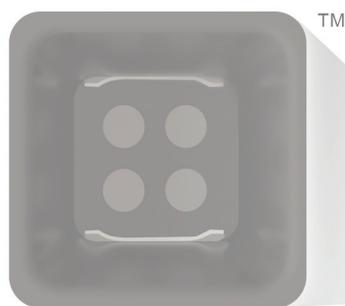


# MSD<sup>®</sup> MULTI-SPOT Assay System

## Human Cytokine Assays Tissue Culture Kit (384-well)

**T**-PLEX<sup>®</sup>



# MSD Cytokine Assays

## Human Cytokine Assays

### Tissue Culture Kit (384-well)

*This package insert must be read in its entirety before using this product.*

**FOR RESEARCH USE ONLY.**

**NOT FOR USE IN DIAGNOSTIC PROCEDURES.**

## MESO SCALE DISCOVERY®

A division of Meso Scale Diagnostics, LLC.

1601 Research Boulevard

Rockville, MD 20850-3173 USA

[www.mesoscale.com](http://www.mesoscale.com)

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, mesoscale.com, www.mesoscale.com, methodicalmind.com, www.methodicalmind.com, DISCOVERY WORKBENCH, InstrumentLink, MESO, MesoAccess, MesoAdvantage, MesoSphere, Methodical Mind, Methodical Mind Enterprise, MSD GOLD, MULTI-ARRAY, MULTI-SPOT, QuickPlex, Parsec, ProductLink, SECTOR, SECTOR PR, SECTOR HTS, SULFO-TAG, TeamLink, TrueSensitivity, TURBO-BOOST, TURBO-TAG, N-PLEX, R-PLEX, S-PLEX, T-PLEX, U-PLEX, V-PLEX, MSD (design), MSD (luminous design), Methodical Mind (design), Parsec (design), 96 WELL SMALL-SPOT (design), 96 WELL 1-, 4-, 7-, 9-, & 10-SPOT (designs), 384 WELL 1- & 4-SPOT (designs), N-PLEX (design), R-PLEX (design), S-PLEX (design), T-PLEX (design), U-PLEX (design), V-PLEX (design), It's All About U, SPOT THE DIFFERENCE, The Biomarker Company, and The Methodical Mind Experience are trademarks and/or service marks owned by or licensed to Meso Scale Diagnostics, LLC. All other trademarks and service marks are the property of their respective owners.  
©2009-2017, 2020 Meso Scale Diagnostics, LLC. All rights reserved.

# Table of Contents

Introduction .....	4
Principle of the Assay .....	4
Reagents Supplied.....	5
Additional Materials and Equipment.....	6
Safety .....	6
Sample Preparation .....	7
Reagent Preparation .....	7
Assay Protocol .....	10
Assay Notes.....	11
Specificity.....	12
Summary Protocol .....	13
Appendix .....	14
Plate Diagram .....	15

## Contact Information

### MSD Customer Service

Phone: 1-240-314-2795  
Fax: 1-301-990-2776  
Email: [CustomerService@mesoscale.com](mailto:CustomerService@mesoscale.com)

### MSD Scientific Support

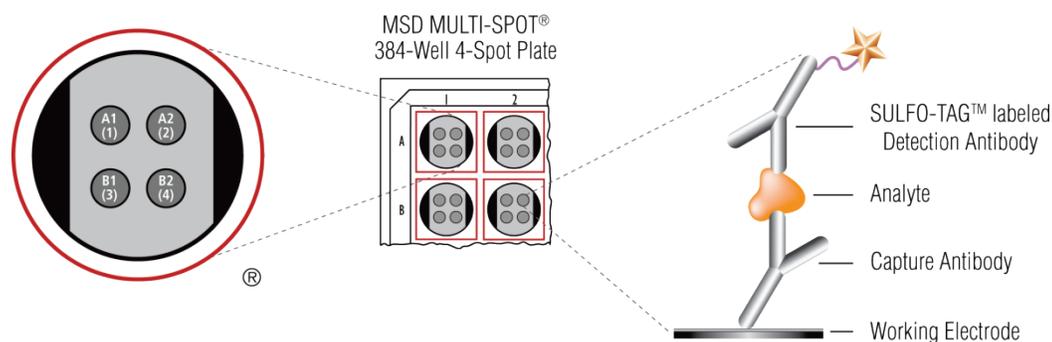
Phone: 1-240-314-2798  
Fax: 1-240-632-2219 attn: Scientific Support  
Email: [ScientificSupport@mesoscale.com](mailto:ScientificSupport@mesoscale.com)

