

U-PLEX[®] Biomarker Group 1 (NHP)

Singleplex Assays



MSD U-PLEX Platform

U-PLEX Biomarker Group 1 (NHP) Singleplex Assays

For use with serum, EDTA plasma, and cell culture supernatants.

Catalog numbers of U-PLEX Biomarker Group 1 (NHP) Singleplex Assays are provided in Table 8 on page 15.

This package insert should 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 Blvd.

Rockville, MD 20850 USA

www.mesoscale.com

MESO SCALE DISCOVERY, Meso Scale Diagnostics, MSD, mesoscale.com, www.mesoscale.com, methodicalmind.com, www.methodicalmind.com, DISCOVERY WORKBENCH, InstrumentLink, MESO, MesoSphere, Methodical Mind, MSD GOLD, MULTI-ARRAY, MULTI-SPOT, QuickPlex, ProductLink, SECTOR, SECTOR HTS, SECTOR PR, 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 (head logo), 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.

©2015-2018, 2020-2021, 2023 Meso Scale Diagnostics, LLC. All rights reserved.

Table of Contents

Introduction	4
Principle of the Assay	5
Components.....	6
Instrument Compatibility.....	8
Additional Materials and Equipment	8
Safety	8
Best Practices	9
Reagent Preparation	10
Assay Protocols.....	13
Assay Performance	14
Appendix	15
Summary Protocols.....	18
Plate Diagrams.....	20

Contact Information

MSD Customer Service

Phone:1-240-314-2795

Fax: 1-301-990-2776

Email: CustomerService@mesoscale.com

MSD Scientific Support

Phone:1-240-314-2798

Fax: 1-240-632-2219 Attn: Scientific Support

Email: ScientificSupport@mesoscale.com

Introduction

The MSD U-PLEX platform combines high sensitivity and a rapid read time (less than 2 minutes) with the flexibility to easily design and build custom assays and efficiently transition from singleplex to multiplex assays. U-PLEX Singleplex assays have high sensitivity, provide up to 5 logs of linear dynamic range, and use minimal sample volume.

The U-PLEX Biomarker Group 1 (NHP) contains 59 analytes (Table 1) that are important in many biological processes.

A representative data set for each of the assays in U-PLEX Biomarker Group 1 (NHP) is presented in the product-specific datasheets available at www.mesoscale.com/support/datasheets.

Table 1: Assays included in U-PLEX Biomarker Group 1 (NHP)

Assays			
CTACK (CCL27)	IL-1RA	IL-17A/F	MIF
Eotaxin (CCL11)	IL-2	IL-17B	MIP-1 α (CCL3)
Eotaxin-2 (CCL24)	IL-4	IL-17C	MIP-1 β (CCL4)
Eotaxin-3 (CCL26)	IL-5	IL-17D	MIP-3 α (CCL20)
ENA-78 (CXCL5)	IL-6	IL-17F	MIP-3 β (CCL19)
FLT3L	IL-7	IL-18	MIP-5 (CCL15)
Fractalkine (CX3CL1)	IL-8 (CXCL8)	IL-22	SDF-1 α (CXCL12)
G-CSF (CSF2)	IL-9	IL-23	TARC (CCL17)
GM-CSF (CSF3)	IL-10	IP-10 (CXCL10)	TNF- α
GRO- α (CXCL1)	IL-12/IL-23p40	I-TAC (CXCL11)	TNF- β
I-309 (CCL1)	IL-12p70	MCP-1 (CCL2)	TPO
IFN- α 2a	IL-13	MCP-2 (CCL8)	TRAIL
IFN- γ	IL-15	MCP-4 (CCL13)	VEGF-A
IL-1 α	IL-16	M-CSF (CSF1)	YKL-40
IL-1 β	IL-17A	MDC (CCL22)	

Principle of the Assay

Singleplex assays are supplied on MSD GOLD™ Small Spot Streptavidin 96-well or MSD Streptavidin 384-well plates (Figure 1). These plates provide high sensitivity, consistent performance, and excellent inter- and intralot uniformity.

Each singleplex assay is supplied with a biotinylated capture antibody that binds to streptavidin on the plate surface. Analytes in the sample bind to the capture reagents. Detection antibodies conjugated with electrochemiluminescent labels (MSD GOLD SULFO-TAG™) bind to the analytes to complete the sandwich immunoassay. Once the immunoassay is complete, the plate is loaded 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 (which is proportional to the amount of analyte present in the sample) and provides a quantitative measure of each analyte in the sample.

