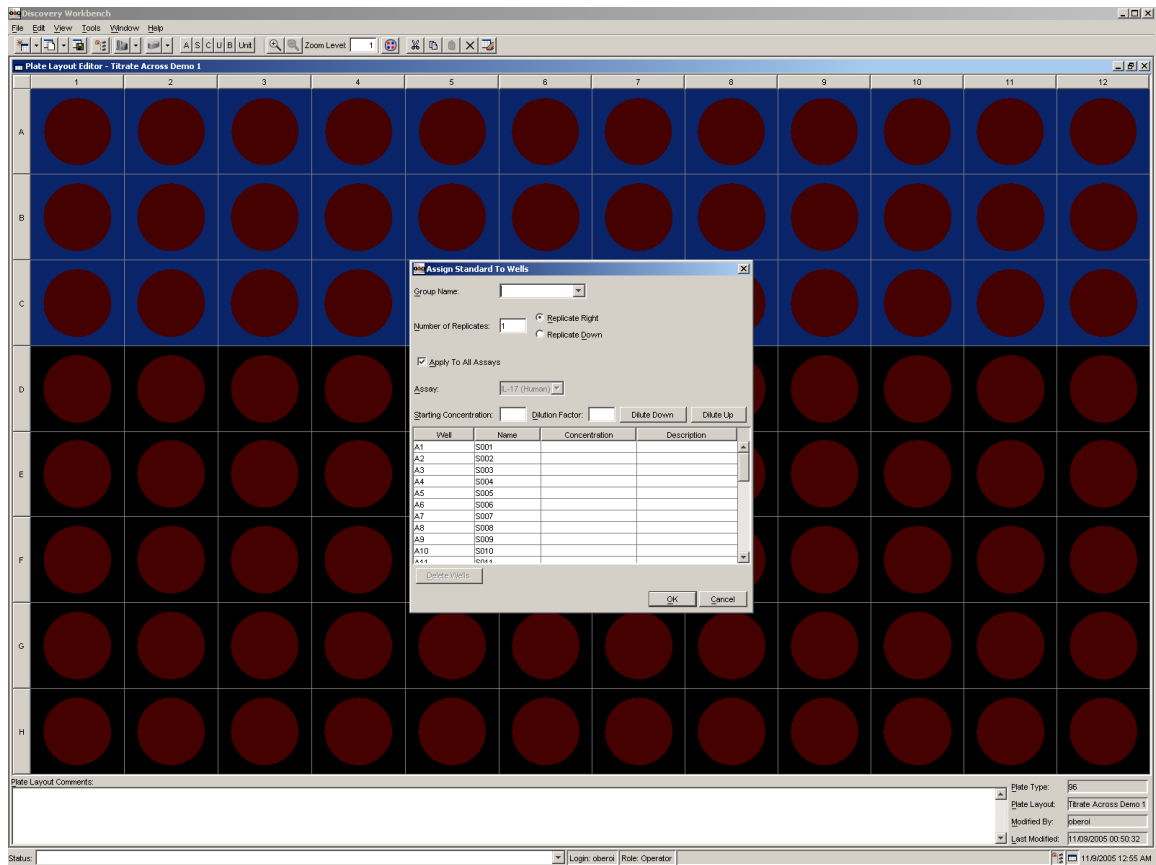
After assigning the assay, click OK and you're ready to move on to assigning standards, controls, unknowns and blanks. Notice that other buttons on the toolbar **S**, **C**, **U**, **B**, and **Unit** are now enabled.

In the case of a multispot plate, follow the procedure outlined above for each spot. The legend and colors represent which assays are assigned to which spots.

Step 2.3 Assign Standards (S)

The plate map is selectable like a table in Excel. You can select individual wells, rows, columns, or use the control and shift keys to select interesting pattern.

- For this example, select rows A, B, C, and D. Then click on the **S** button in the tool bar to open the **Assign Standards to Wells** window. A representative screen shot is shown below:



The **Assign Standard to Wells** window should look like the following:

Well	Name	Concentration	Description
A1	S001		
A2	S002		
A3	S003		
A4	S004		
A5	S005		
A6	S006		
A7	S007		
A8	S008		
A9	S009		
A10	S010		
A11	S011		

- In the *Group Name* type in: *Std Curve*. Since this example has 3 replicates, change the number of replicates to 3 and select *Replicate Down*. The wells should now be grouped A1, B1, C1 and A2, B2, C2... Since the standard curve starts at 10,000 pg/ml and we diluted across by a factor of 4, enter the starting concentration at 10,000 and dilution factor of 4, then click on the Dilute Down button. Scroll down to S012 and overwrite the concentration to be 0. Note that **Name**, **Concentration**, and **Description** can be modified in the table as well. The final window should look like the following:

Assign Standard To Wells

Group Name: Standard Curve

Number of Replicates: 4 Replicate Right Replicate Down

Apply To All Assays

Assay: mKC (Mouse)

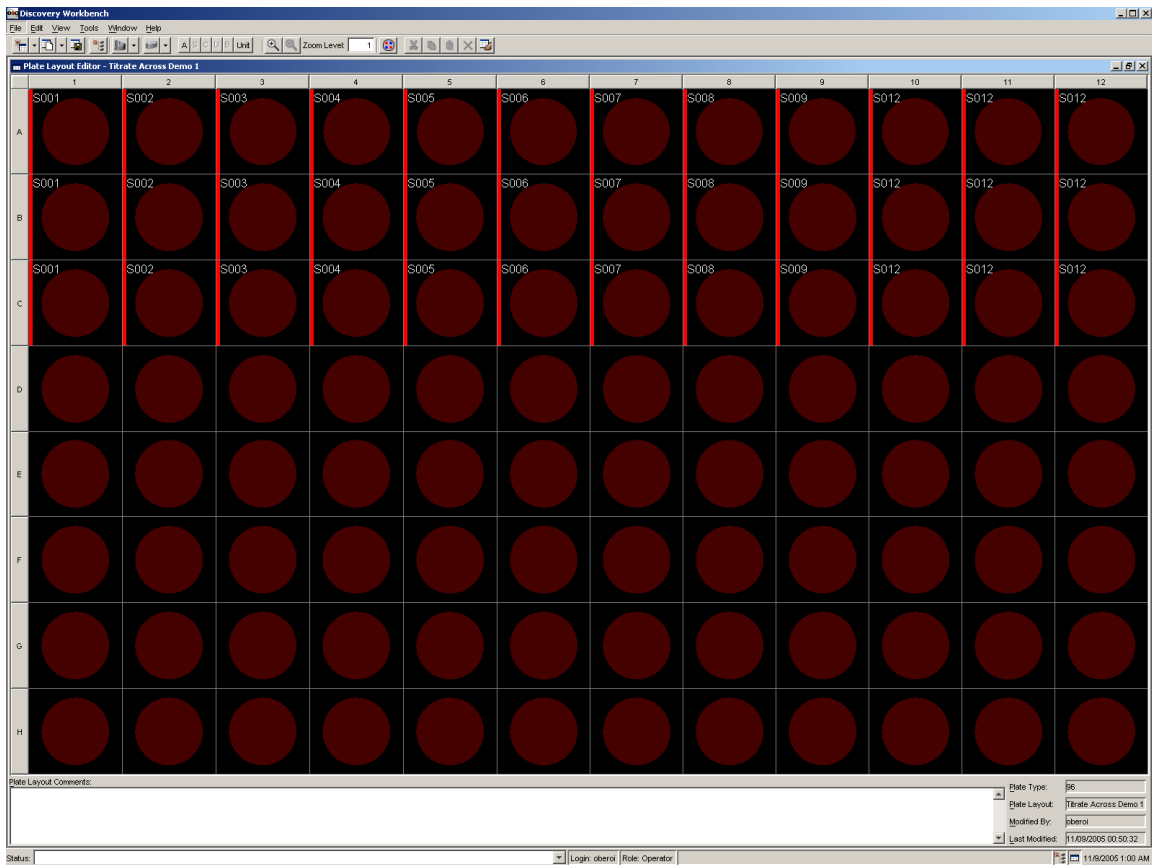
Starting Concentration: 10000 Dilution Factor: 4

Well	Name	Concentration	Description
A1, B1, C1, D1	S001	10000.0	
A2, B2, C2, D2	S002	2500.0	
A3, B3, C3, D3	S003	625.0	
A4, B4, C4, D4	S004	156.25	
A5, B5, C5, D5	S005	39.0625	
A6, B6, C6, D6	S006	9.765625	
A7, B7, C7, D7	S007	2.44140625	
A8, B8, C8, D8	S008	0.6103515625	
A9, B9, C9, D9	S009	0.15258789062	
A10, B10, C10,...	S010	0.03814697266	
A11, B11, C11	S011	0.00953674316	

- Click OK and the Plate layout should now have a color description with the sample names in the well.

We deliberately made a mistake in this example. The problem was that the standards in A10, B10, C10, A11, B11, and C11 were also zero concentration. How can we correct this?

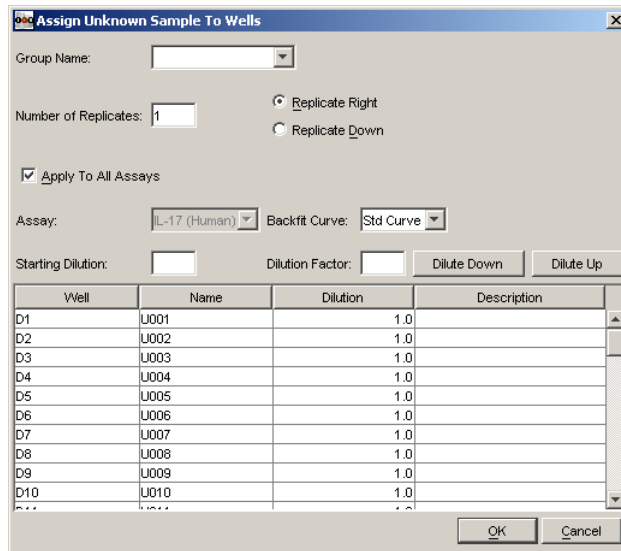
- First select wells A10, B10, C10, A11, B11, C11 in the plate view and press the **DELETE** button on your keyboard. Now select A12, B12, and C12 as a group. Press **Ctrl-C** to copy those wells. The 3 wells can be copied as a group. Select A10 and press **Ctrl-V** to paste S12 into those wells and repeat for A11. The grid should appear as shown below with 9 replicates of S12.



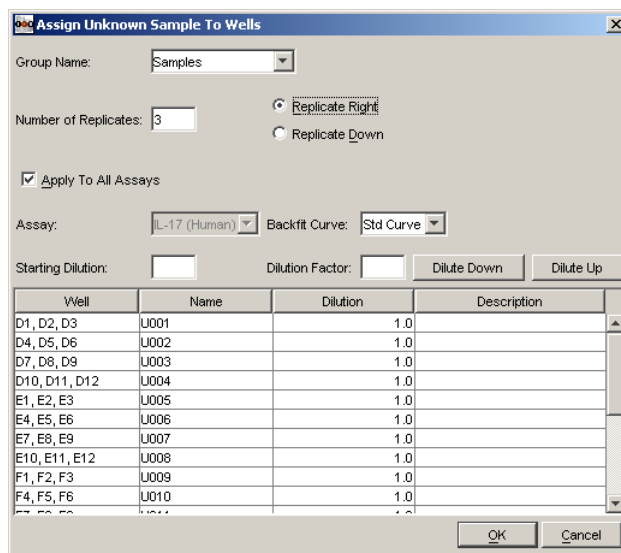
Step 2.4 Assign Unknowns (**U**)

The bottom 5 rows contain unknowns that are assayed in triplicate.

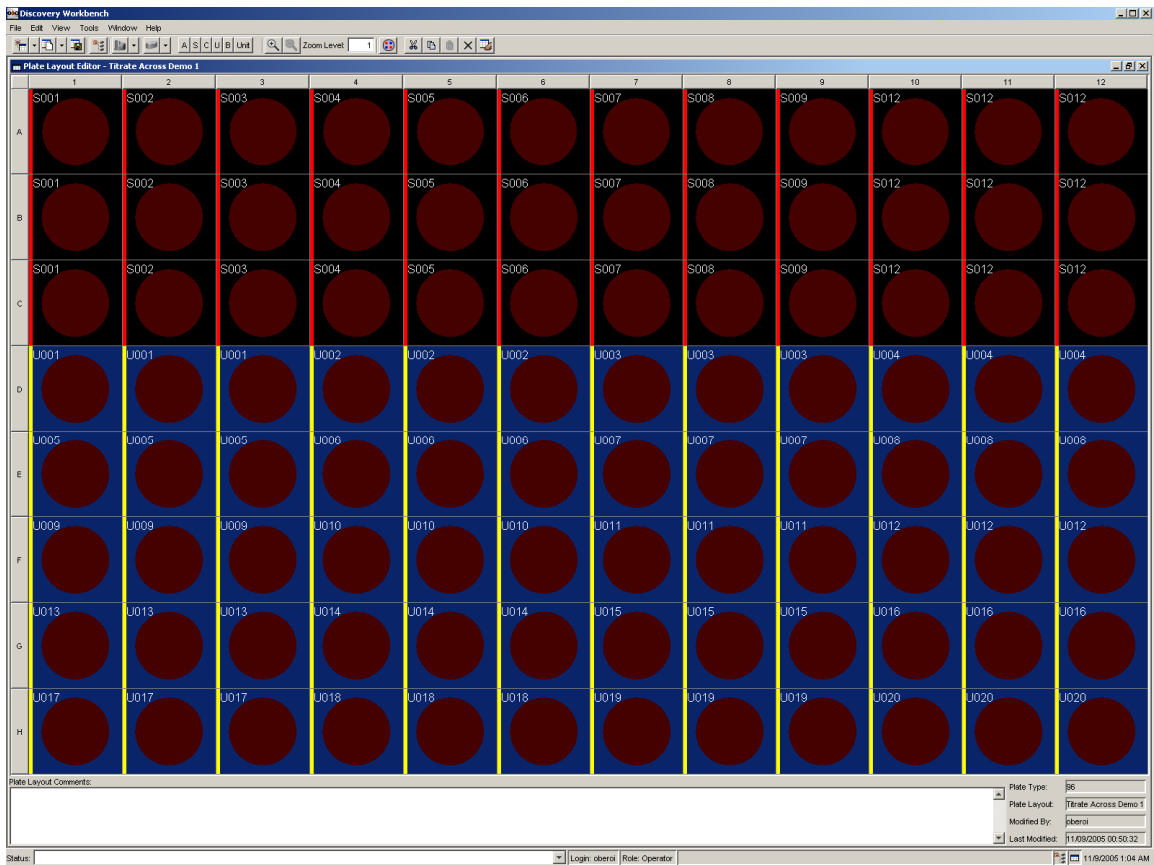
- Select rows D, E, F, G, and H and then click on the **U** button in the tool bar to open the Assign Unknowns to Wells window. This step is similar to assigning the standards, with a few exceptions. First, one can select which backfit curve to fit the unknowns data to if there are multiple standard curves on the plate. Second, one can assign the dilution rather than the concentration. The calculated concentrations will take into account the dilution factor.