# Introduction

MSD offers a broad range of human cytokine assays in 1-Spot MULTI-ARRAY® and 4-Spot MULTI-SPOT® 384-well plate formats. This product insert outlines an assay protocol recommended for tissue culture samples. This insert also describes ways the user can modify these protocols to meet specific workflow or performance requirements.

## Principle of the Assay

MSD 384-well Cytokine Assays measure one to four cytokines in an MSD 384-well MULTI-ARRAY or MULTI-SPOT plate. The assays employ a sandwich immunoassay (Figure 1). MSD provides a plate precoated with capture antibodies. The user adds the sample and a solution containing detection antibodies conjugated with electrochemiluminescent (MSD SULFO-TAG™) labels throughout one or more incubation periods. Analytes in the sample bind to capture antibodies immobilized on the working electrode surface (Figure 2); recruitment of the detection antibodies by the bound analytes completes the sandwich. The user adds an MSD buffer that provides the appropriate chemical environment for electrochemiluminescence (ECL) and loads the plate into an MSD® instrument where a voltage applied to the plate electrodes causes the captured labels to emit light. The instrument measures the intensity of emitted light to provide a quantitative measure of analytes in the sample.



**Figure 1.** A cytokine capture antibody is precoated onto specific spots of a 4-Spot MSD MULTI-SPOT plate. Calibrator solutions or samples are incubated in the MULTI-SPOT plate, and each cytokine binds to its corresponding capture antibody spot. Cytokine levels are quantitated using a cytokine-specific detection antibody conjugated with MSD SULFO-TAG reagent to stimulate the production of ECL when voltage is applied to the plate electrodes.

# Reagents Supplied

Necessary components (Table 1) are supplied with the Human Cytokine Assay Tissue Culture Kit (96-well) kit.

**Table 1.** Components supplied

Product Description	Storage
<b>MULTI-SPOT or MULTI-ARRAY Plate</b> 384-well MSD plate spotted with specific anticytokine capture antibodies	2–8 °C
<b>Detection Antibody Blend</b> Anticytokine antibodies labeled with MSD SULFO-TAG reagent and premixed to provide a 50X stock solution of each antibody	2–8 °C
<b>Cytokine Calibrators</b> <b>Human Tissue Culture Calibrator 2:</b> Provided for singleplex or multiplex combinations including GM-CSF, IL-1 $\beta$ , IL-2, IL-4, IL-5, IL-6, IL-8, IL-10, IL-12p70, IL-13, and TNF- $\alpha$ assays.* <b>Human IFN-<math>\gamma</math> Calibrator:</b> Provided for singleplex or multiplex assays including IFN- $\gamma$ assay.* For the lot-specific concentration of Calibrators mentioned above, please refer to the Certificate of Analysis (COA) supplied with the kit. The COAs are also available at <a href="http://www.mesoscale.com">www.mesoscale.com</a> . <b>Other Calibrators:</b> For all other cytokine assays, individual stock Calibrators are provided at 1 $\mu$ g/mL concentration.	$\leq$ -70 °C
<b>Diluent 1</b> RPMI-based medium for dilution of Calibrators	2–8 °C
<b>Diluent 100</b> Contains blocking and stabilizing agents	2–8 °C
<b>Blocker B</b> (included with multiplexed assays containing IFN- $\gamma$ or IL-12p40) Contains blocking and stabilizing agents	RT
<b>Read Buffer T (4X)</b> 4X Read Buffer T with surfactant	RT

\*For the lot-specific concentration of these Calibrators, please refer to the COA supplied with the kit. The COAs are also available at [www.mesoscale.com](http://www.mesoscale.com).