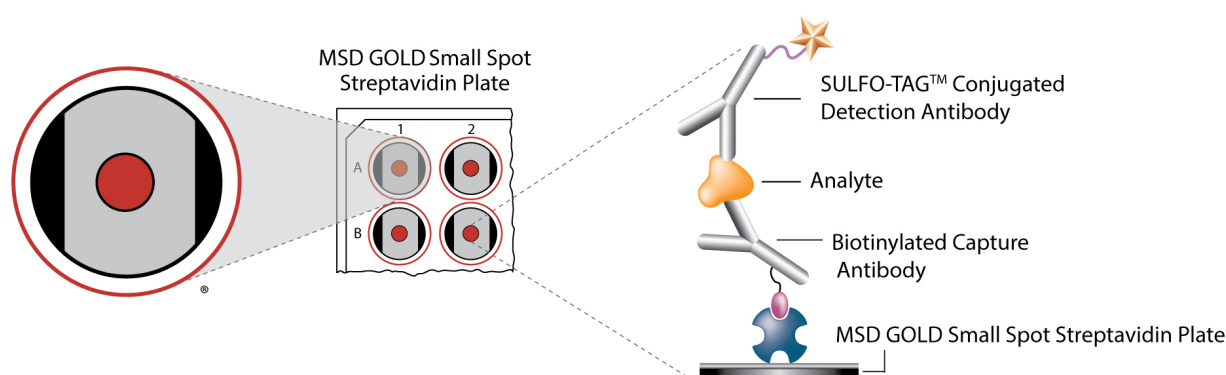


Figure 1: U-PLEX Singleplex Assay on an MSD GOLD 96-well Small Spot Streptavidin Plate. The 384-well assay is similar.

Components

Tables 2, 3, and 4 list the components provided with U-PLEX Biomarker Group 1 (NHP) Singleplex Assays. U-PLEX Singleplex Assays are available with either SECTOR™ or QuickPlex® plates or SECTOR 384-well plates (Table 6).

Reagents Supplied With All U-PLEX Singleplex Assays

Table 2: Reagents that are supplied with all U-PLEX biomarker Group 1 (NHP) 96-well Singleplex Assays

Reagent	Storage	Catalog No.	Size	Quantity Supplied			Description
				1 Plate	5 Plates	25 Plates	
MSD GOLD 96-Well Small Spot Streptavidin SECTOR Plate	2–8 °C	L45SA-1	1-spot	1 plate	5 plates	25 plates	96-well plate, foil sealed, with desiccant
MSD GOLD 96-Well Small Spot Streptavidin QuickPlex Plate		L4BSA-1					
Diluent 100	2–8 °C	R50AA-4	50 mL	1 bottle	1 bottle	5 bottles	Diluent for biotinylated capture antibody
Diluent 57	≤–10 °C	R50BZ-1	10 mL	1 bottle	—	—	Diluent for samples and Calibrator
		R50BZ-2	50 mL	—	1 bottle	5 bottles	
Diluent 3	≤–10 °C	R50AP-1	8 mL	1 bottle	—	—	Diluent for detection antibody
		R50AP-2	40 mL	—	1 bottle	5 bottles	
MSD GOLD Read Buffer B	RT	R60AM-1	18 mL	1 bottle	—	—	Buffer to catalyze the electrochemiluminescent reaction
		R60AM-2	90 mL	—	1 bottle	5 bottles	

RT = room temperature

Dash (—) = not applicable

Table 3: Reagents that are supplied with all U-PLEX Biomarker Group 1 (NHP) 384-well Singleplex Assays

Reagent	Storage	Catalog No.	Size	Quantity Supplied		Description
				5 Plates	25 Plates	
MSD 384-well Streptavidin SECTOR Plate	2–8 °C	L21SA-1	—	5 plates	25 plates	384-well plate, foil sealed, with desiccant
Diluent 100	2–8 °C	R50AA-4	50 mL	2 bottles	10 bottles	Diluent for biotinylated capture antibody
Diluent 57	≤–10 °C	R50BZ-2	50 mL	2 bottles	10 bottles	Diluent for samples and Calibrators
Diluent 3	≤–10 °C	R50AP-2	40 mL	2 bottles	10 bottles	Diluent for detection antibody
MSD GOLD Read Buffer B	RT	R60AM-2	90 mL	1 bottle	5 bottles	Buffer to catalyze the electrochemiluminescent reaction

RT = room temperature

Dash (—) = not applicable

Assay-Specific Reagents

U-PLEX Antibody Set

Based upon the analyte selected, you will receive a U-PLEX Antibody Set containing a biotinylated capture antibody and SULFO-TAG conjugated detection antibody. A complete list of all Antibody Sets available for U-PLEX Biomarker Group 1 (NHP) and their respective catalog numbers is provided in the Appendix (Table 9).

Table 4. Contents of U-PLEX Antibody Set

Name	Storage	Size	Quantity Supplied			Description
			1 Plate	5 Plates	25 Plates	
U-PLEX NHP Analyte-Specific Antibody Set	2–8 °C	1-Plate	1	—	—	Set containing biotinylated capture antibody and SULFO-TAG conjugated detection antibody
		5-Plate	—	1	5	

Dash (—) = not applicable

Calibrators

Calibrators are multianalyte blends, each containing multiple human proteins lyophilized in a buffered diluent.