- Create a name for this group called Samples. Since we have triplicates going across rather than down, enter 3 for the number of replicates and then select Replicate Right. All the dilutions should remain at 1. The window should now appear as shown below:



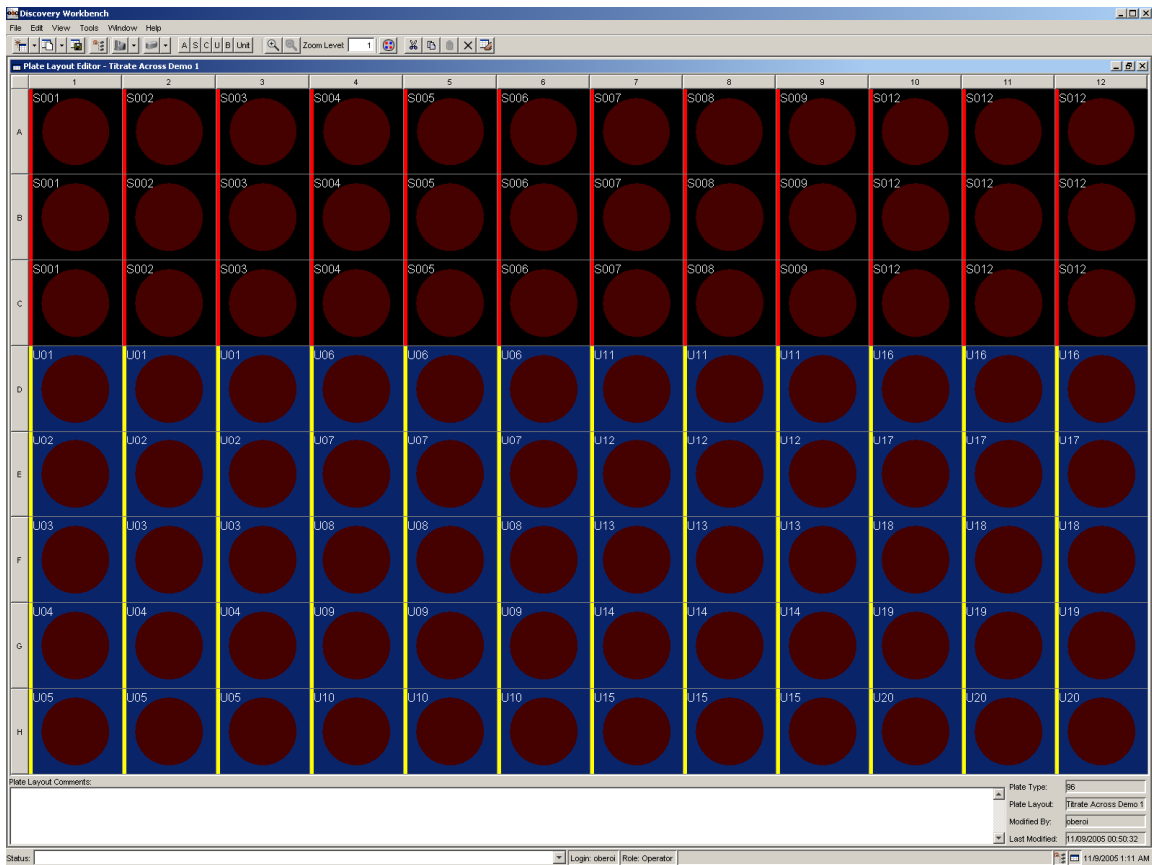
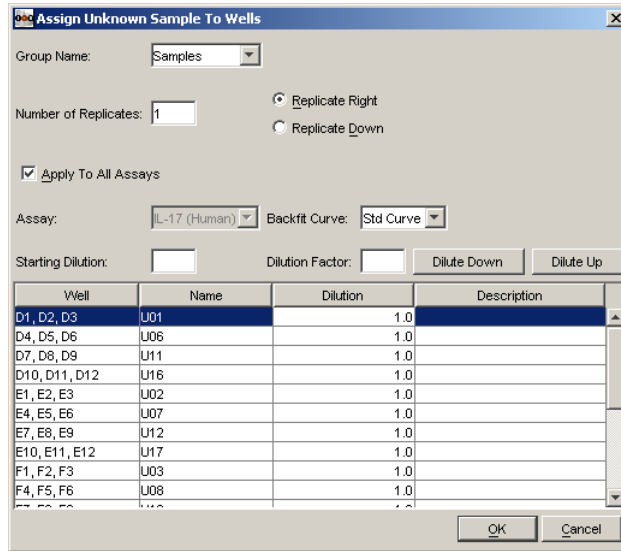
After clicking OK, the plate view should look like the following:



If however, we wanted the samples numbered going down rather than going across. For example, D1 should be U001, E1 should be U002 not U005.

There are two ways to change sample numbering.

First, new names can be assigned from the Assign Unknown Sample to Well Menu. Highlight the rows and select **U** from the tool bar. (Note that while the triplicates are shown, the window shows number of replicates as being 1. This is a bug in the current version of the program. If you need to re-assign replicates, delete the well assignment before creating the new assignment). From this window you can assign the new names by selecting the names in the table. See final view below.

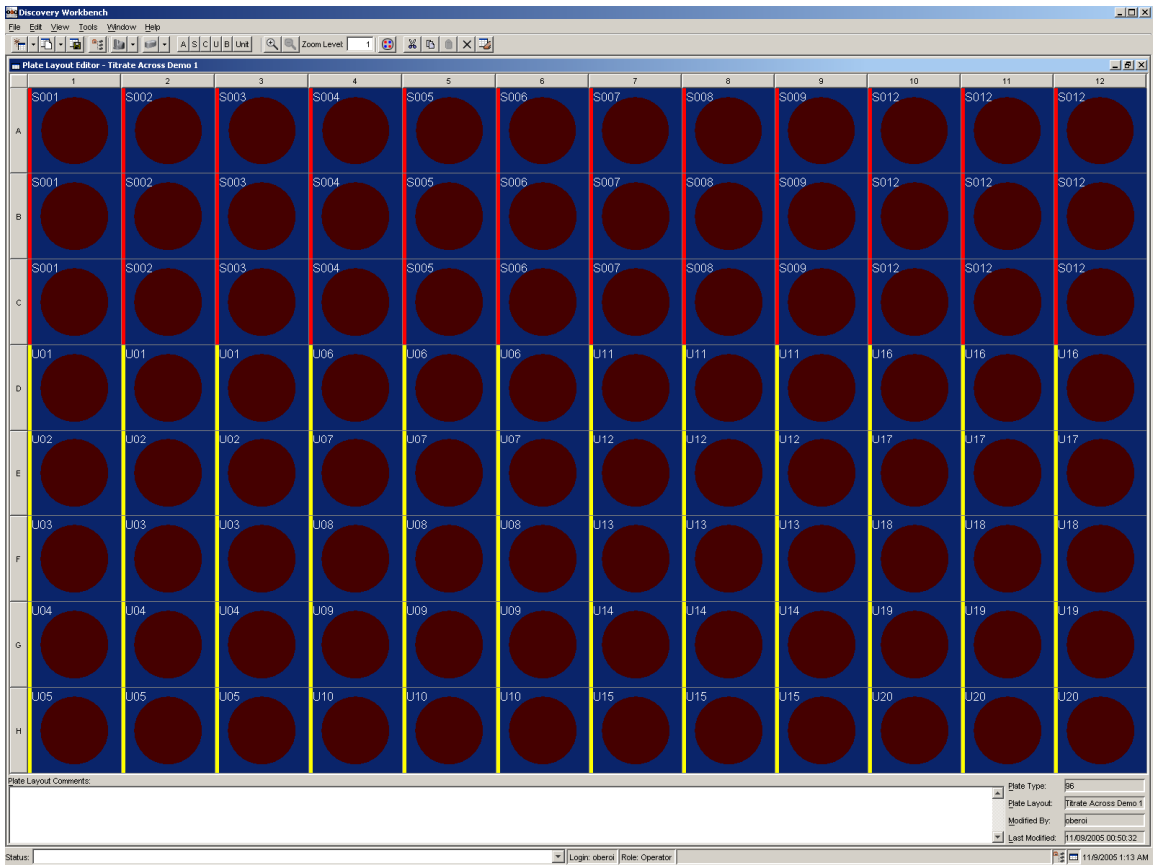


Alternatively, we can select the D1 through H3 and enter 3 for the number of replicates and then select Replicate Right then repeat for D4 through H6, D7 through H9 and D10 through H12. The results would be the same.

- Use one of these options to change the names to match the ones in the table shown above.

Step 2.5 Save the Plate Layout

The final Plate Layout should look like the following:

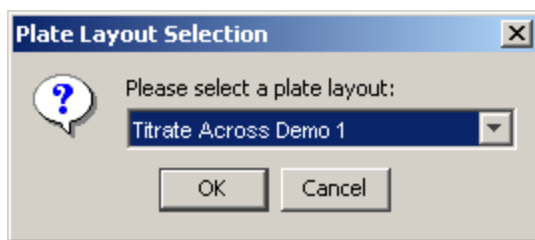


- Save the Plate Layout by going to **File->Save Plate Layout in Library** and then close the window by selecting the x in the upper right part of the plate view window. This will save the plate layout in the local Library of layouts.

4.2.3. Step 3. Create an Experiment Using the Plate Layout

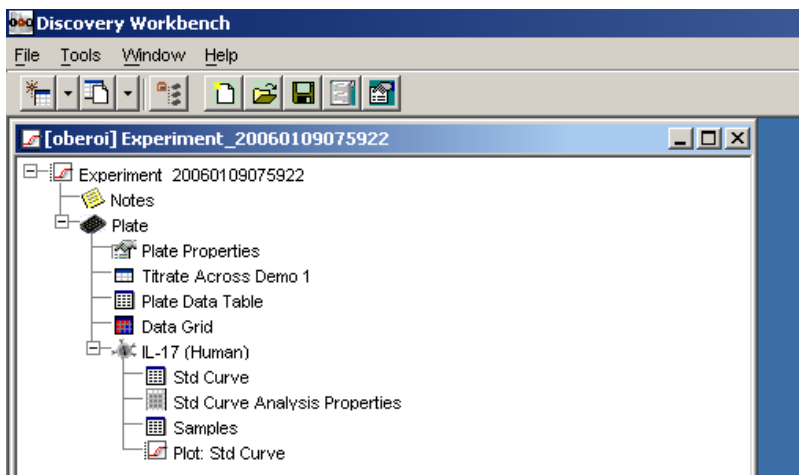
Now that you have the plate layout in the Library, the plate can be analyzed.

- Highlight the plate to be analyzed in the Plate Data History (For this exercise the plate is: *10/06/2005 15:23:27*). Select **File->Analyze Plates**. You may also right click on a highlighted plate, then select Analyze Plates.
- A pull down menu with the available Plates Layouts for that plate type will be shown.



Since we have only created one Layout for this type of plate, only one option will appear.

- Select *Titrates Across Demo 1* as the plate layout. Click OK, and an Experiment is automatically created as shown below. Expand the tree by clicking on the +.

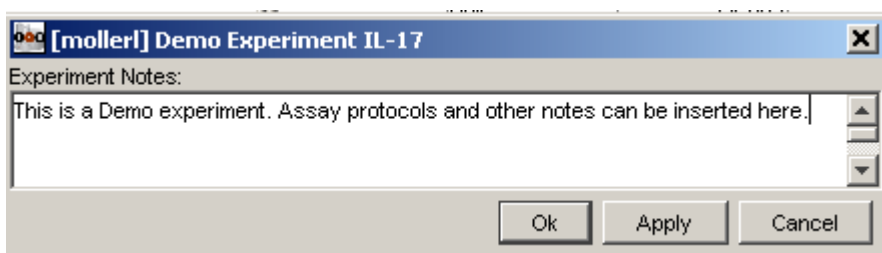


- Rename the Experiment by right clicking on: Experiment_20051109011512. Rename the Experiment to **Demo Experiment IL-17**. Then save the experiment using Ctrl-S or **File->Save Experiment**.

4.2.4. Step 4. Manipulating and Configuring the Experiment.

By double clicking on any of the items in the Experiment Tree, a window for that branch will open up. Since you can have a lot of windows open, it is generally good practice to close windows when they are no longer required for viewing.

When double clicking on the **Notes**, a Notes window will open up where you can record your notes. Click the apply button to save Notes and close the Notes window.



Viewing Results and Plots

Under the tree there will be a single plate. You can expand the branch by clicking on the +. Under the plate, there is the *Plate Properties*, the plate layout called *Titrate Across Demo 1*, *Plate Data Table*, *Data Grid*, and the assay.

- When you expand the IL-17 (Human) assay, 4 branches will appear:
 1. Standard: This is the table of data from the Standards.
 2. Standard Analysis Properties: This table contains the analysis statistics and detection limits.
 3. Unknown: This is a table of the unknowns.
 4. Plot: Standard: This is a plot of the standard curve.

The names of the groups are shown (Standard and Unknown). If more groups were created, more branches would be present.