RT = room temperature

# Additional Materials and Equipment

- Deionized water for diluting concentrated buffers
- 50-mL tubes for reagent preparation
- 15-mL tubes for reagent preparation
- Microcentrifuge tubes for preparing serial dilutions
- Phosphate-buffered saline (PBS) plus 0.05% Tween-20 for plate washing or MSD Wash Buffer (Catalog No. R61AA-1)
- Appropriate liquid handling equipment for the desired throughput, capable of accurately dispensing 10  $\mu$ L to 40  $\mu$ L into a 384-well microtiter plate
- Plate washing equipment: automated plate washer or multichannel pipette
- Adhesive plate seals
- Microtiter plate shaker

## Safety

Use safe laboratory practices and wear gloves, safety glasses, and lab coats when handling kit components. Handle and dispose of all hazardous samples properly in accordance with local, state, and federal guidelines.

Additional safety information is available in the Product Safety Data Sheet (SDS), which can be obtained from MSD Customer Service or at [www.mesoscale.com](http://www.mesoscale.com).

# Sample Preparation

This section provides a general guide for the preparation of tissue culture supernatant for use in tissue culture assays. Safe laboratory practices and personal protective equipment such as gloves, lab coat, and safety glasses should be used at all times when handling samples. All samples of a potentially infectious or hazardous origin should be handled in the manner outlined by the Centers for Disease Control and the Occupational Safety and Health Administration for blood-borne pathogens and human- or animal-sourced materials. When analyzing some samples, it may be necessary to dilute the sample by a factor of 2 to 10 to achieve the most accurate quantitation in the MSD Human Cytokine Assays. Please contact MSD Customer Service with any questions.

## Tissue Culture Supernatant

Most tissue culture supernatant samples generally do not require any dilution before being used in the MSD Human Cytokine Assays. Samples from experimental conditions with extremely high levels of cytokines may require dilution.

# Reagent Preparation

Bring all reagents to room temperature. Thaw the stock calibrator on ice.

## Prepare Blocker B Solution

Blocker B should be made fresh prior to each run. Blocker B will appear milky and not completely in solution, but can still be used. Blocker B does not require filtration.

For multiplexed assays containing IFN- $\gamma$ :

Prepare a 1% (w/v) solution of Blocker B in PBS (20 mL per plate) by adding 200 mg Blocker B to 20 mL PBS. Vortex to mix and use.

For multiplexed assays containing IL-12p40:

Prepare a 0.1% (w/v) solution of Blocker B in PBS (20 mL per plate) by adding 20 mg Blocker B to 20 mL PBS. Vortex to mix and use.

## Prepare Calibrator and Control Solutions

Dilute Calibrators in Diluent 1: The calibration curve preparation instructions listed below will generate a standard curve ranging from 10,000 pg/mL to 2.4 pg/mL. The curve should be adjusted as necessary to provide the proper range for test samples.

Diluent 1 is a standard tissue culture growth medium with 10% serum. If tissue culture samples are in a different medium, use that medium for calibration curve preparation. However, please note that if using a serum-free medium, the presence of some carrier protein in solution is necessary to prevent loss of analyte to the labware.

## Prepare Calibration Curve

The Human Tissue Culture Calibrator 2 should be used for singleplex and multiplex assays that include the following assays: GM-CSF, IL-1 $\beta$ , IL-2, IL-4, IL-5, IL-6, IL-8, IL-10, IL-12p70, IL-13, and TNF- $\alpha$ .

The Human IFN- $\gamma$  Calibrator should be used for singleplex and multiplex assays that include IFN- $\gamma$ .

**Note:** For the calibrators mentioned above, please refer to the COA supplied with each kit for the measured concentration values and top of the curve concentrations. Alternatively, continue to use the nominal top of curve concentration value of 10,000 pg/mL in the protocol and data analysis for convenience.

For the remaining cytokine assays, individual stock Calibrators are provided at 1 µg/mL concentration, and 10,000 pg/mL should be used as the nominal top of curve concentration value.

The following protocol (see Figure 2) should be used to prepare the calibration curve.