Individual analyte concentrations are provided in the lot-specific certificates of analysis (COA). Based on the analyte selected, you will receive one of the following calibrators.

Table 5. Analytes included in the calibrator blends available for U-PLEX Biomarker Group 1 (NHP)

Name	Storage	Catalog No.	Size	Quantity Supplied			Analytes
				1 Plate	5 Plates	25 Plates	
Calibrator 1	2–8 °C	C0060-2	1 vial	1 vial	5 vials	25 vials	GM-CSF, IFN- γ , IL-1 β , IL-2, IL-4, IL-5, IL-6, IL-8, IL-10, IL-12p70, IL-17A, TNF- α , VEGF-A
Calibrator 2	2–8 °C	C0061-2	1 vial	1 vial	5 vials	25 vials	Eotaxin, Eotaxin-3, IP-10, MCP-1, MCP-4, MDC, MIP-1 α , MIP-1 β , TARC
Calibrator 3	2–8 °C	C0062-2	1 vial	1 vial	5 vials	25 vials	G-CSF, IL-1 α , IL-7, IL-12/IL-23p40, IL-15, IL-16, IL-18, TNF- β , TPO
Calibrator 4	2–8 °C	C0063-2	1 vial	1 vial	5 vials	25 vials	CTACK, Fractalkine, I-TAC, MIP-3 β , SDF-1 α
Calibrator 6	2–8 °C	C0072-2	1 vial	1 vial	5 vials	25 vials	IL-17A/F, IL-17E/IL-25, IL-17F, IL-21, IL-22, IL-23, IL-27, IL-29/IFN- λ 1, IL-31, IL-33, TSLP
Calibrator 9	2–8 °C	C0090-2	1 vial	1 vial	5 vials	25 vials	EPO, FLT3L, IFN- β , IL-1RA, IL-2R α , IL-3, IL-9, IL-17B, IL-17C, IL-17D
Calibrator 10	2–8 °C	C0091-2	1 vial	1 vial	5 vials	25 vials	Eotaxin-2, GRO- α , I-309, MCP-2, MCP-3, M-CSF, MIF, MIP-5, TRAIL, YKL-40

Instrument Compatibility

MSD offers U-PLEX Singleplex Assays designed for use on specific instrument platforms (Table 6).

Table 6. Instrument compatibility

Instrument	Assays on 96-well SECTOR Plates	Assays on 96-well QuickPlex Plates	Assays on 384-well SECTOR Plates
MESO® QuickPlex SQ 120	Y	—	—
MESO QuickPlex® SQ 120MM	Y	—	—
MESO SECTOR® S 600	Y	—	Y
MESO SECTOR S 600MM	Y	—	Y
MESO QuickPlex Q 60MM	—	Y	—

Dash (—) = not applicable

Additional Materials and Equipment

- ☐ Appropriately sized tubes for reagent preparation
- ☐ Polypropylene microcentrifuge tubes for preparing dilutions
- ☐ Liquid-handling equipment suitable for dispensing 10 to 150 µL/well into 96-well or 384-well microtiter plates
- ☐ Plate-washing equipment: automated plate washer or multichannel pipette
- ☐ Microtiter plate shaker (rotary) capable of shaking at 500–1,000 rpm (1,500 rpm for 384-well plates)
- ☐ MSD® Wash Buffer (20X, 100 mL, catalog number R61AA-1) for plate washing. The standard protocol uses a minimum of 415 mL of 1X Wash Buffer for a 384-well plate and 130 mL for a 96-well plate. Automated plate washers may need overage added to these volumes.
- ☐ Adhesive plate seals
- ☐ Deionized water
- ☐ Vortex mixer

Safety

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

Additional product-specific safety information is available in the applicable safety data sheet (SDS), which can be obtained from MSD Customer Service or at the www.mesoscale.com® website.