The default table for standards is shown below:

Well	Concentration	Signal	Mean	CV	Calc. Conc. Mean	Calc. Conc. CV	% Recovery Mean
A01		928,794					
B01	10,000	915,146	916,742	1.24	10020	1.38	100
C01		906,286					
A02		244,055					
B02	2,500	241,550	241,098.667	1.33	2479	1.34	99.2
C02		237,891					
A03		64,993					
B03	625	58,560	60,521.333	6.41	631	6.29	101
C03		58,011					
A04		15,489					
B04	156.25	14,086	14,517.667	5.81	156	5.68	100
C04		13,978					
A05		3,352					
B05	39.062	3,340	3,380.667	1.78	37.1	1.78	95
C05		3,450					
A06		953					
B06	9.766	920	958.333	4.31	10.1	4.64	104
C06		1,002					
A07		307					
B07	2.441	362	351	11.3	3.11	15.1	127
C07		384					
A08		123					
B08	0.61	128	122	5.37	0.343	24.2	56.2
C08		115					
A09		96					
B09	0.153	102	106	11.8	0.137	119	89.5
C09		120					
A10		124					
B10		114					
C10		126					
A11		125					
B11	0	151	107.889	29.2	0.272	87.6	N/A
C11		123					
A12		66					
B12		86					
C12		56					

Notice that there are a lot of columns. The following columns are available:

Assay: Name of the Assay

Sample: Name of the Sample

Well: Well ID (A01, A02, ...)

Signal: Electrochemiluminescent Signal

Concentration: Concentration for standard, or expected concentration for controls

Mean: Mean signal for replicates

Std. Deviation: Standard deviation of the Signal

CV: Coefficient of variation for the Signal -- (std. deviation/mean)*100

Calc. Concentration: Backfit calculated concentration from the curve fit.

Calc. Concentration Mean: Mean of Calc. Conc.

Calc. Concentration Std. Deviation: Standard deviation of Calc. Conc.

Calc. Concentration CV: Coefficient of variation

% Recovery: For standards and controls percentage of expected concentration (Calc. Concentration/Concentration)*100

% Recovery Mean: Mean of % recovery for replicates

Detection Range: Sample in Detection Range, below or above detection range, above or below lowest/highest standard.

To add or remove columns, right click anywhere on the column titles then select or deselect columns from the menu. Columns can be moved around by dragging them with the left mouse button.

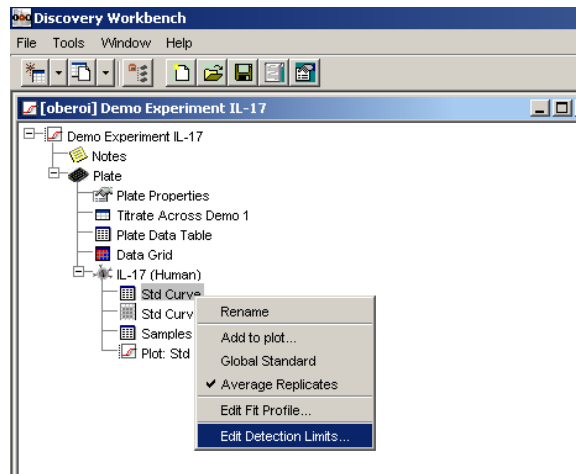
- Try to edit the table by selecting and deselecting appropriate columns such that the following columns are displayed: Well, Concentration, Signal, Mean, CV, Calc. Conc. Mean, Calc. Conc. CV and % Recovery mean.

Well	Concentration	Signal	Mean	CV	Calc. Conc. Mean	Calc. Conc. CV	% Recovery Mean
A01		928,794					
B01	10,000	915,146	916,742	1.24	10020	1.38	100
C01		906,286					
A02		244,055					
B02	2,500	241,550	241,098.667	1.33	2479	1.34	99.2
C02		237,691					
A03		64,993					
B03	625	58,560	60,521.333	6.41	631	6.29	101
C03		58,011					
A04		15,489					
B04	156.25	14,086	14,517.667	5.81	156	5.68	100
C04		13,978					
A05		3,352					
B05	39.062	3,340	3,380.667	1.78	37.1	1.78	95
C05		3,450					
A06		953					
B06	9.766	920	958.333	4.31	10.1	4.64	104
C06		1,002					
A07		307					
B07	2.441	362	351	11.3	3.11	15.1	127
C07		384					
A08		123					
B08	0.61	128	122	5.37	0.343	24.2	56.2
C08		115					
A09		96					
B09	0.153	102	106	11.8	0.137	119	89.5
C09		120					
A10		124					
B10		114					
C10		126					
A11		125					
B11	0	151	107.889	29.2	0.272	87.6	N/A
C11		123					
A12		66					
B12		86					
C12		56					

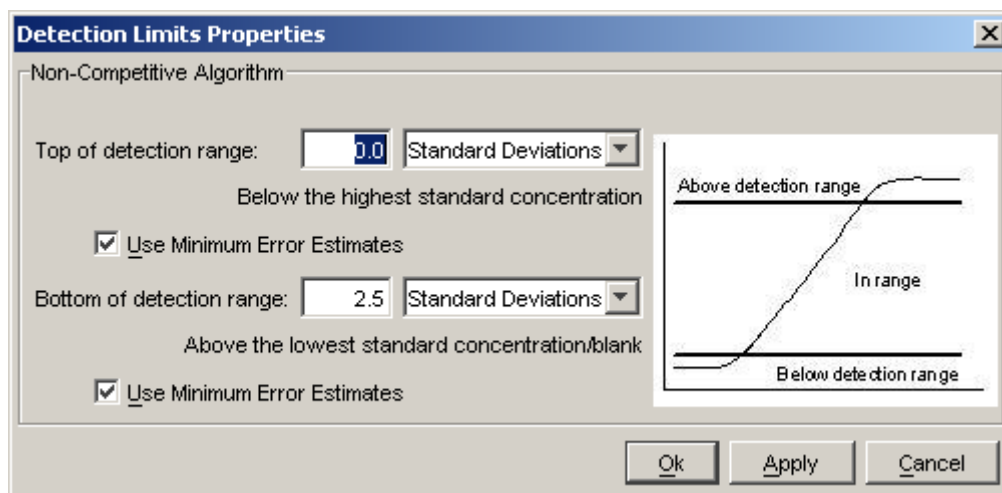
- Open the Sample Table and edit it such that the following columns are displayed: Sample, Detection Range, Well, Calc. Concentration, Calc. Conc. Mean and Calc. Conc. CV (as shown below):

Sample	Detection Range	Well	Calc. Concentration	Calc. Conc. Mean	Calc. Conc. CV
U01	In Detection Range	D01	542	523	3.31
	In Detection Range	D02	517		
	In Detection Range	D03	509		
U02	In Detection Range	E01	137	130	5.62
	In Detection Range	E02	131		
	In Detection Range	E03	122		
U03	In Detection Range	F01	37	35.7	3.05
	In Detection Range	F02	35.1		
	In Detection Range	F03	35.2		
U04	In Detection Range	G01	10.1	12.4	25.9
	In Detection Range	G02	11		
	In Detection Range	G03	16		
U05	Below Detection Range	H01	0.216	0.135	87.1
	Below Detection Range	H02	0.19		
	Below Fit Curve Range	H03	0		
U06	In Detection Range	D04	563	520	7.07
	In Detection Range	D05	501		
	In Detection Range	D06	497		
U07	In Detection Range	E04	145	135	6.15
	In Detection Range	E05	131		
	In Detection Range	E06	130		
U08	In Detection Range	F04	39	36.6	6.11
	In Detection Range	F05	36.1		
	In Detection Range	F06	34.7		
U09	In Detection Range	G04	10.7	10.2	4.75
	In Detection Range	G05	9.9		
	In Detection Range	G06	9.89		
U10	Below Detection Range	H04	0.229	0.19	29.6
	Below Detection Range	H05	0.216		
	Below Detection Range	H06	0.126		
U11	In Detection Range	D07	535	491	7.61
	In Detection Range	D08	471		
	In Detection Range	D09	469		
U12	In Detection Range	E07	126	125	1.81
	In Detection Range	E08	124		
	In Detection Range	E09	124		
U13	In Detection Range	F07	37.3	34.5	7.08
	In Detection Range	F08	33.3		
	In Detection Range	F09	32.8		
U14	In Detection Range	G07	10.2	9.43	9.19
	In Detection Range	G08	9.57		
	In Detection Range	G09	8.5		
U15	Below Detection Range	H07	0.033	0.095	78.8
	Below Detection Range	H08	0.073		
	Below Detection Range	H09	0.178		
U16	In Detection Range	D10	575	545	5.65
	In Detection Range	D11	547		
	In Detection Range	D12	514		
U17	In Detection Range	E10	151	144	4.33
	In Detection Range	E11	142		
	In Detection Range	E12	139		
U18	In Detection Range	F10	40.1	38.5	3.78
	In Detection Range	F11	38		
	In Detection Range	F12	37.4		
U19	In Detection Range	G10	11.4	11.1	1.83
	In Detection Range	G11	11		
	In Detection Range	G12	11.1		
U20	Below Detection Range	H10	0.343	0.347	45.6
	Below Detection Range	H11	0.19		
	Below Detection Range	H12	0.506		

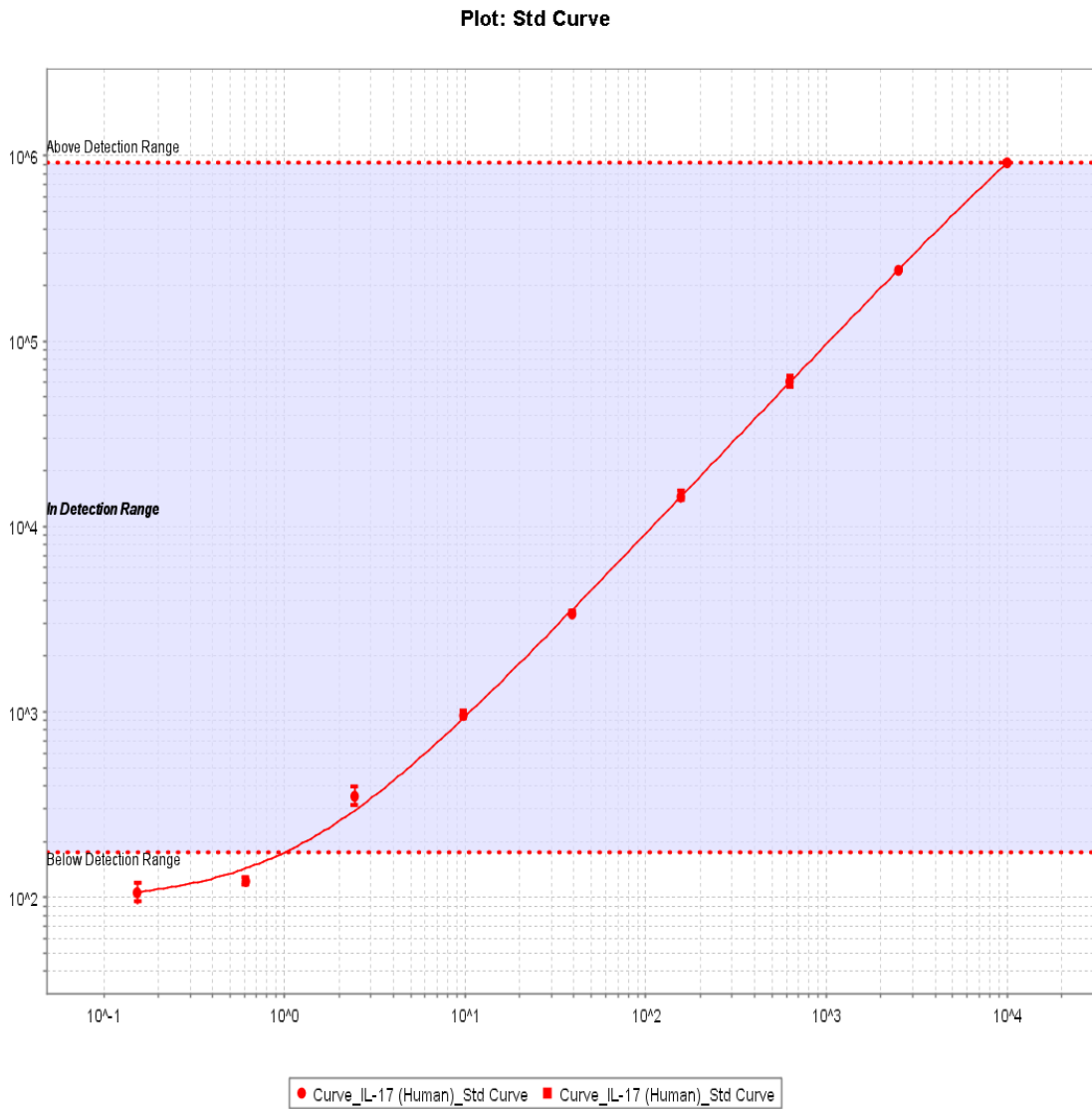
Notice the new column shown called Detection Range. The detection limits can be set by selecting the Std Curve branch in the tree and right clicking to open a pull down menu and selecting **Edit Detection Limits**.



The upper and lower limits of detection can be set as a number of standard deviations over the bottom and the top of the curve in the Edit Detection Limits window. The default is 2.5 standard deviations from the bottom and the top calibrator.