CAL-1: Add 10 µL of each calibrator supplied to Diluent 1 such that the final volume is 1,000 µL. Use this high Calibrator (nominal concentration: 10,000 pg/mL) to prepare the standard curve following a 1:4 dilution series.

CAL-2: Add 50 µL of 10,000 pg/mL combined high Calibrator to 150 µL of Diluent 1.

CAL-3: Add 50 µL of 2,500 pg/mL Calibrator to 150 µL of Diluent 1.

CAL-4: Add 50 µL of 625 pg/mL Calibrator to 150 µL of Diluent 1.

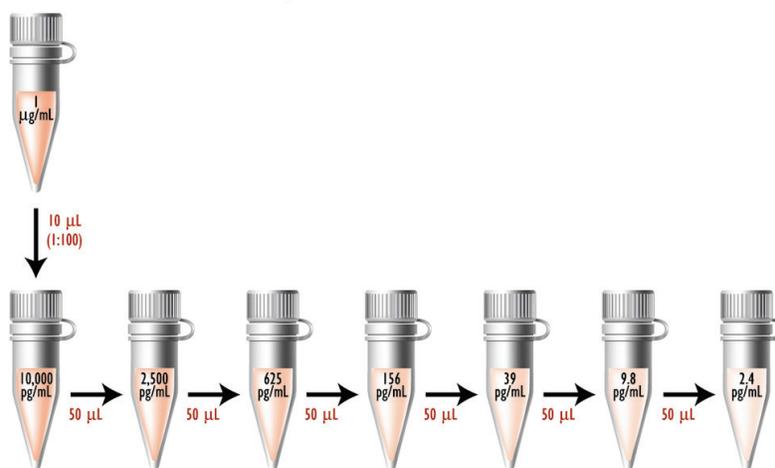
CAL-5: Add 50 µL of 156 pg/mL Calibrator to 150 µL of Diluent 1.

CAL-6: Add 50 µL of 39 pg/mL Calibrator to 150 µL of Diluent 1.

CAL-7: Add 50 µL of 9.8 pg/mL Calibrator to 150 µL of Diluent 1.

CAL-8: 150 µL of Diluent 1 (zero Calibrator blank).

Alternatively, Calibrators can be prepared in the sample matrix or diluent of choice to verify acceptable performance in these matrices. In general, the presence of some protein in the sample matrix helps prevent loss of analyte by adsorption onto the sides of tubes, pipette tips, and other surfaces.



**Figure 2.** Calibration curve preparation from 1 µg/mL stock solution.

We anticipate that most customers will continue to use the nominal (10,000 pg/mL) values in their protocol and data analysis of GM-CSF, IFN-γ, IL-1β, IL-2, IL-4, IL-5, IL-6, IL-8, IL-10, IL-12p70, IL-13, and TNF-α assays. Many applications do not require absolute quantitation of analytes relative to a fixed standard, and the nominal values allow convenient use and consistent data analysis. Dose-response studies, IC-50 measurements, and experiments where the relative abundance of the analyte is sufficient are examples of situations where using nominal values for the Calibrators is consistent with good practice.

Applications that require accurate, absolute quantitation of analytes relative to a fixed standard will benefit from using the measured value for the Calibrators. There will be no change in the protocol for the preparation of Calibrator solutions, but users will need to enter the measured value of the Calibrators into their data analysis template. In MSD's DISCOVERY WORKBENCH® software, users will simply enter the concentration for the Calibrators at the top of the curve in the plate layout. Each lot of kits will have specific measured values.

## Prepare Detection Antibody Solution

**Note:** Detection antibody solutions should be kept in the dark because some antibodies may be light sensitive.

MSD provides detection antibodies as a 50X stock solutions (some vials may be labeled 50 µg/mL). The working detection antibody solution is 1X (or 1.0 µg/mL).

For one plate, combine:

- 160 µL of 50X Detection Antibody Blend
- 7.84 mL of Diluent 100

## Prepare Wash Buffer

MSD provides Wash Buffer as a 20X stock solution. The working solution is 1X. PBS + 0.05% Tween-20 can be used instead.