Best Practices

- Bring frozen diluents to room temperature in a 20–26 °C water bath before use. If a controlled water bath is not available, thaw at room temperature. Diluents may also be thawed overnight at 2–8°C.
- Ensure that diluents, wash buffer, and Read Buffer are equilibrated to room temperature before use. Mix well before use. Plates should be brought to room temperature before opening the foil packet.
- To avoid cross-contamination between vials, open vials for one protocol step at a time. Use filtered pipette tips, and use a fresh pipette tip for each reagent addition.
- MSD assays are tested and characterized between 21–26 °C; testing outside this temperature range may result in increased variability.
- Prepare calibrators, samples, and controls in a polypropylene container of sufficient volume.
- Avoid prolonged exposure of detection antibody (stock or diluted) to light. During the antibody incubation step, plates should not be exposed to direct sunlight.
- To ensure that all lyophilized powder is reconstituted, it is recommended that vials be inverted 3 times to distribute the diluent inside the vial. Then vortex the vial with 3 short pulses (upright, inverted, upright) after the solution sits at room temperature for the recommended amount of time in the product protocol.
- Ensure that all reagents are within their expiration date at the time of the test.
- For additional accuracy and precision, pre-wet pipette tips before transferring reagents and samples. Avoid pipetting bubbles while doing so.
- Plate shaking should be vigorous, with a rotary motion between 500 and 1,000 rpm for 96-well plates and 1,000–1,500 rpm for 384-well plates. Binding reactions may reach equilibrium sooner if you use shaking at the middle of the range or above. For long-term studies, the shaking speed and shaker model be kept consistent.
- Tap the plate on a paper towel after washing to ensure the removal of residual fluid.
- Consistent incubation times will improve the reproducibility of test results.
- Ensure that all necessary instruments, equipment, and reagents for the next step are prepared before washing the plates to prevent the plates from drying out.
- Avoid excessive drying of the plate during washing steps, especially if working inside a laminar flow hood or another high-airflow environment. Cover the plate with a new plate seal immediately after washing to protect from airflow and add solutions to the plate as soon as possible.
- Use a new adhesive plate seal for all incubation steps. Avoid re-using plate seals.
- Avoid creating bubbles in wells during all pipetting steps as they may lead to variable results.
- Use reverse pipetting when necessary and do not blow out residual liquid to avoid the introduction of bubbles. For empty wells, pipette gently to the bottom corner.
- Dispense reagents and wash fluids at the side of the well towards the bottom corner away from the coated spots.
- Protect plates from sources of heat such as vents, sunlight, etc. which may introduce variability across the plate surface. Some models of shakers generate heat that may affect plates on the platform.
- Ensure that all equipment is serviced and calibrated on a routine basis.
- Remove the plate seal before reading the plate.
- Read Buffer should be at room temperature (20–26 °C) before adding it to the plate.
- Keep time intervals consistent between the addition of Read Buffer and reading the plate to improve inter-plate precision. It is recommended that an MSD instrument be prepared to read a plate before adding Read Buffer. Unless otherwise directed, read the plate as soon as possible after adding Read Buffer.

- Do not shake the plate after adding Read Buffer.
- Do not obscure or damage the plate barcode; it is required for the plate reader.
- Only use the Read Buffer and wash buffer recommended for use with this kit.
- For 384-well assays, the protocol assumes the use of automated plate washers that can begin to aspirate before the total 90 μL is dispensed. If this ability is not present, reduce the wash volume to 80 μL to avoid overflowing the wells.

Reagent Preparation

Bring all reagents to room temperature and refer to the Best Practices section (page 9) before beginning the protocol.

Important: Upon the first thaw, aliquot diluents into suitable volumes before refreezing.

To prepare MSD Wash Buffer and other supplemental reagents, please refer to the Additional Materials and Equipment section (page 8).

Coat 96-well Plate

- ☐ Add 200 μL of biotinylated capture antibody to 3.3 mL of Diluent 100. Mix by vortexing.
- ☐ Add 25 μL of the above solution to each well of the provided MSD GOLD Small Spot Streptavidin Plate. Tap the plate gently on all sides. Seal the plate with an adhesive plate seal and incubate with shaking at room temperature for 1 hour.
- ☐ Wash the plate 3 times with at least 150 μL /well of 1X MSD Wash Buffer. The plate is now coated and ready for use. Plates may be sealed and stored overnight at 4 $^{\circ}\text{C}$.

Coat 384-well Plate

- ☐ Add 240 μL of biotinylated capture antibody to 11.76 mL of Diluent 100. Mix by vortexing.
- ☐ Add 25 μL of the above solution to each well of the provided plate. Tap the plate gently on all sides. Seal the plate with an adhesive plate seal and incubate with shaking at room temperature for 2 hours.
- ☐ Wash the plate 3 times with 90 μL /well of 1X MSD Wash Buffer. The plate is now coated and ready for use. Plates may be sealed and stored overnight at 4 $^{\circ}\text{C}$.

Prepare Calibrator Standards

Bring the calibrator vial(s) to room temperature. Reconstitute each vial of calibrator by adding 250 μL of Assay Diluent to the glass vial. This will result in a 10X concentrated stock of the calibrator, which will be diluted 5-fold (per the instructions given below) to generate the highest point in the standard curve (i.e., Calibrator Standard 1). A 2-fold dilution in the assay plate completes the 10-fold dilution. Invert the reconstituted calibrator at least 3 times. Do not vortex. Let the reconstituted solution equilibrate at room temperature for 15–30 minutes and then vortex briefly. The calibrator is now ready for use. Keep the dilutions at room temperature.

Note: We recommend that reconstituted or thawed calibrators be used immediately. If storage is necessary, divide calibrators into suitably sized aliquots (60 μL aliquots are recommended) and store immediately at $\leq -70^{\circ}\text{C}$.

The following instructions will enable you to prepare 7 Calibrator Standard solutions plus a zero Calibrator Standard for up to 6 replicates (see Figure 2; Table 7).