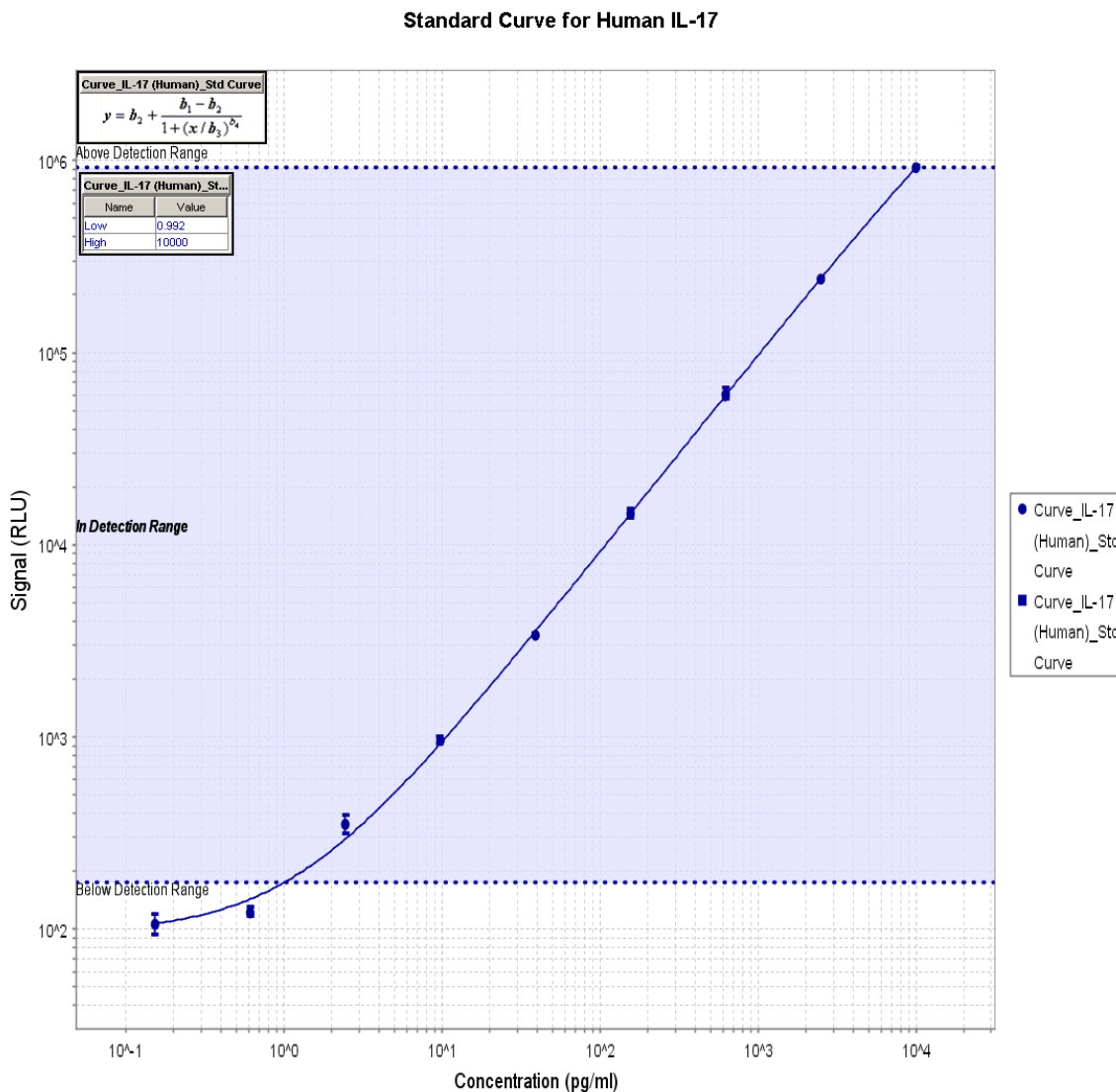


- Double click on the plot icon in the tree, and a plot window should appear, as shown below:



Right clicking in the plot gives a large pulldown menu. From this menu, one can edit the colors of the graphs by selecting *Edit Plot*. The axis can be labeled and scaled by selecting *Edit Plot Settings*. The plot can be copied to the clipboard for pasting into other documents. The image can be saved. Annotations in the graph and the analysis parameters can be added to the graph.

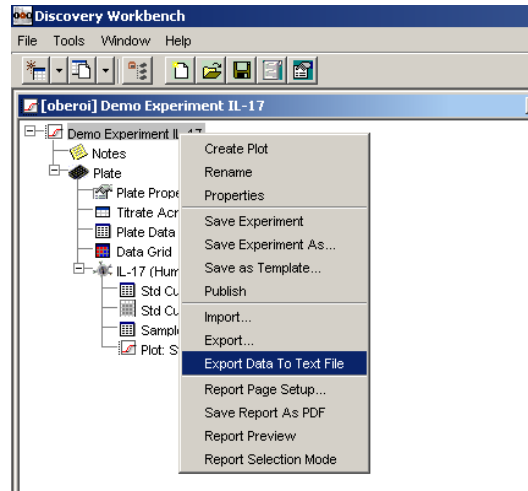
- Try to create the following plot by changing the color of the curve and symbols, moving the legend, editing the title, naming the axes and displaying the detection limits and fitting function:



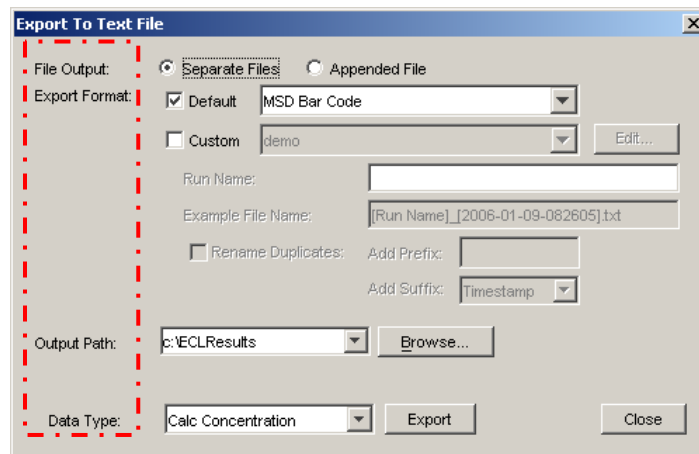
4.2.5. Step 5. Saving, reporting, and copying plots and tables.

The calculated concentrations for the plate can be exported in formats similar to the text files that are generated by the instrument in signal space.

- To create a text file with the concentrations rather than signals, right click on the experiment in the tree and select **Export Data To Text File**.



This will open a window where you can select the **Data Type** to be exported, **Export Format**, **Output Path**, and whether multiple plates should be appended or as separate files. Click on the Export button to export the data.

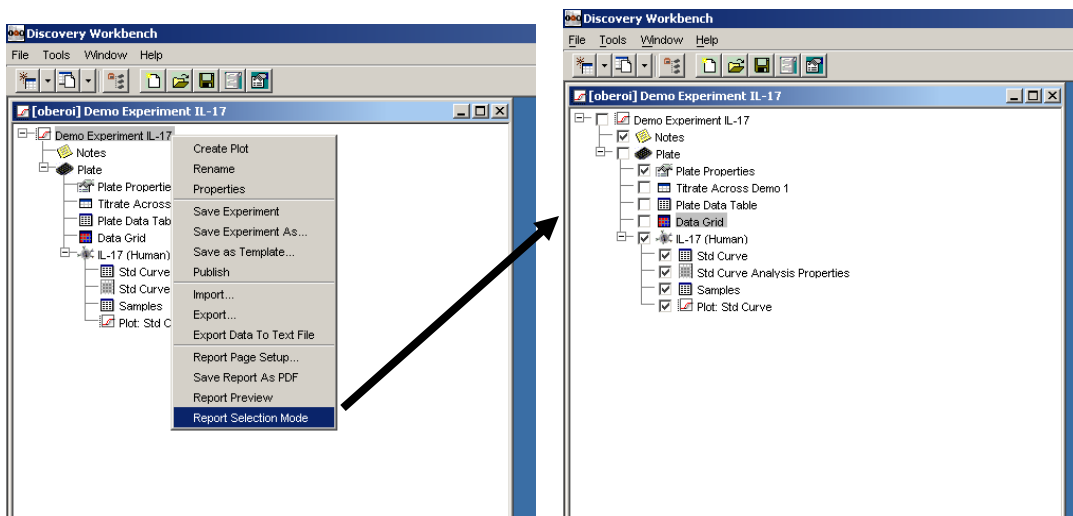


- Open the text file generated by the export function. It should display concentrations rather than the signals.

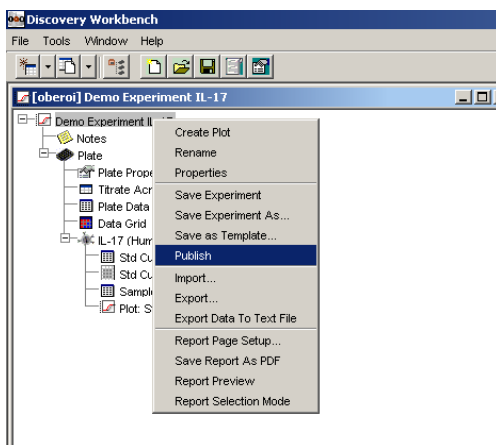
A PDF report of the Experiment can be saved for printing purposes. The number of pages may be very large if the whole experiment is printed, so the user can select the parts of the experiment to be exported.

- For this experiment, try saving the PDF in two ways. First select **File -> Save Report as PDF** and save the PDF as **Full_IL-17Report.PDF**. The report should be 17 pages long.

To print selected portions of the experiment, right click on the experiment and select **Report Selection Mode**. Check boxes will now appear in the tree structure. Select the portions of the experiment to be printed. For example in the right part of the figure below, the Notes, Plate Properties, Data tables, Analysis Properties and Plot are printed. Using this example, a 9 page report will be generated.



Experiments can be published by selecting Publish from the right click menu on the Experiment. Once published, an experiment cannot be modified. The experiment can be opened and saved as a new experiment in the Plate Data History, but a new name must be assigned to the experiment.



Data can be copied from tables to other Windows programs by cutting and pasting. For example, the entire data set for the plate is present in the Plate Data Table. Open this table and select **Edit -> Copy Table to Clipboard**. This will copy that entire table to the Windows Clipboard. The data table can be pasted into another Windows program

such as Excel or Word. The data can be manipulated in these other programs to create different tables or summary tables. The image of the table can be copied by selecting ***Edit ->Copy Image to Clipboard***. When this is copied to another windows program, the data cannot be manipulated and is just an image of the table. Plots can be copied the same way. This function is useful for making presentation and putting tables and graphs in reports.

4.2.6. Advanced Functions: Excluding Data

Data can be excluded from the curve fit at multiple different screens: the plot, the table, or the Data Grid. Excluded data on a plot will show up as an open symbol. If the symbol is a dot, it will appear as an open circle. Excluded data in a table will appear in a red font. Excluded data in the Data Grid will appear grayed out.

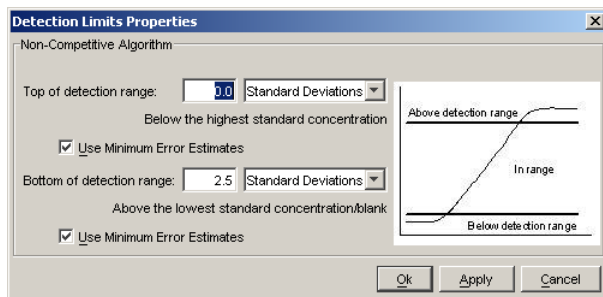
Data can be excluded from a plot by right clicking on a point on the curve and selecting **Exclude Data Point(s)**. The data symbol will change and the curve should automatically adjust to the exclusion.

- From the plot, exclude the data point at 2.441 pg/ml. Before excluding, notice the detection limits from the Analysis Properties (right) after exclusion, the detection limits are slightly higher (left).

Std Curve Analysis Properties		Std Curve Analysis Properties	
Name	Value	Name	Value
Algorithm Parameters		Algorithm Parameters	
Initial Top	925909	Initial Top	925909
Initial Bottom	95.4	Initial Bottom	95.4
Initial MidPoint	4912	Initial MidPoint	4912
Initial HillSlope	1	Initial HillSlope	1
Weighting	1/y^2	Weighting	1/y^2
Max Iteration	500	Max Iteration	500
Fit Statistics		Fit Statistics	
RSquared	1	RSquared	1
Calculated Parameters		Calculated Parameters	
Top	7135571	Top	4561776
Bottom	95.6	Bottom	96.3
MidPoint	64245	MidPoint	37607
HillSlope	1.03	HillSlope	1.05
Detection Range Parameters		Detection Range Parameters	
Low	0.992	Low	1.11
High	10000	High	10000
Equation		Equation	
FourPL	$y = b_2 + \frac{b_1 - b_2}{1 + (x / b_3)^{b_4}}$	FourPL	$y = b_2 + \frac{b_1 - b_2}{1 + (x / b_3)^{b_4}}$

4.2.7. Advanced Functions: Detection Limits

Detection limits are calculated from the curve fit of the standard curve. The detection limits are by default to be 2.5 standard deviations from the bottom of the curve and the top is the top calibrator. These limits can be modified by right clicking on the **Standards Table**. The Detection Limits Properties box allows one to select the number of standard deviations from the highest and lowest standards to compute the top and bottom of the detection range.



When graphs are generated, they automatically have the detection range shaded in grey to be able to see the points that fall into the detection range. A column can be added in the tables to show whether the data is:

- a) **Below Fit Curve Range** – Signal is below the bottom of the bottom of the curve fit. No concentration is given.
- b) **Below Detection Range** – Signal is above the bottom of the curve, but below the detection limit as defined by the Detection Limit Properties.
- c) **Above Detection Range** – Signal is above the top of detection range as defined by the Detection Limit Properties.
- d) **Above Fit Curve Range** – Signal is above the top of the curve fit. No concentration is given.
- e) **In Detection Range** – Signal is within the top and bottom detection limits and a concentration is provided.

(See step 4 for a review of how to insert columns in a table).

Average of Replicates

By default, the curve fitting is done using the average of the replicates of each standard. To change to using individual points for the curve fitting, right click on the standards table and uncheck Average Replicates. Individual points will be displayed for the standards on the plots when the Average Replicates option is not selected.