For one plate, combine:

- 30 mL of MSD Wash Buffer (20X)
- 570 mL of deionized water

## Prepare Read Buffer T

MSD provides Read Buffer T as a 4X stock solution. The working solution is 2X.

For one plate, combine:

- 8.5 mL of Read Buffer T (4X)
- 8.5 mL of deionized water

Diluted read buffer can be prepared in advance and stored at room temperature in a tightly sealed container for up to one month.

## Prepare MSD Plate

MSD plates are precoated with capture antibodies (Figure 1) and exposed to a proprietary stabilizing treatment to ensure the integrity and stability of the immobilized antibodies. Plates can be used as delivered; no additional preparation (e.g., prewetting) is required. The plate has not been preblocked, and a discrete blocking step is not generally required for MSD cytokine assays. For some cytokines, however, a blocking step may improve the assay sensitivity.

# Assay Protocol

A suggested plate layout for calibrating Cytokine Assay Kits is shown in Figure 3.

For MULTI-SPOT Assays containing IFN- $\gamma$ , a blocking step is required before beginning the assay to achieve the best performance. Dispense 20  $\mu$ L of the 1% (w/v) Blocker B Solution into each well of the MSD plate. Seal the plate with an adhesive plate seal and incubate with vigorous shaking for 1 hour at room temperature. Wash the plate three times with at least 90  $\mu$ L/well of 1X MSD Wash Buffer or PBS + 0.05% Tween-20 and proceed with the assay protocol at Step 1.

For MULTI-SPOT Assays containing IL-12p40, a blocking step is required before beginning the assay to achieve the best performance. Dispense 20  $\mu$ L of the 0.1% (w/v) Blocker B Solution into each well of the MSD plate. Seal the plate with an adhesive plate seal, and incubate with vigorous shaking for 1 hour at room temperature. Wash the plate three times with at least 90  $\mu$ L/well of 1X MSD Wash Buffer or PBS + 0.05% Tween-20, and proceed with the assay protocol at Step 1.

	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	
A	10,000 pg/mL																								
B																									
C	2500 pg/mL																								
D																									
E	625 pg/mL																								
F																									
G	2.4 pg/mL																								
H																									
I	39 pg/mL																								
J																									
K	9.8 pg/mL																								
L																									
M	2.4 pg/mL																								
N																									
O	0 pg/mL																								
P																									

**Figure 3.** Suggested plate setup for calibrating MSD Cytokine Assay Kits.

In Figure 3, columns 1 and 2 provide a standard curve in duplicate. The remaining wells are available for samples. Some assays may require lower limits of detection, and in these cases, the 625 pg/mL point can be removed and a 0.6 pg/mL point added between 2.4 pg/mL and 0 pg/mL to provide an additional lower calibration point. An alternative setup that contains a larger dynamic range for the standard curve can be prepared by running Calibrators in a 12-point titration in duplicate across the top or bottom of the plate. The concentrations of Calibrators run may be adjusted depending on the desired dynamic range for the experiment. The data from this calibration curve can be analyzed using any standard data analysis package.

## STEP 1: Add Sample or Calibrator Standard

- Dispense 10  $\mu$ L of each Calibrator or Sample Solution into a separate well of the MSD plate. Figure 3 illustrates one plate arrangement of Calibrator solutions that can be used to evaluate the performance of the assay.
- Seal the plate with an adhesive plate seal, and incubate for 1 to 2 hours with vigorous shaking (500–1,000 rpm) at room temperature.

**Note:** In general, shaking the plate results in better mixing and more rapid binding of the sample to the capture antibodies. If a protocol without shaking is preferred, a longer incubation time (4 hours or longer) may be required to achieve the same sensitivity.

## STEP 2: Add Detection Antibody Solution

- ❑ Dispense 10  $\mu$ L of the 1X detection antibody solution into each well of the MSD plate.
- ❑ Seal the plate and incubate for 1 to 2 hours with vigorous shaking at room temperature. If a protocol without shaking is preferred, a longer incubation time may be required to achieve the same sensitivity.