- ❑ Prepare Calibrator Standard 1 by adding 50 μL of the reconstituted or thawed calibrator to 200 μL of Assay Diluent. Mix by vortexing.
- ❑ For Calibrator Standard 2, add 75 μL of Calibrator Standard 1 to 225 μL of Assay Diluent. Mix by vortexing.
- ❑ Repeat 4-fold serial dilutions 5 additional times to generate a total of 7 Calibrator Standards. Mix by vortexing between each serial dilution.
- ❑ Use Assay Diluent as Calibrator Standard 8 (zero calibrator/blank).

Note: For the lot-specific concentration of calibrators in the blend, refer to the COA supplied with the assay pack. You can also find a copy of the COA at www.mesoscale.com.

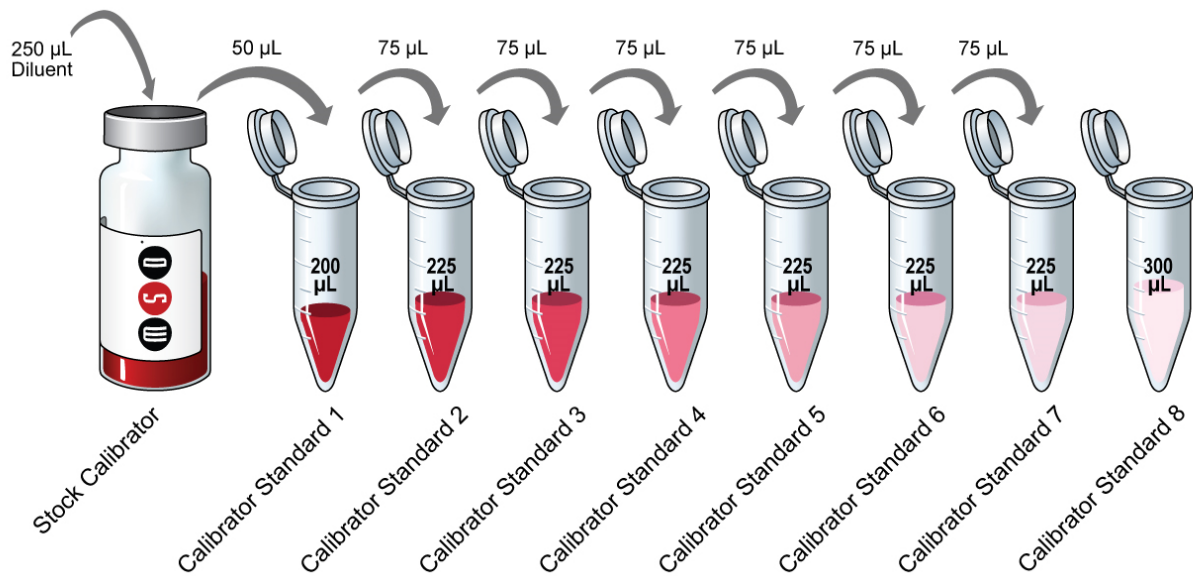


Figure 2. Dilution schema for Calibrator Standards for U-PLEX Biomarker Group 1 (NHP) Singleplex Assays

Table 7. Serial dilution to generate the standard curve

Calibrator Standard No.	Tube No.	Source of Calibrator	Volume of Reconstituted Calibrator (μL)	Assay Diluent (μL)	Total Volume (μL)
1	1	Stock Calibrator vial	50	200	250
2	2	From tube 1	75	225	300
3	3	From tube 2	75	225	300
4	4	From tube 3	75	225	300
5	5	From tube 4	75	225	300
6	6	From tube 5	75	225	300
7	7	From tube 6	75	225	300
8 (zero Calibrator)	8	—	0	300	300

Dash (—) = not applicable

Dilute Samples

Depending on the sample set under investigation, dilution may be necessary. Assay Diluent may be used for sample dilution. The dilution factor for the given sample type may need to be optimized.

Note: For MIF and YKL-40, the concentrations in normal serum and EDTA plasma may exceed the standard working range of the assays. Preassay dilution of samples may be required to generate optimal results. Refer to the product-specific datasheets for additional information. Diluent 100 may be used in place of Assay Diluent for samples that require high dilution.

Prepare Detection Antibody Solution

The detection antibody is provided as a 100X stock solution. The working solution is 1X for 96-well plates and 0.5X for 384-well plates. Prepare the detection antibody solution immediately before use.

For one plate, combine:

- ☐ 60 μ L of the supplied 100X detection antibody
- ☐ 5,940 μ L of Diluent 3 (11.94 mL for 384-well assays)

Read Buffer

MSD provides MSD GOLD Read Buffer B ready for use. Do not dilute.

Prepare Wash Buffer

Prepare a 1X working solution by diluting the 20X stock with deionized water. 1X MSD Wash Buffer can be stored at room temperature for up to two weeks.

Assay Protocols

96-well Plate Assays

Note: Follow Reagent Preparation before beginning this assay protocol.