4.3. Example 3: 10-spot Mouse Cytokine Panel

The data file: Example3.txt is a titration of the mouse cytokine with unknown samples below. The standard curve consists of a 1:4 dilution series of calibrators from 10,000 pg/ml to 0.4 pg/ml across the top 4 rows in quadruplicate. The spots are:

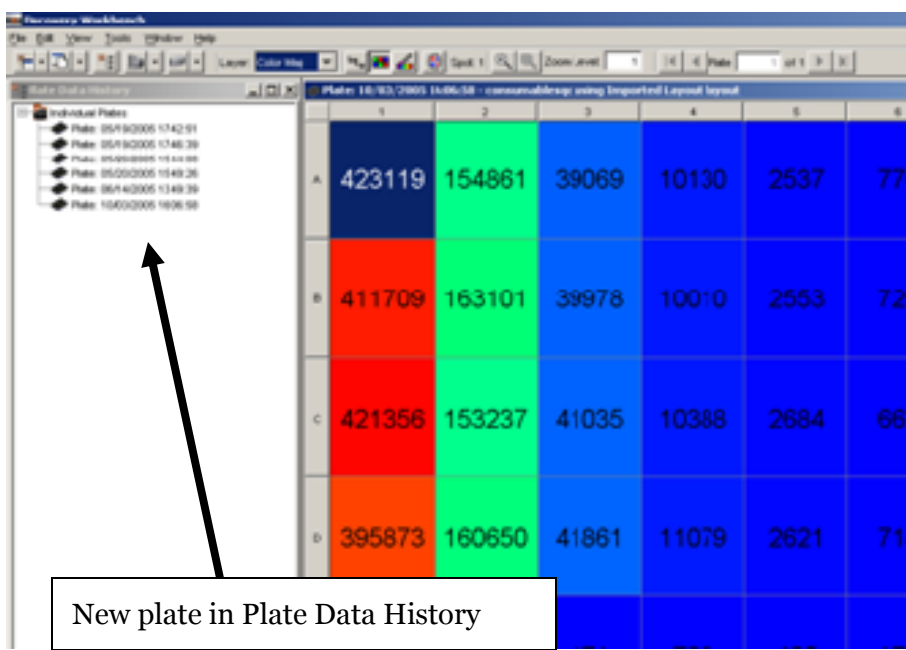
- 1: mIFNg 6: mKC
- 2: mIL-1b 7: mIL-10
- 3: mIL-2 8: blank
- 4: mIL-4 9: blank
- 5: mIL-5 10:mTNFa.

The unknowns are assayed in quadruplicate (replicated down) in the bottom 4 rows.

4.3.1. Step 1: Import the data file into the Plate Data History.

Open MSD Discovery Workbench.

Open the Plate Data History by selecting the icon, or selecting it from the **Tools** menu at the top. Once the Plate Data History is open, select **File->Import->Plate Data File**. Then select the file *Example3.txt*. A new plate should open up and the data window should open and a plate created on **10/03/2005 16:06:58** should be created in the Plate Data History. The plate is a Sector Imager 6000, 10-spot plate.

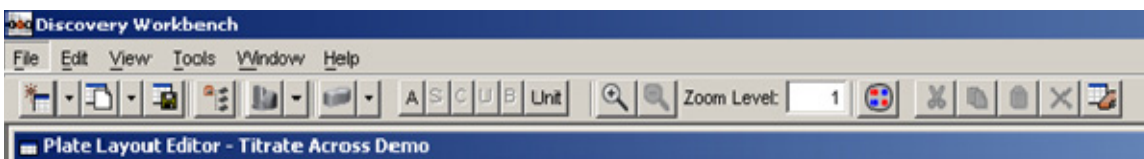


4.3.2. Step 2: Create Plate Layout

Close the plate window before continuing. Create a new plate layout for this plate and add it to the library of plate layouts and experiment templates in the Discovery Workbench.

Step 2.1: Select Plate Type

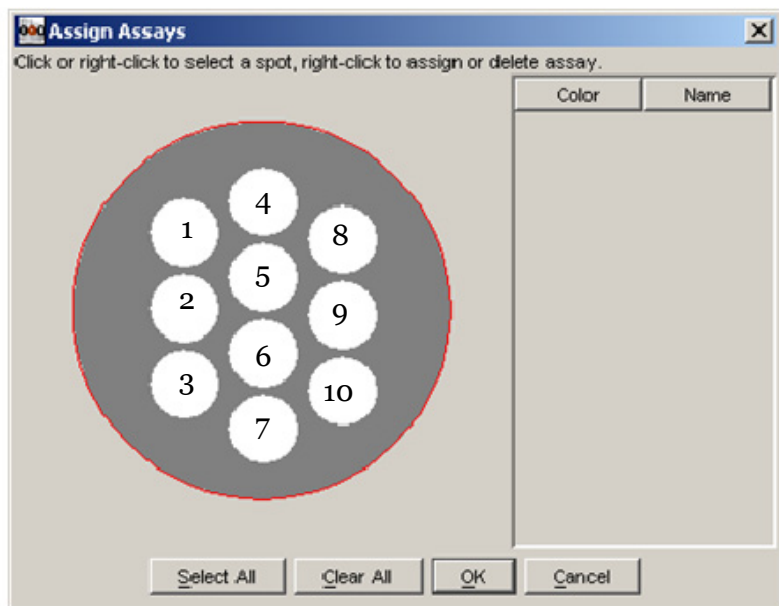
- From the **File** menu, select **File->New Plate Layout in Library->Sector Imager 6000**. Select *96 Multi-Spot 10* from the **Plate Type** dropdown menu. For the name, type in: *Titrate Across Demo*. Click on the New Button. This should open a 10-spot plate view with all of the spots being white. The Toolbar shown below should appear on the top of the window:



Step 2.2 Select Assays (A)

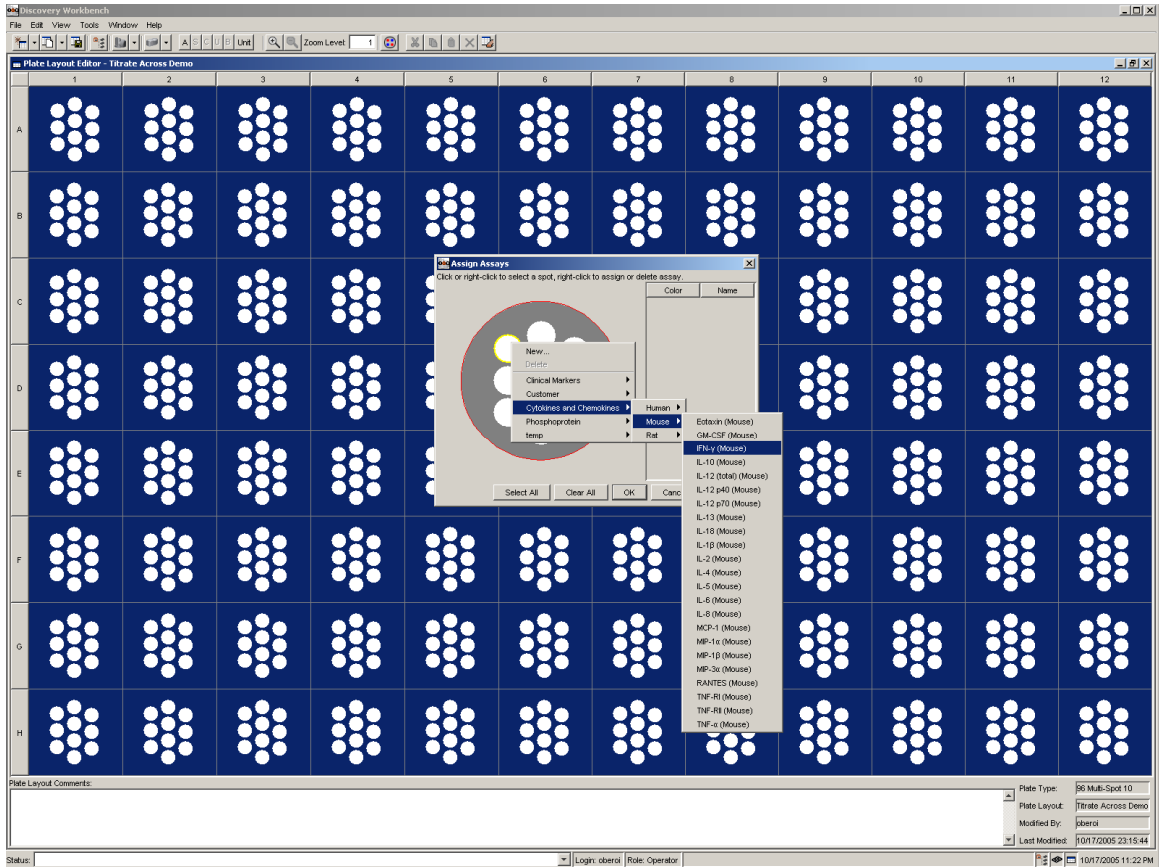
The **A** button in the toolbar is used to assign the assays for the spots.

- Left click on the **A**, and the **Assign Assays** window shown below should appear:



Note: there are no numbers shown in the **Assign Assays** window, you can find the spot assignment on the foil package of the MSD plate used in the assay.

- To assign a spot, right click on the spot and a drop down menu will open up. Right click on spot 1 (top left) and from the drop down menu, select **Cytokines and Chemokines->Mouse->IFN- γ (Mouse)**.

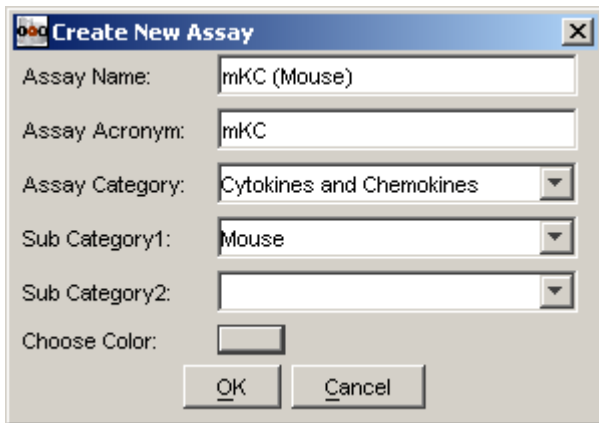


Repeat to select the other spots.

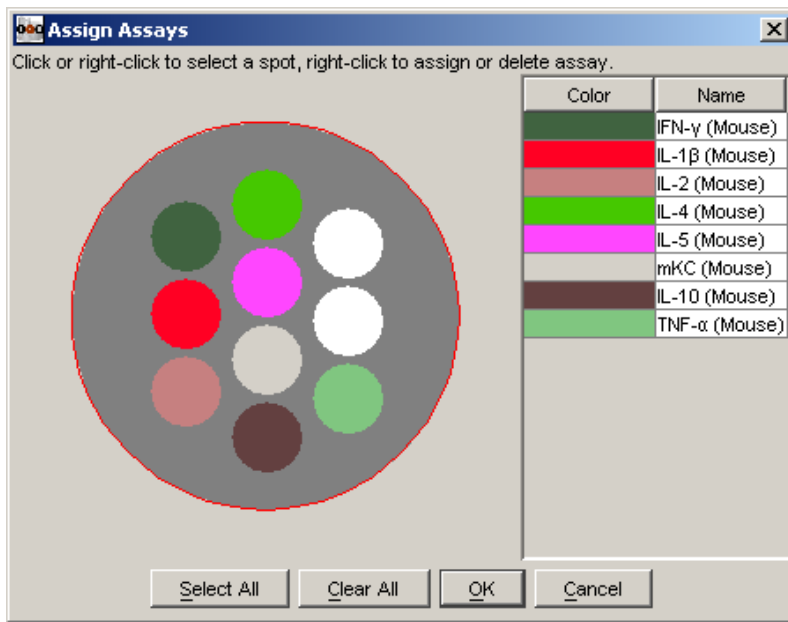
The spots are: 1: mIFNg 6: mKC
 2: mIL-1b 7: mIL-10
 3: mIL-2 8: blank
 4: mIL-4 9: blank
 5: mIL-5 10:mTNFa.

For the blank spots, do not assign any assays.

Notice that there is no mKC assay. To create a new assay, select a New when clicking on the spot. Create a new mouse KC assay as shown in the window below.



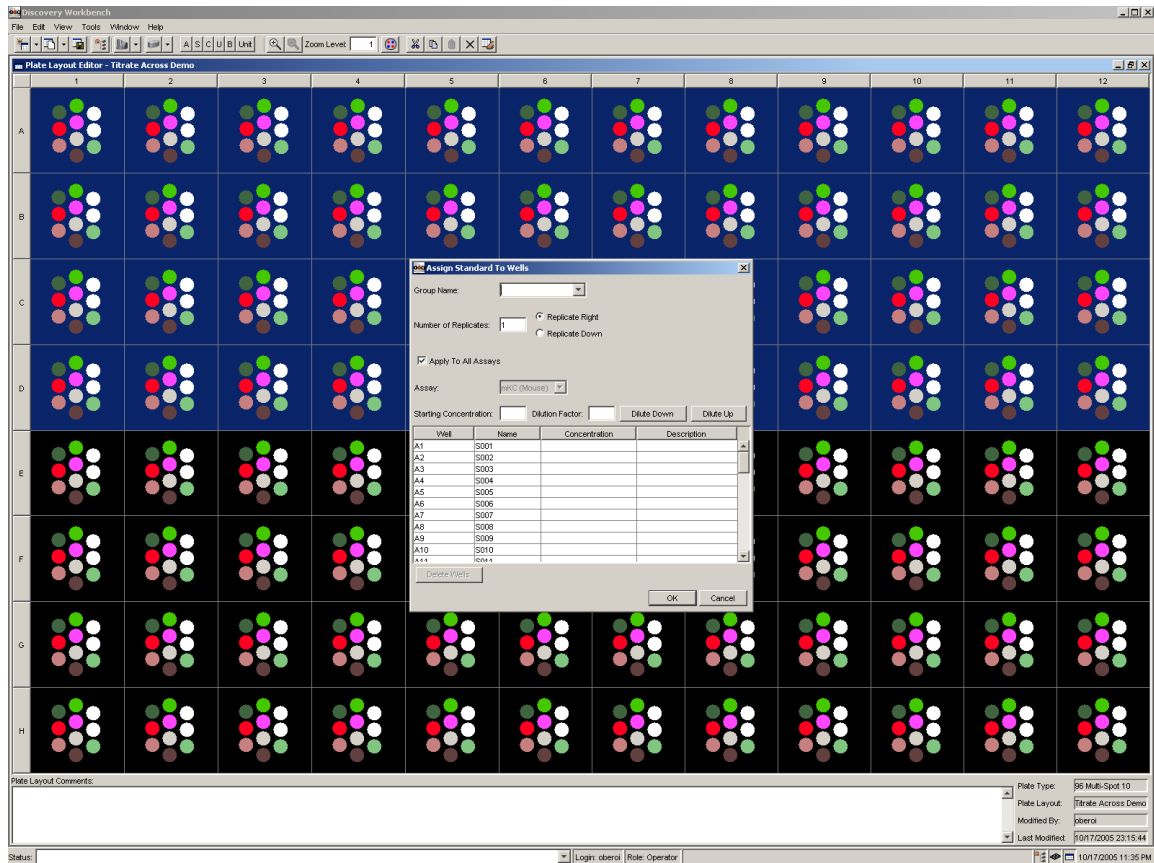
After you have assigned all the spots, the **Assign Assays** window should look like the one below. Notice the legend and colors represent which assays are assigned to which spots.