## STEP 3: Wash and Read

- ❑ Wash the plate three times with at least 90  $\mu$ L/well of 1X MSD Wash Buffer or PBS + 0.05% Tween-20.
- ❑ Add 35  $\mu$ L of 2X Read Buffer T to each well of the MSD plate. Analyze the plate on the MSD instrument. Incubation in Read Buffer T is not required before reading the plate.

**Note:** Bubbles in the fluid will interfere with a reliable reading of the MULTI-SPOT plate. Use reverse pipetting (Appendix) techniques to ensure bubbles are not created when dispensing the Read Buffer T.

# Assay Notes

- ❑ Unwashed Assay: The protocol can be converted to an unwashed assay by eliminating the wash step following the detection antibody incubation, adding 15  $\mu$ L 2X Read Buffer T, and analyzing the plate on the MSD instrument. This method may result in a loss of sensitivity, the effect of which will vary for each cytokine.
- ❑ Sample Matrices: In general, these plates have been found to work well to measure cytokine levels in cell culture supernatants and simple buffers. Sample dilution and/or spike recovery studies in the sample matrix of interest should be carried out to verify acceptable performance in the matrix. Use the Human Cytokine Ultra-Sensitive Assays or V-PLEX<sup>®</sup> Human Cytokine Assays available on [www.mesoscale.com](http://www.mesoscale.com) for testing serum and plasma samples.
- ❑ Combining the Sample and Detection Antibody Addition Steps: The protocols described above call for incubating the sample in the wells before the addition of the detection antibody (sequential incubations). This procedure is used because in some selected assays using a polyclonal detection, the detection antibody itself may include antibodies that compete with the immobilized capture antibody. In some assays, however, MSD has found that the sample and detection antibodies may be added concurrently (simultaneous incubation) with little or no loss in performance.
- ❑ Stability of Assay in Read Buffer T: The plates do not need to be read immediately after the addition of Read Buffer T. In the washed assay, the loss of signal is typically less than 20% over a 1-hour incubation in Read Buffer T, and it is usually stabilized by 2 hours. The observed loss is due to the reestablishment of equilibrium in the well.
- ❑ Optimizing Assay Sensitivity: The wash protocols described above were developed to provide excellent assay performance with a minimum number of steps, minimal sample volume requirement, and a rapid assay completion time. One or more of the following assay modifications may be used to further improve the assay sensitivity.
  - Increasing incubation times: The incubation time with the detection antibody can be increased to 4 hours with shaking (or at least 12 hours without shaking).
  - Increasing sample volumes up to 20  $\mu$ : Efficient shaking and/or longer incubation times become more important at higher volumes. If a larger sample volume is used, the concentration of the detection antibody must be adjusted in the detection antibody solution to compensate for a change in volume.
  - Washing sample from the well before the addition of the detection antibody mix: This may be especially useful for large sample volumes to avoid dilution of the detection antibodies. Not all assays will benefit from the addition of this step, and some assays may experience decreased sensitivity.

**Note:** See the Appendix for additional notes.

# Specificity

The capture and detection antibody pairs used in MSD cytokine assays have been selected by an optimization process that is designed to minimize cross-reactivity with other cytokine assays. In addition, the chosen antibodies are specific for the particular human cytokine of interest and show minimal cross-reactivity with other species.

# Summary Protocol

## MSD 384-Well MULTI-ARRAY and MULTI-SPOT Human Cytokine Assays: Tissue Culture Kit

MSD provides this summary protocol for convenience.

Please read the entire detailed protocol before performing the MSD Human Cytokine Assay.

### Sample and Reagent Preparation

- Prepare Blocker B Solution.
  - For MULTI-SPOT Assays with IFN- $\gamma$ , prepare a solution of 1% Blocker B in PBS.
  - For MULTI-SPOT Assays with IL-12p40, prepare a solution of 0.1% Blocker B in PBS.
- Prepare Calibrator solutions and calibration curves.
  - Use the 1  $\mu\text{g/mL}$  Calibrator stock to prepare an 8-point calibration curve of 10,000, 2,500, 625, 156, 39, 9.8, 2.4, and 0  $\text{pg/mL}$ .
  - The calibration curve can be modified as necessary to meet specific assay requirements.
- Prepare detection antibody solution by diluting stock detection antibody to 1X (or 1  $\mu\text{g/mL}$ ) in 8.0 mL (per plate) of Diluent 100.
- Prepare 17 mL of 2X Read Buffer T by diluting 4X Read Buffer T 2-fold with deionized water.
- Prepare 1X MSD Wash Buffer by diluting the 20X stock solution to 1X with deionized water. Use at least 90  $\mu\text{L/well}$  for plate washing.