STEP 1: Add Samples and Calibrators

- ☐ Add 25 μ L of Assay Diluent to each well. Tap the plate gently on all sides.
- ☐ Add 25 μ L of the prepared Calibrator Standard or sample to each well. Seal the plate with an adhesive plate seal. Incubate at room temperature with shaking for 1 hour.

STEP 2: Wash and Add Detection Antibody Solution

- ☐ Wash the plate 3 times with at least 150 μ L/well of 1X MSD Wash Buffer.
- ☐ Add 50 μ L of detection antibody solution to each well. Seal the plate with an adhesive plate seal. Incubate at room temperature with shaking for 1 hour.

STEP 3: Wash and Read

- ☐ Wash the plate 3 times with at least 150 μ L/well of 1X MSD Wash Buffer.
- ☐ Add 150 μ L of MSD GOLD Read Buffer B to each well. Analyze the plate on an MSD instrument. Incubation in Read Buffer is not required before reading the plate.

384-well Plate Assays

STEP 1: Add Samples and Calibrators

- ☐ Wash the plate 3 times with 90 μ L/well of 1X MSD Wash Buffer.
- ☐ Add 25 μ L of the prepared Calibrator Standard or sample to each well. Seal the plate with an adhesive plate seal. Incubate at room temperature with shaking for 2 hours.

STEP 2: Wash and Add Detection Antibody Solution

- ☐ Wash the plate 3 times with 90 μ L/well of 1X MSD Wash Buffer.
- ☐ Add 25 μ L of detection antibody solution to each well. Seal the plate with an adhesive plate seal. Incubate at room temperature with shaking for 2 hours.

STEP 3: Wash and Read

- ☐ Wash the plate 3 times with 90 μ L/well of 1X MSD Wash Buffer.
- ☐ Add 40 μ L of MSD GOLD Read Buffer B to each well. Analyze the plate on an MSD instrument. Incubation in Read Buffer is not required before reading the plate.

Alternate Protocols

The suggestions below may be useful for simplifying the protocol.

- ❑ **Alternate Protocol 1, Co-incubate:** Co-incubating samples and detection antibody solution may improve the sensitivity for some assays. Note that the use of the co-incubation protocol may result in sample concentrations that vary from concentrations obtained with the standard protocol. If this protocol is chosen, we recommend that this protocol be used for the entirety of the research project.
- ❑ **Alternate Protocol 2, Shortened Incubation:** Some 384-well assays may achieve acceptable performance with shorter incubations. Consider reducing the incubation time of samples in the plate and of detection antibody each to 1 hour.
- ❑ **Alternate Protocol 3, Reduced Wash:** For cell culture supernatants, you may simplify the protocol by eliminating one of the wash steps. After incubating the Calibrator Standard or sample, add detection antibody solution to the plate without decanting or washing the plate.

Assay Performance

A representative data set for each assay is presented in the product-specific datasheets available at www.mesoscale.com/U-PLEX-documents. The data represent the performance of the assay tested in multiplex format on U-PLEX plates. The data were generated during the development of the assay and do not represent the product specifications. Under your experimental conditions, the assay may perform differently than the representative data shown.

Appendix

U-PLEX Singleplex Assays

Assays include Antibody Sets, plates, Diluents, calibrators, and MSD GOLD Read Buffer B (Table 8).