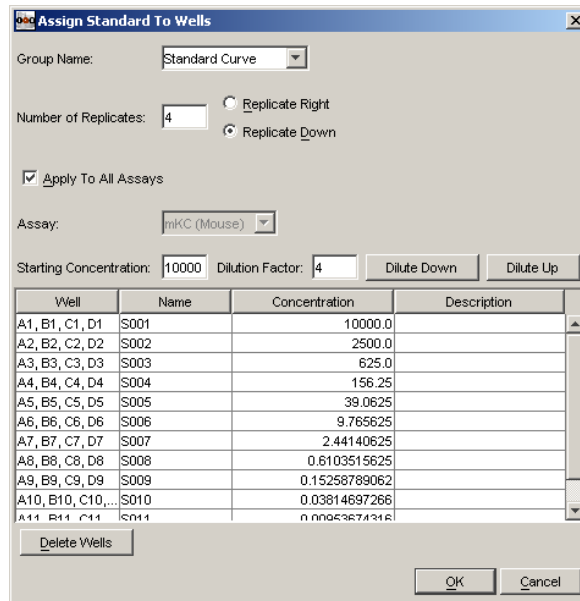
After assigning the assays, click OK and you're ready to move on to assigning standards, controls, unknowns and blanks.

Step 2.3 Assign Standards (S)

For this example, select rows A, B, C, and D. Then click on the **S** button in the tool bar to open the **Assign Standards to Wells** window. A representative screen shot is shown below:



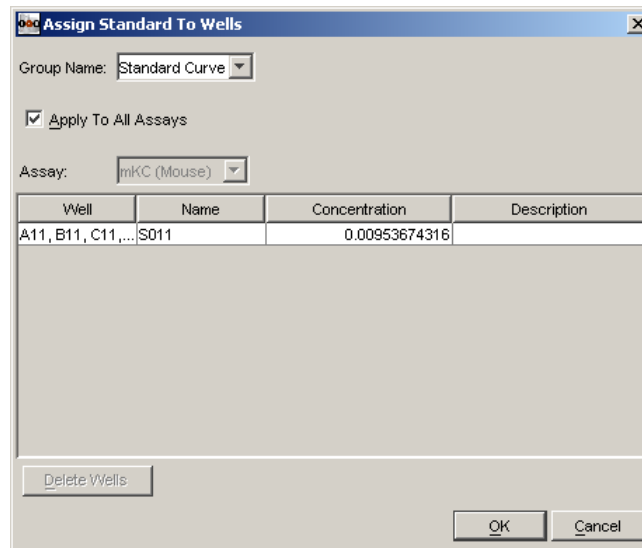
- Complete the **Assign Standard to Wells** window. In the *Group Name* type in: *Standard Curve*. Since this example has 4 replicates, change the number of replicates to 4 and select *Replicate Down*. The wells should now be grouped A1,B1,C1,D1 and A2,B2,C2,D2,... Since the standard curve starts at 10,000 pg/ml and we diluted across by a factor of 4, put the starting concentration at 10,000 and dilution factor of 4, then click on the Dilute Down button. Scroll down to S012 and overwrite the concentration to be 0. Note that **Name**, **Concentration**, and **Description** can be modified in the table as well. The final window should look like the following:



Click OK and the Plate layout should now have a color description with the sample names in the well.

You've just learned that the concentration of calibrator in column 11 is zero.

- How can we change the concentration of A11, B11, C11, and D11 to zero? Select just A11 and click on the **S** button to open the Assign Standard to Wells window. Notice that the window opened all the wells with the S011 name:



- Overwrite the concentration to be 0 and Click OK. The problem is that now we have 2 different standards with 0 concentration. This is OK, but then we will not get the appropriate CVs and standards curves for the zero. So we want to copy the

S011 standard to wells A12, B12, C12, and D12. To do this, select A11 through D11 as a block. Then copy with *Ctrl-C* or Copy from the Edit menu. Then select A12 through D12 as a block and paste using *Ctrl-V* or Paste from the Edit menu. Now there are 8 wells with the 0 standard.

Step 2.4 Assign Unknowns (U)

The bottom 4 rows contain unknowns assayed in quadruplicate. Select rows E, F, G, and H and then click on the **U** button in the tool bar to open the Assign Unknowns to Wells window. This step is similar to assigning the standards, with a few exceptions. First, one can select which backfit curve to fit the unknowns to if there are multiple standard curves on the plate. Second, one can assign the dilution rather than the concentration. The calculated concentrations will take into account the dilution factor.

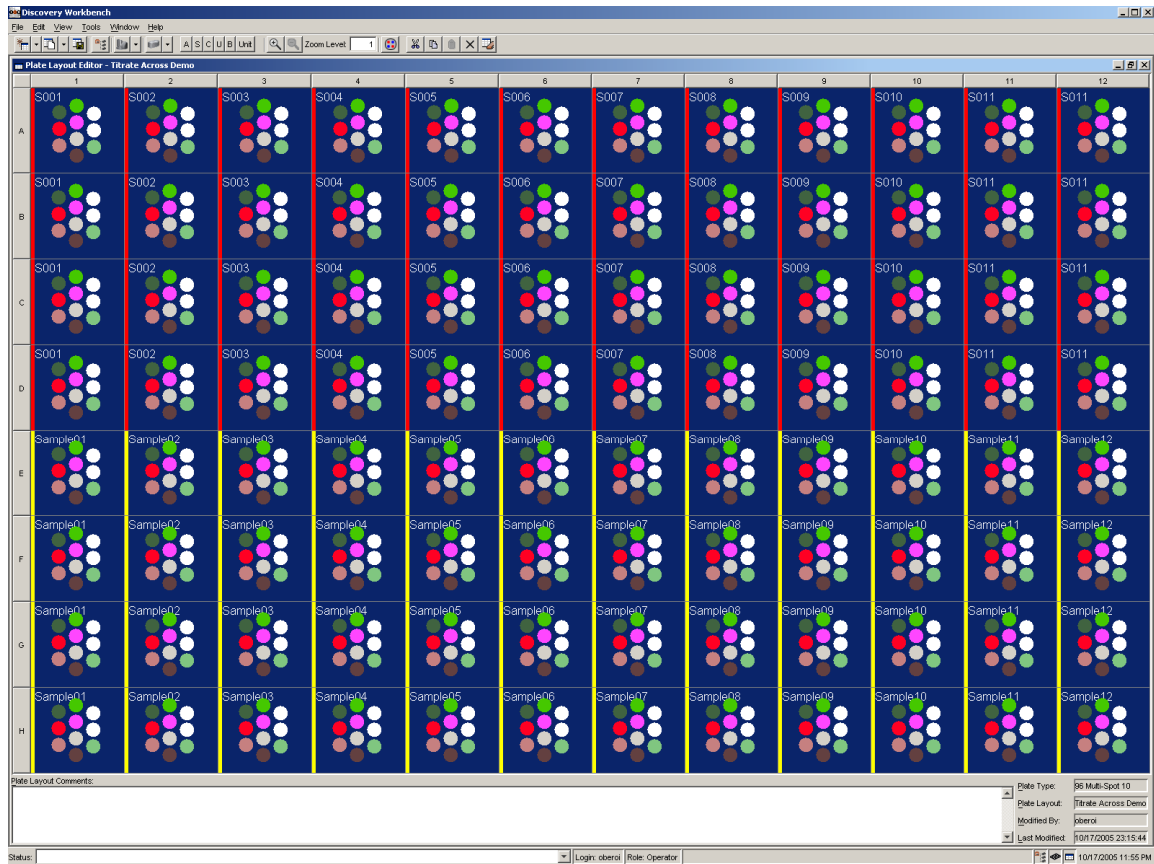
Try to generate the following on your own:

Well	Name	Dilution	Description
E1, F1, G1, H1	U001	1.0	
E2, F2, G2, H2	U002	1.0	
E3, F3, G3, H3	U003	1.0	
E4, F4, G4, H4	U004	1.0	
E5, F5, G5, H5	U005	1.0	
E6, F6, G6, H6	U006	1.0	
E7, F7, G7, H7	U007	1.0	
E8, F8, G8, H8	U008	1.0	
E9, F9, G9, H9	U009	1.0	
E10, F10, G10, H10	U010	1.0	

Let's assume for this example that the samples have names. We can change the names from U001, U002, ... U012 to Sample01, Sample02...Sample12 by typing in the Name in the table. Note that we pad the number with a 0 so that the samples can be sorted appropriately.

Step 2.5 Save the Plate Layout

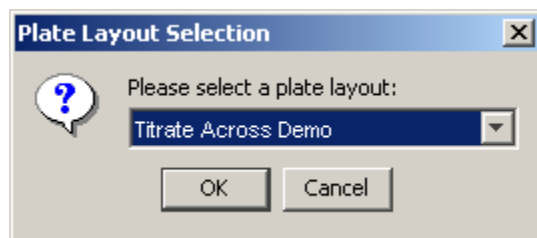
The final Plate Layout should look like the following:



Save the Plate Layout: **File->Save Plate Layout in Library** and then close the window.

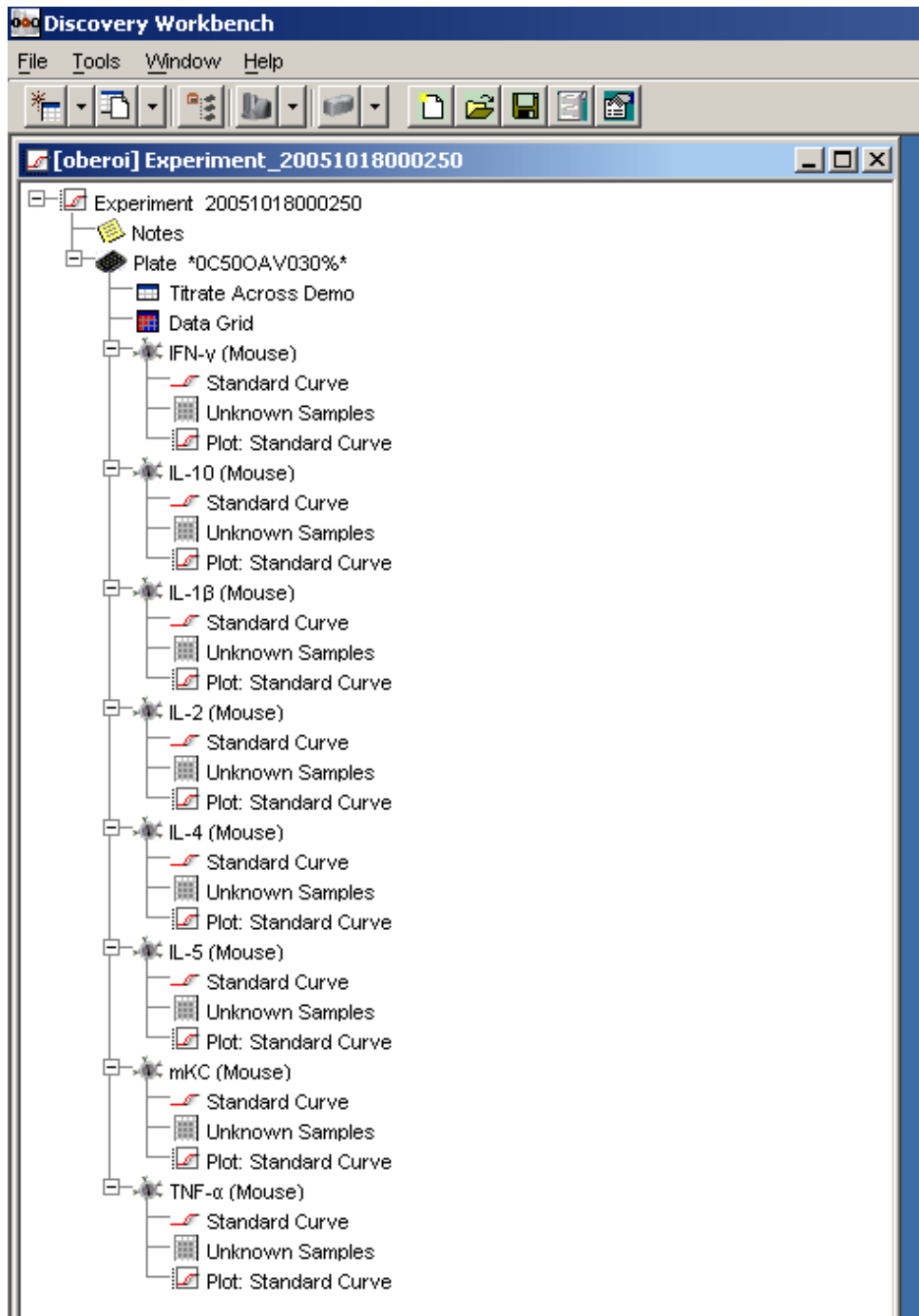
4.3.3. Step 3. Create an Experiment Using the Plate Layout

Now that you have the plate layout in the Library, the plate can be analyzed. Highlight the plate to be analyzed in the Plate Data History (**10/03/2005 16:06:58**). Select **File->Analyze Plates**. A pull down menu with the available Plates Layouts for that plate type will be shown.



Since we have only created one Layout for this type of plate, only one option will appear.

- Select *Titrates Across Demo* as the plate layout. Click OK, and an Experiment is automatically created as shown below. Expand the tree by clicking on the +.



- Rename the experiment by right clicking on: Experiment 20051018000250. Rename the experiment to ***Demo Experiment 10 Spot***. Then save the experiment by Ctrl-S or **File->Save Experiment**.