### STEP 1: For MULTI-SPOT Assays with IFN- $\gamma$ or IL-12p40 only:

- Dispense 20  $\mu\text{L/well}$  1% Blocker B Solution for IFN- $\gamma$  or 0.1% Blocker B Solution for IL-12p40.
- Incubate at room temperature with vigorous shaking (500–1,000 rpm) for 1 hour.
- Wash the plate three times with at least 90  $\mu\text{L/well}$  of 1X MSD Wash Buffer or PBS+ 0.05% Tween-20.
- Proceed with sample/Calibrator addition.

### STEP 2: Add Sample or Calibrator

- Dispense 10  $\mu\text{L/well}$  Calibrator or sample.
- Incubate at room temperature with vigorous shaking for 1 to 2 hours.

### STEP 3: Add Detection Antibody Solution

- Add 10  $\mu\text{L/well}$  of 1X detection antibody solution.
- Incubate at room temperature with vigorous shaking for 1 to 2 hours.

### STEP 4: Wash and Read Plate

- Wash the plate three times with at least 90  $\mu\text{L/well}$  of 1X MSD Wash Buffer or PBS + 0.05% Tween-20.
- Add 35  $\mu\text{L/well}$  of 2X Read Buffer T.
- Analyze plate on the MSD instrument.

# Appendix

## Background signal and negative signals

The output signal produced by electrochemiluminescence assays is in units of counts of light measured by a charge-coupled device (CCD) camera or photodiode. As with any measurement technique, there is a certain amount of normal variation in the signal (instrument noise) which sets the threshold for the lowest levels of signal that can be measured (noise floor). This variation is different depending upon the size of the working electrode with typical values of about 10 counts for 384-well small spot and 50 counts for 384-well 4-spot plates. When the background signal of an assay approaches the noise floor (i.e., the mean signal of negative controls or sample blanks is close to zero), it is possible to observe negative counts for some wells.

## Signal Levels

The camera system is linear over nearly a 6-log dynamic range. The highest achievable signals on the MSD instruments are between 1.0 and 2.0 million counts. If the signals from the highest point on the calibration curve are not approaching 1.0 million counts, the high end of the calibration curve may be extended. The lowest observed signals using Read Buffer T (2X) are between 10 and 50. Negative signal values may occur due to instrument noise, omission or usage of the incorrect read buffer, or incorrect amount of detection antibody.

## Curve Fitting

MSD DISCOVERY WORKBENCH software uses least-squares fitting algorithms to generate the standard curve that will be used to calculate the concentration of analytes in the samples. The assays have a wide dynamic range (3–4 logs) that allows accurate quantification without the need for dilution in many cases. By default, the software uses a 4-parameter logistic model (or sigmoidal dose-response) and includes a  $1/Y^2$  weighting function. The weighting function is important because it provides a better fit of data over a wide dynamic range, particularly at the low end of the standard curve.

## Reverse Pipetting

Most manual hand pipets have two plunger positions for pipetting liquids. The first position is calibrated to allow aspiration and dispensing of user-specified amounts of liquid and the second (blowout) position enables the user to expel any residual liquid after the pipet has been pushed to the first position. When a pipet is used to dispense liquid by moving the plunger to the first position followed by the second (blowout) position, bubbles may be created in the dispensed liquid. The reverse pipetting technique is designed to allow precise pipetting while avoiding the creation of bubbles. The technique is to push the pipet plunger past the first position to the second position before aspirating liquid into the tip, thereby aspirating slightly more liquid than the desired volume (overdraw). To dispense the liquid from the tip, the pipet plunger is pushed to the first position only. This allows precise dispensing without the introduction of bubbles. When using the reverse pipetting technique, it is important not to overdraw excess liquid into the pipetting instrument.