Table 8. Catalog numbers of U-PLEX Biomarker Group 1 (NHP) Singleplex Assays

Product	96-Well SECTOR Assays (1/5/25 Plates)	96-Well QuickPlex Assays (1/5/25 Plates)	384-Well SECTOR Assays (5/25 Plates)
U-PLEX NHP CTACK Assay	K156VDK-1/-2/-4	K156VDK-21/-22/-24	K256VDK-2/-4
U-PLEX NHP Eotaxin Assay	K156UDK-1/-2/-4	K156UDK-21/-22/-24	K256UDK-2/-4
U-PLEX NHP Eotaxin-2 Assay	K156XQK-1/-2/-4	K156XQK-21/-22/-24	K256XQK-2/-4
U-PLEX NHP Eotaxin-3 Assay	K156UEK-1/-2/-4	K156UEK-21/-22/-24	K256UEK-2/-4
U-PLEX NHP ENA-78 Assay	K156VEK-1/-2/-4	K156VEK-21/-22/-24	K256VEK-2/-4
U-PLEX NHP FLT3L Assay	K156XFK-1/-2/-4	K156XFK-21/-22/-24	K256XFK-2/-4
U-PLEX NHP Fractalkine Assay	K156VCK-1/-2/-4	K156VCK-21/-22/-24	K256VCK-2/-4
U-PLEX NHP G-CSF Assay	K156VGK-1/-2/-4	K156VGK-21/-22/-24	K256VGK-2/-4
U-PLEX NHP GM-CSF Assay	K156UMK-1/-2/-4	K156UMK-21/-22/-24	K256UMK-2/-4
U-PLEX NHP GRO- α Assay	K156UXK-1/-2/-4	K156UXK-21/-22/-24	K256UXK-2/-4
U-PLEX NHP I-309 Assay	K156UYK-1/-2/-4	K156UYK-21/-22/-24	K256UYK-2/-4
U-PLEX NHP IFN- α 2a Assay	K156VHK-1/-2/-4	K156VHK-21/-22/-24	K256VHK-2/-4
U-PLEX NHP IFN- γ Assay	K156TTK-1/-2/-4	K156TTK-21/-22/-24	K256TTK-2/-4
U-PLEX NHP IL-1 α Assay	K156UNK-1/-2/-4	K156UNK-21/-22/-24	K256UNK-2/-4
U-PLEX NHP IL-1 β Assay	K156TUK-1/-2/-4	K156TUK-21/-22/-24	K256TUK-2/-4
U-PLEX NHP IL-1RA Assay	K156XPK-1/-2/-4	K156XPK-21/-22/-24	K256XPK-2/-4
U-PLEX NHP IL-2 Assay	K156TVK-1/-2/-4	K156TVK-21/-22/-24	K256TVK-2/-4
U-PLEX NHP IL-4 Assay	K156TWK-1/-2/-4	K156TWK-21/-22/-24	K256TWK-2/-4
U-PLEX NHP IL-5 Assay	K156UOK-1/-2/-4	K156UOK-21/-22/-24	K256UOK-2/-4
U-PLEX NHP IL-6 Assay	K156TXK-1/-2/-4	K156TXK-21/-22/-24	K256TXK-2/-4
U-PLEX NHP IL-7 Assay	K156UPK-1/-2/-4	K156UPK-21/-22/-24	K256UPK-2/-4
U-PLEX NHP IL-8 Assay	K156TYK-1/-2/-4	K156TYK-21/-22/-24	K256TYK-2/-4
U-PLEX NHP IL-9 Assay	K156XKK-1/-2/-4	K156XKK-21/-22/-24	K256XKK-2/-4
U-PLEX NHP IL-10 Assay	K156TZK-1/-2/-4	K156TZK-21/-22/-24	K256TZK-2/-4
U-PLEX NHP IL-12/IL-23p40 Assay	K156UQK-1/-2/-4	K156UQK-21/-22/-24	K256UQK-2/-4
U-PLEX NHP IL-12p70 Assay	K156UAK-1/-2/-4	K156UAK-21/-22/-24	K256UAK-2/-4

Table 8 (continued). Catalog numbers of U-PLEX Biomarker Group 1 (NHP) Singleplex Assays