4.3.4. Step 4. Manipulating and Configuring the Experiment.

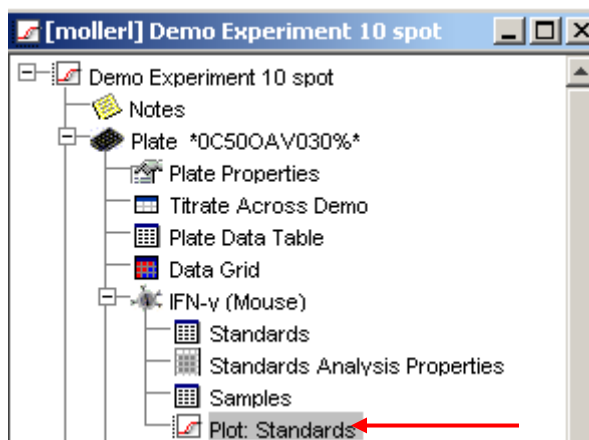
By Double clicking on any of the items in the Experiment Tree, a window for that branch will open up. Remember to limit the number of windows open at the same time, as having too many windows open could cause the computer to perform sluggishly.

4.1 Displaying samples on a plot

The distribution of samples relative to the standards can quickly be visualized on a plot.

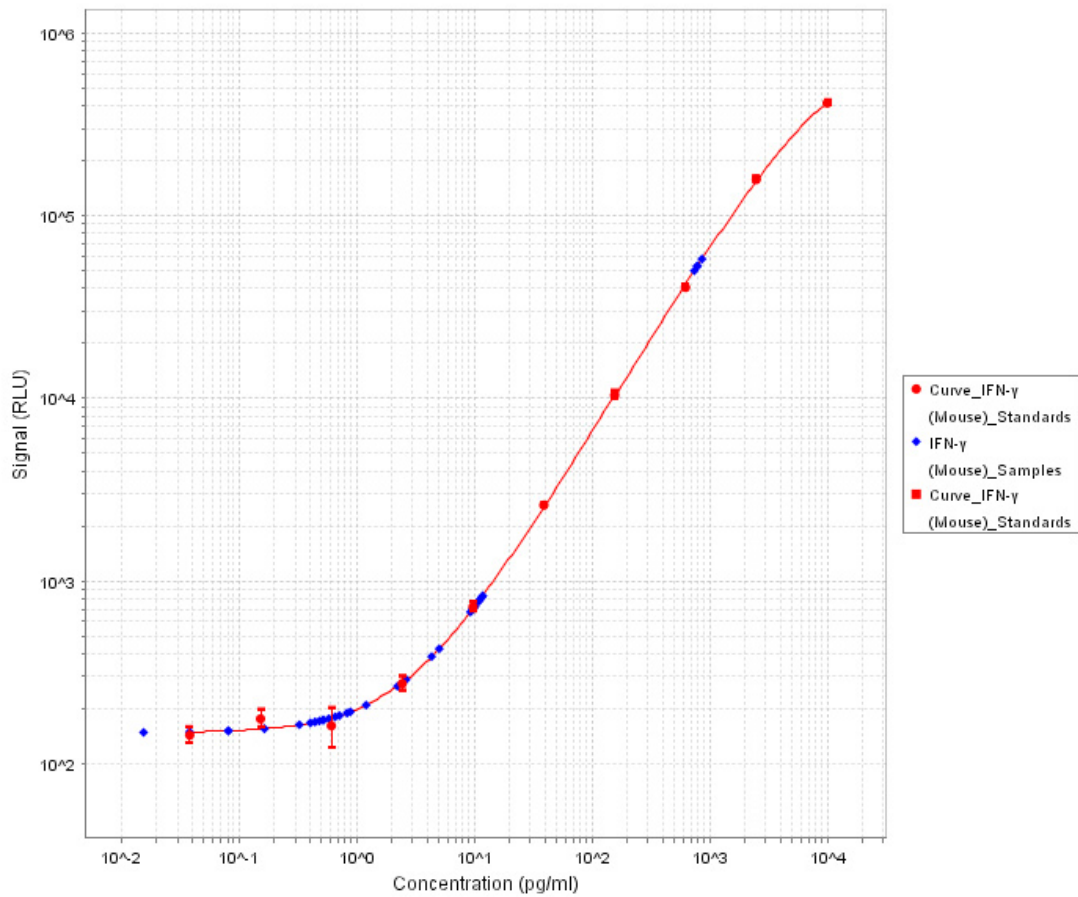
We will use the IFN-g assay as an example.

- Open the IFN-g plot by double clicking on Plot: Standards



- Highlight the Samples table for IFN-g and with the left mouse button drag the table onto the plot window. The samples should now appear on the plot as shown below.

Multiple groups (if available) can be dragged and dropped into the plot window.

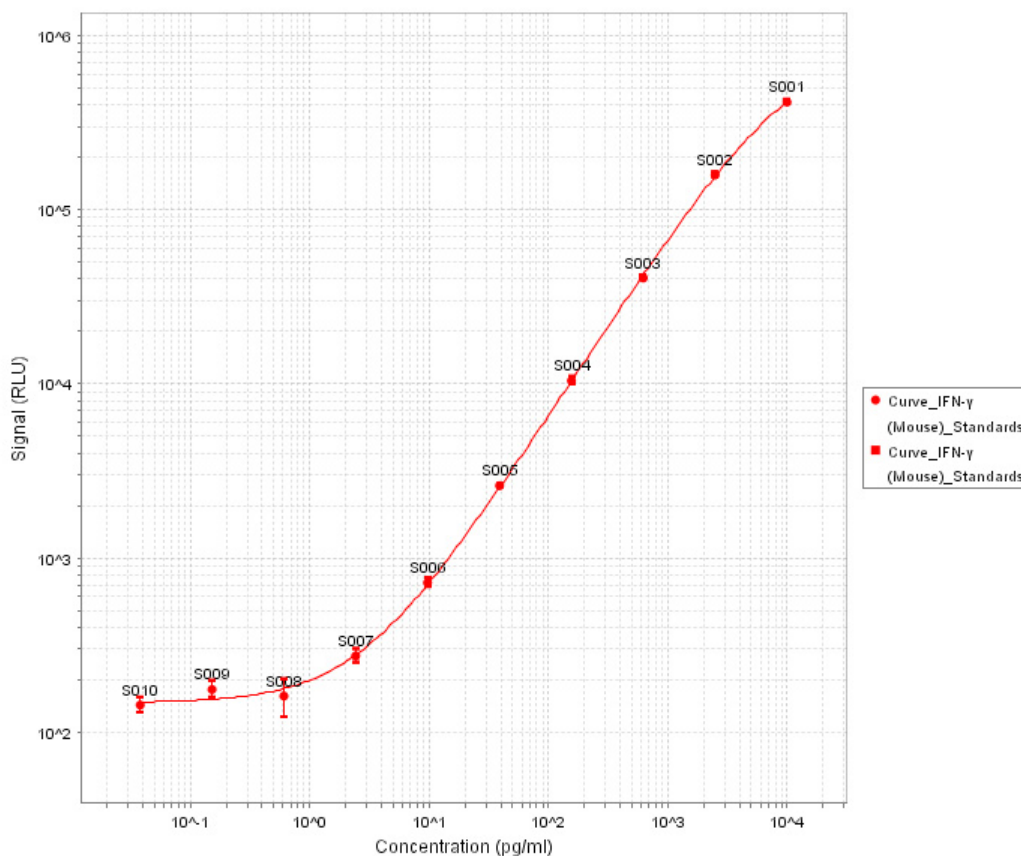


To remove items from a plot, either right click the plot and select the *Remove From Plot* option from the menu or go to *Edit > Remove From Plot*.

- Remove the samples from the plot

4.2 Displaying sample names

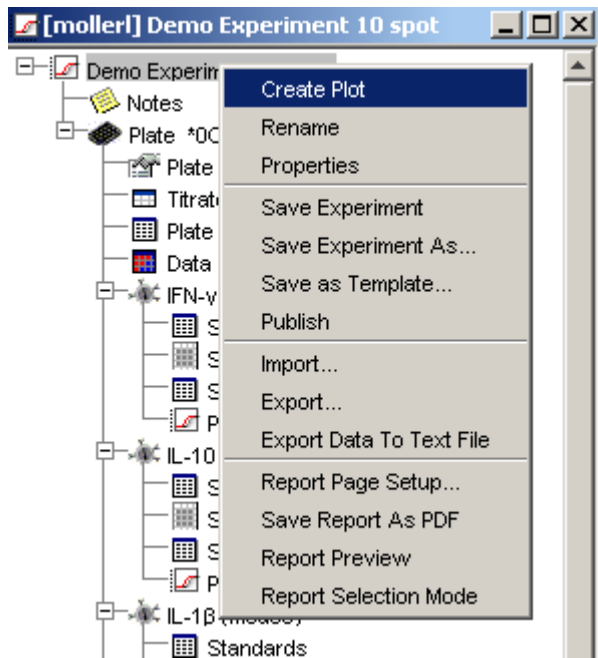
- To display sample names in a plot go to *View -> Sample Annotation*. Add the names of the standards to the IFN-g plot as shown below.



4.3 Creating plots

In addition to the plots available for each assay, custom plots can easily be created.

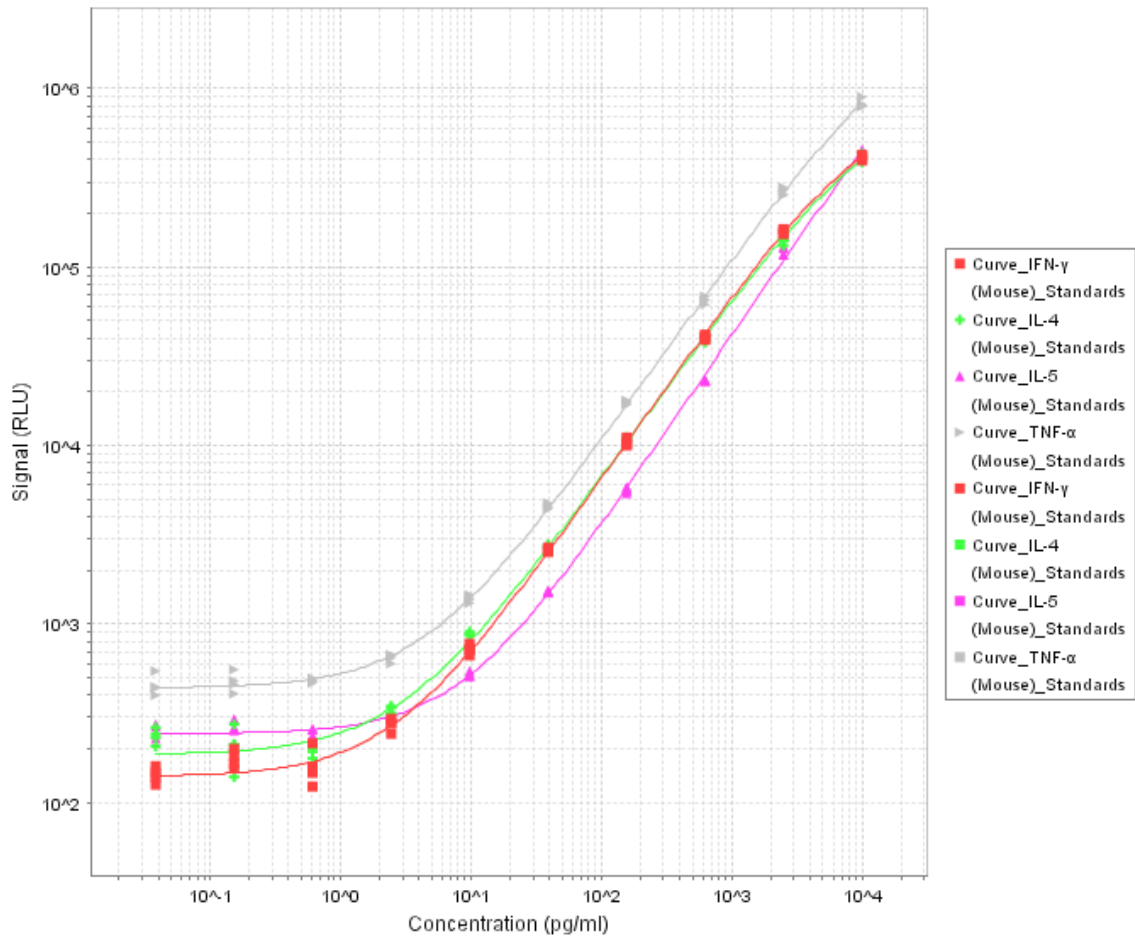
- To create a new plot, highlight the experiment and select *Create Plot* from the right click menu. A blank plot window will appear.



We will create a plot of the standard curves for IFN-g, IL4, IL5 and TNF-a.

- Highlight the Standards table for IFN-g and with the left mouse button drag the table onto the empty plot window. The IFN-g calibration curve should now appear on the plot. Repeat this step for IL4, IL5 and TNF-a.
- Try to edit the plot in the following manner:
 - Remove the detection ranges
 - Use individual replicates instead of the average of replicates
 - Editing the title and axes so that it looks like the one shown below

Standard Curves



5. Frequently Asked Questions

5.1. Layouts

5.1.1. How do I assign assays?

- 1) To assign the same assay pattern to all wells in the plate, right click on a layout in the layout editor and select Assay Assignment.
- 2) To assign an assay pattern to less than the whole plate, select the wells that have the same pattern and choose assign assay to selected wells. NOTE: the selected wells do not have to be contiguous. By holding down the shift key when clicking, ranges can be selected and by using the control key disjoint wells can be selected. Additionally the row and columns can be selected by clicking on the row and column headers.
- 3) An assay map will appear. Right click on a spot and select an assay from the popup menu of assays.
- 4) Assign assays to all spots in a well except a blank or BSA spot. Click OK when done.
- 5) The layout has now been saved.

5.1.2. How do I use an assay that is not in the list?

In the assay assignment dialog, one of the choices in the right click popup menu is to create a new assay. You can create and organize the assays as needed from this popup menu.

5.1.3. How do I assign replicates?

In the *assign* dialogs, you can choose to set replicates. Pick how many and in which direction. The table in the dialog will update if the values are entered. If you choose OK, your choices will overwrite the old data that may have been there.