# Plate Diagram

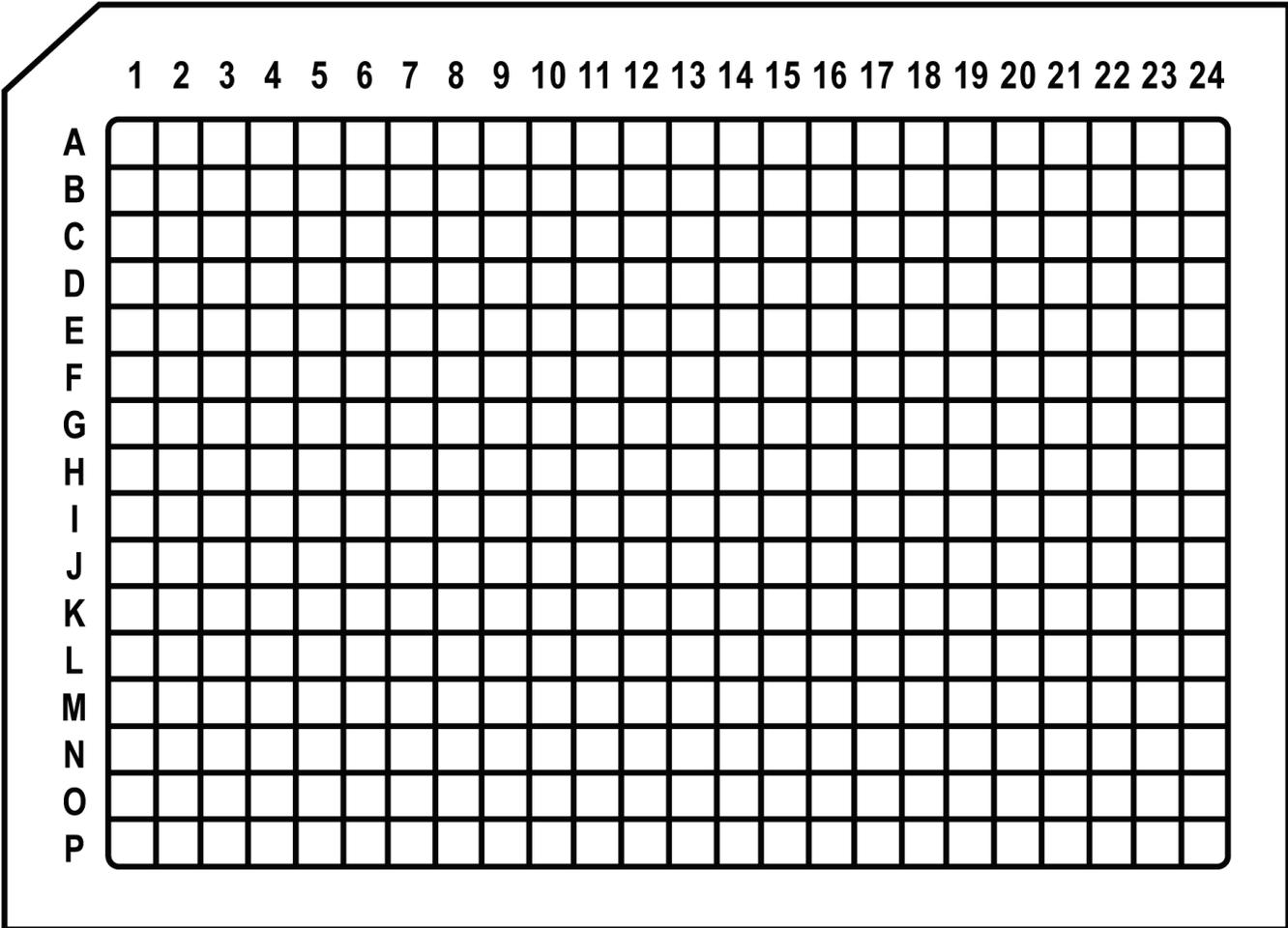


Figure 4. Plate diagram. A similar plate layout can be created in Excel and easily imported into DISCOVERY WORKBENCH software.