5.1.4. How do I enter unusual dilution patterns?

There are several autofill options available but if these are not enough, you can hand enter the values in the table by directly editing it.

5.1.5. How do I name my samples?

The *assign* dialogs creates names by default but they can be changed in the table in the bottom of the dialog window by directly editing the fields.

5.2. Experiments

5.2.1. How do I get my data into a new experiment?

Before data is imported, a layout needs to be available. While it is possible to edit the layout after the plate has been imported it is easier to prepare the layout first.

Data is imported three ways:

- 1) Use the mouse to drag a plate listed in the Plate Data History onto the experiment. You will be prompted for the layout that you want to use with the plate.
- 2) Use the file menu to import plate. NOTE: This option is the only choice if the plate was run on different machine than the one where the analysis is being done. A standard formatted export report for the desired plate(s) needs to be available. Browse to the plate report. It will show up in the Plate Data History after it has been imported and then can be dragged into the experiment. If the report was generated using a custom format, use the Plate Data History on the instrument to re-export in the default format.
- 3) Pick a template when running a plate. On the Run dialog, there is a choice to create an experiment using a template. When the plates are finished the template is used to put the plates into an experiment. The template can be created by saving one from an existing experiment or by saving a new experiment as a template.

5.2.2. I have an experiment that I like, how do I reuse it?

The easiest way is to store the experiment as a template. Templates belong to the person who created them, so only you can edit the template. However anyone can make their own copy and then edit it to suit their own needs. The template will save the experiment with your preferences for what items are displayed without the actual data. All of these choices can then be used by creating another experiment from the template.

5.2.3. How do I share the templates/experiments with other users?

Experiments and templates are visible to everyone that uses the machine. Only you can change the items that you have created. However others can view the experiments and save a copy for themselves as desired. For a group wanting to use the same format in an experiment, creating a template that everyone uses is a great way to start.

Both experiments and templates can be exported and imported. This allows the work developing an experiment to be shared across all the instruments.

5.2.4. How do I create an experiment template?

The experiment template is created from an experiment. We create and adjust the experiment and then save it as a template. It is not necessary to save the experiment

used to generate the template, if so desired. To create an experiment template from scratch.

- 1) Create a new experiment.
- 2) Create the layout that you will want to use if it does not already exist.
- 3) Import data from the desired plate type. This data is just a placeholder so we can see what the table/plot options are doing and will not be saved in the template.
- 4) Adjust the plots and tables as desired.
- 5) Right click on the experiment (the top folder) in the experiment window.
- 6) Select “save as template.”
- 7) The experiment is not saved at this point and can be discarded if desired.

To create an experiment template from an existing experiment:

- 1) Open the experiment.
- 2) Right click on the experiment in the experiment window (top level).
- 3) Select “save as template” and create a name for the template.

5.3. Miscellaneous

5.3.1. Where did the menu item go?

Many options in the software are context sensitive so that they are not in your way when working on other parts of the software. Try making sure the item that you are going to manipulate is selected to see if the options reappear.

5.4. Data Analysis

5.4.1. How do I exclude data?

Occasionally, there are values in the imported data that we know are not correct (for experimental reasons) and need to be excluded from the data analysis. In many of the views of the data including tables, plots and the data grid, a right click brings up a popup menu that has a choice for exclude data. A few things to consider:

- 1) Data is never removed or altered. The data is simply not included in calculated values.
- 2) Excluded values are highlighted so you know they are not included in the calculations.

- 3) What is excluded depends on the context of the view. For example, if you are viewing single data points for each spot, only that spot's data is excluded. If on the other hand the view is of the well then all spots in that well are excluded.
- 4) Excluded values can be restored.

5.4.2. How do I control detection limits?

Options for the detection limits can be found by right clicking on the standard group and selecting the detection limits option.

5.4.3. How do I control the fitting?

The fitting can be adjusted for each standard. Open the experiment and right click on the standards in the plate to select "Edit Fit Profile". The dialog that appears allows you to control the type of fitting and the parameters for the fitting.

5.5. Reporting and Working with Other Applications

5.5.1. How do I print out the experiment?

The experiment can be printed to a PDF file. This in turn can be printed. A print preview can be used to make sure the content of the PDF contains the items you want to print.

5.5.2. How do I remove items from the printed output?

- 1) Right click the experiment in the tree view of the experiment and select "Report Selection Mode."
- 2) Select the "show report selection" option.
- 3) The tree now has check boxes next to each item. Items that are checked are included in the report.
- 4) Deselect "show report selection"

5.5.3. Can I get a text file containing the data?

Yes...

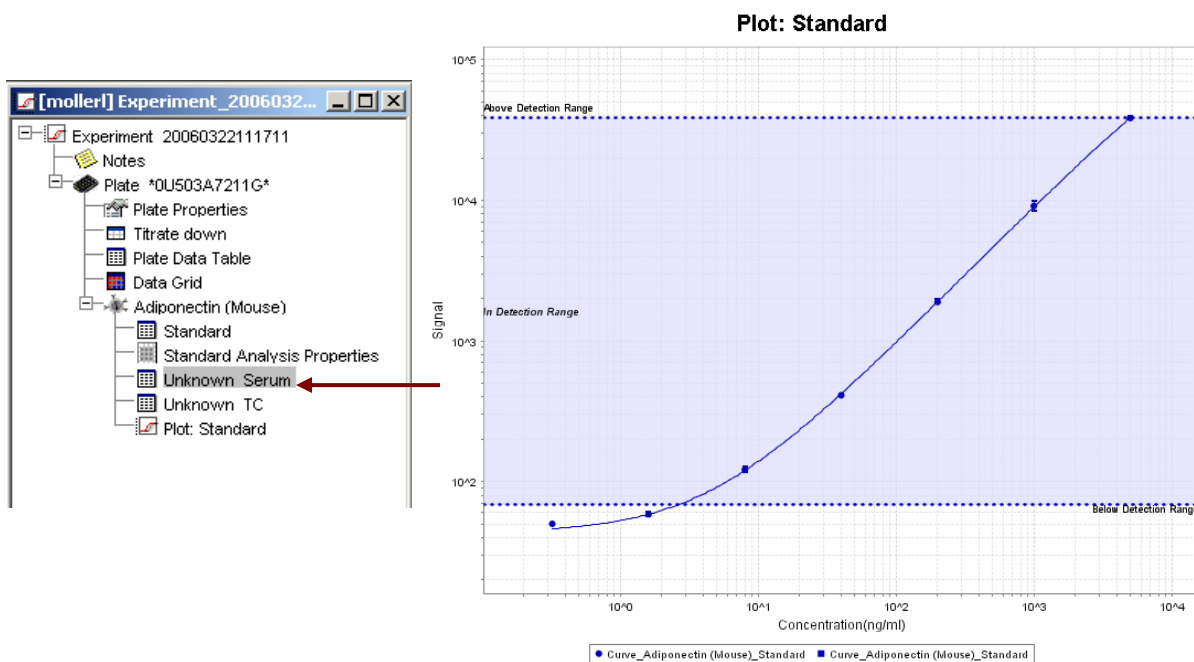
- 1) Go to the experiment tree.
- 2) Right click on the experiment to see "export data to text file"
- 3) The formatting options are the same as those for the ECL data exports.
- 4) Pick which piece of data you wish to export from the pull down box.
- 5) Click on the Export button.

5.5.4. How do I get a plot/table into MS Word or MS Powerpoint?

The software supports cut and paste of many items into other applications using standard clipboard operations. Right click the plot or table and select *copy plot / table* or go to *EDIT > Copy plot / table* then paste in the application of choice.

6. Appendix - Fitting functions

The Plot: Standard window displays the standard curve for an analyte within an experiment. To open a plot, double click on the Plot icon within the Experiment tree menu.



6.1. Four Parameter Logistic Fit

The default fitting function is a weighted four parameter logistic fit (4PL) which best describes the standard curve for most assays run on MSD instruments. Compared to linear regression, four parameter logistic regression provides accurate quantitation over a wider range of analyte concentrations

The minimum number of non-zero calibrators for 4PL regression is five however it is recommended that at least eight standard concentrations, including a zero concentration standard be used to generate the standard curve. Recommended standard concentrations are listed in the protocol insert of kitted assays that are provided with calibrators.

The 4PL equation is:

$$y = b_2 + \frac{b_1 - b_2}{1 + (x / b_3)^{b_4}}$$

Where:

y = ECL counts (response)

x = concentration

b1 = minimum response plateau

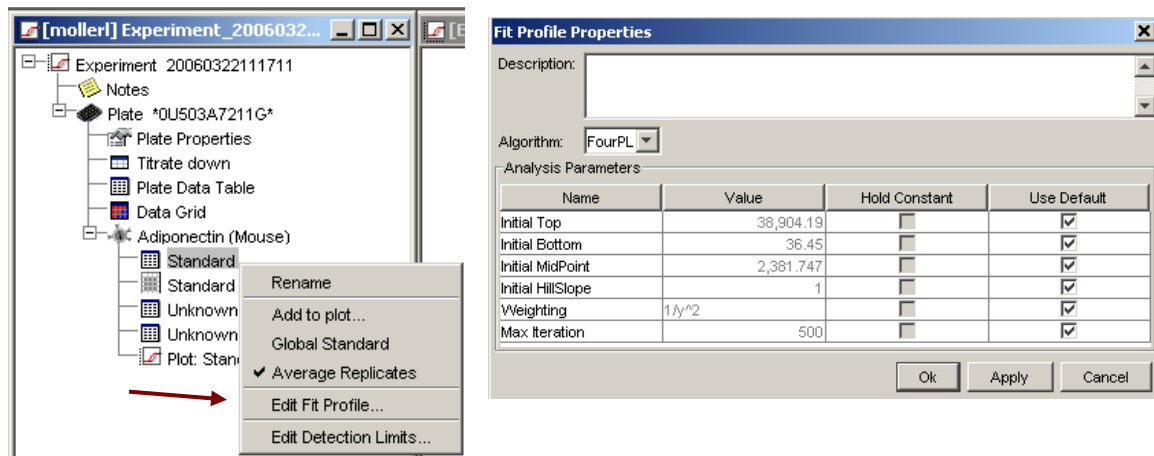
b2 = maximum response plateau

b3 = the concentration at which 50% of the maximal response is observed

b4 = the "slope" or shape parameter

Curve fitting is an iterative process that optimizes parameter values by minimizing the sum of the squared residuals. Residuals are the differences between the experimental data and the fitted curve at each point. The curve fitting algorithm uses a weighting function ($1/y^2$) to compensate for the differences in magnitude of the residuals at low and high analyte concentrations. The residuals are typically lower for standards with the smallest responses and higher for standards with the largest responses. By weighting this way, curve fitting is not biased by small or large residuals.

The default initial values for parameters can be modified by right clicking the **Standard table** in the Experiment tree and selecting **Edit Fit Profile**. Initial values can be changed in the **Value** column after un-checking a parameter in the **Use Default** column.

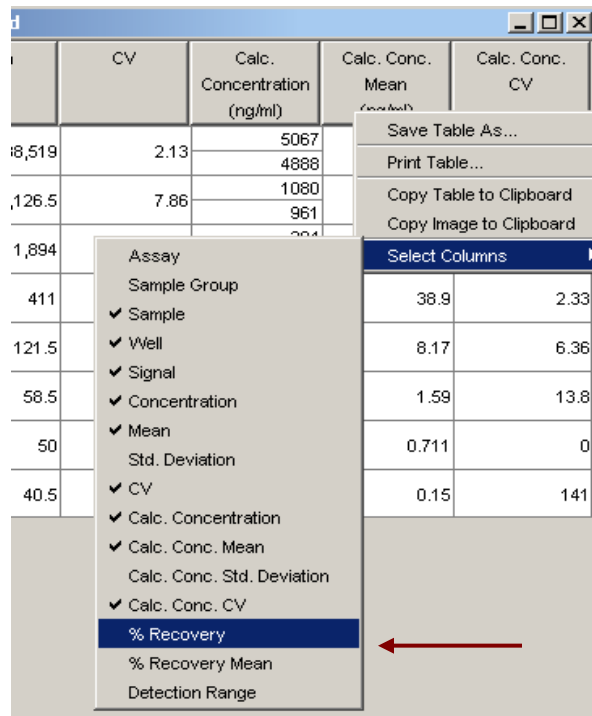


For competitive assays, it is often necessary to try the fitting with and without the weighting function. The weighting may reduce the sensitivity of the assay by fitting the higher calibrators (lower signals) better in the competitive assay.

Evaluating the quality of the fit

A practical means of evaluating the quality of the fit is to compare the input concentration of the standards to the calculated concentration obtained by backfitting to the calibration curve using the formula: $\% \text{ Recovery} = \text{Calculated Concentration} / \text{Expected Concentration} \times 100$. If a curve describes the experimental data well the $\% \text{ Recovery}$ values should fall between 80 to 120 $\%$ of the expected concentration.

To display $\% \text{ recovery}$ in the Standards table, open the table and right click on the grey Table header and select **$\% \text{ Recovery}$** or **$\% \text{ Recovery Mean}$** from the list of available columns in the **Select Columns** menu.



	CV	Calc. Concentration (ng/ml)	Calc. Conc. Mean	Calc. Conc. CV
18,519	2.13	5067		
		4888		
,126.5	7.86	1080		
		961		
1,894		204		
411			38.9	2.33
121.5			8.17	6.36
58.5			1.59	13.8
50			0.711	0
40.5			0.15	141

- Assay
- Sample Group
- ✓ Sample
- ✓ Well
- ✓ Signal
- ✓ Concentration
- ✓ Mean
- Std. Deviation
- ✓ CV
- ✓ Calc. Concentration
- ✓ Calc. Conc. Mean
- Calc. Conc. Std. Deviation
- ✓ Calc. Conc. CV
- % Recovery**
- % Recovery Mean
- Detection Range

7. Technical Support

Should you encounter an error, please document:

- The Error Code and text in the Error Dialog
- How you generated the error
- Is the error reproducible?
- Get a copy of all of the files in the following folder:

C:\Program Files\MSD Discovery Workbench\log

Email this information to CustomerSupport@mesoscale.com.

Please send comments or feedback on the software, including feature requests, or design change requests, to CustomerSupport@mesoscale.com.