Product	96-Well SECTOR Assays (1/5/25 Plates)	96-Well QuickPlex Assays (1/5/25 Plates)	384-Well SECTOR Assays (5/25 plates)
U-PLEX NHP IL-13 Assay	K156UBK-1/-2/-4	K156UBK-21/-22/-24	K256UBK-2/-4
U-PLEX NHP IL-15 Assay	K156URK-1/-2/-4	K156URK-21/-22/-24	K256URK-2/-4
U-PLEX NHP IL-16 Assay	K156USK-1/-2/-4	K156USK-21/-22/-24	K256USK-2/-4
U-PLEX NHP IL-17A Assay	K156UTK-1/-2/-4	K156UTK-21/-22/-24	K256UTK-2/-4
U-PLEX NHP IL-17A/F Assay	K156VYK-1/-2/-4	K156VYK-21/-22/-24	K256VYK-2/-4
U-PLEX NHP IL-17B Assay	K156XNK-1/-2/-4	K156XNK-21/-22/-24	K256XNK-2/-4
U-PLEX NHP IL-17C Assay	K156WJK-1/-2/-4	K156WJK-21/-22/-24	K256WJK-2/-4
U-PLEX NHP IL-17D Assay	K156XOK-1/-2/-4	K156XOK-21/-22/-24	K256XOK-2/-4
U-PLEX NHP IL-17F Assay	K156WAK-1/-2/-4	K156WAK-21/-22/-24	K256WAK-2/-4
U-PLEX NHP IL-18 Assay	K156VJK-1/-2/-4	K156VJK-21/-22/-24	K256VJK-2/-4
U-PLEX NHP IL-22 Assay	K156WIK-1/-2/-4	K156WIK-21/-22/-24	K256WIK-2/-4
U-PLEX NHP IL-23 Assay	K156W GK-1/-2/-4	K156W GK-21/-22/-24	K256W GK-2/-4
U-PLEX NHP IP-10 Assay	K156U FK-1/-2/-4	K156U FK-21/-22/-24	K256U FK-2/-4
U-PLEX NHP I-TAC Assay	K156U WK-1/-2/-4	K156U WK-21/-22/-24	K256U WK-2/-4
U-PLEX NHP MCP-1 Assay	K156U GK-1/-2/-4	K156U GK-21/-22/-24	K256U GK-2/-4
U-PLEX NHP MCP-2 Assay	K156X HK-1/-2/-4	K156X HK-21/-22/-24	K256X HK-2/-4
U-PLEX NHP MCP-4 Assay	K156U HK-1/-2/-4	K156U HK-21/-22/-24	K256U HK-2/-4
U-PLEX NHP M-CSF Assay	K156X RK-1/-2/-4	K156X RK-21/-22/-24	K256X RK-2/-4
U-PLEX NHP MDC Assay	K156U IK-1/-2/-4	K156U IK-21/-22/-24	K256U IK-2/-4
U-PLEX NHP MIF Assay	K156X JK-1/-2/-4	K156X JK-21/-22/-24	K256X JK-2/-4
U-PLEX NHP MIP-1 α Assay	K156U JK-1/-2/-4	K156U JK-21/-22/-24	K256U JK-2/-4
U-PLEX NHP MIP-1 β Assay	K156U KK-1/-2/-4	K156U KK-21/-22/-24	K256U KK-2/-4
U-PLEX NHP MIP-3 α Assay	K156U ZK-1/-2/-4	K156U ZK-21/-22/-24	K256U ZK-2/-4
U-PLEX NHP MIP-3 β Assay	K156V AK-1/-2/-4	K156V AK-21/-22/-24	K256V AK-2/-4
U-PLEX NHP MIP-5 Assay	K156X SK-1/-2/-4	K156X SK-21/-22/-24	K256X SK-2/-4
U-PLEX NHP SDF-1 α Assay	K156V BK-1/-2/-4	K156V BK-21/-22/-24	K256V BK-2/-4
U-PLEX NHP TARC Assay	K156U LK-1/-2/-4	K156U LK-21/-22/-24	K256U LK-2/-4
U-PLEX NHP TNF- α Assay	K156U CK-1/-2/-4	K156U CK-21/-22/-24	K256U CK-2/-4
U-PLEX NHP TNF- β Assay	K156U UK-1/-2/-4	K156U UK-21/-22/-24	K256U UK-2/-4
U-PLEX NHP TPO Assay	K156V KK-1/-2/-4	K156V KK-21/-22/-24	K256V KK-2/-4
U-PLEX NHP TRAIL Assay	K156X TK-1/-2/-4	K156X TK-21/-22/-24	K256X TK-2/-4
U-PLEX NHP VEGF-A Assay	K156U VK-1/-2/-4	K156U VK-21/-22/-24	K256U VK-2/-4
U-PLEX NHP YKL-40 Assay	K156V LK-1/-2/-4	K156V LK-21/-22/-24	K256V LK-2/-4

U-PLEX Antibody Sets

U-PLEX Antibody Sets (Table 9) include a biotinylated capture antibody and SULFO-TAG conjugated detection antibody.

Table 9. Catalog numbers of Antibody Sets available for U-PLEX Biomarker Group 1 (NHP)

Product	Cat. Nos. (-1/-5 Plate Sizes)	Product	Cat. Nos. (-1/-5 Plate Sizes)
U-PLEX Human CTACK Antibody Set	B21VD-2/-3	U-PLEX Human IL-17A/F Antibody Set	B21VY-2/-3
U-PLEX NHP Eotaxin Antibody Set	B26UD-2/-3	U-PLEX Human IL-17B Antibody Set	B21XN-2/-3
U-PLEX Human Eotaxin-2 Antibody Set	B21XQ-2/-3	U-PLEX Human IL-17C Antibody Set	B21WJ-2/-3
U-PLEX Human Eotaxin-3 Antibody Set	B21UE-2/-3	U-PLEX Human IL-17D Antibody Set	B21XO-2/-3
U-PLEX NHP ENA-78 Antibody Set	B26VE-2/-3	U-PLEX Human IL-17F Antibody Set	B21WA-2/-3
U-PLEX Human FLT3L Antibody Set	B21XF-2/-3	U-PLEX Human IL-18 Antibody Set	B21VJ-2/-3
U-PLEX Human Fractalkine Antibody Set	B21VC-2/-3	U-PLEX Human IL-22 Antibody Set	B21WI-2/-3
U-PLEX Human G-CSF Antibody Set	B21VG-2/-3	U-PLEX Human IL-23 Antibody Set	B21WG-2/-3
U-PLEX Human GM-CSF Antibody Set	B21UM-2/-3	U-PLEX Human IP-10 Antibody Set	B21UF-2/-3
U-PLEX Human GRO- α Antibody Set	B21UX-2/-3	U-PLEX Human I-TAC Antibody Set	B21UW-2/-3
U-PLEX Human I-309 Antibody Set	B21UY-2/-3	U-PLEX Human MCP-1 Antibody Set	B21UG-2/-3
U-PLEX NHP IFN- α 2a Antibody Set	B26VH-2/-3	U-PLEX Human MCP-2 Antibody Set	B21XH-2/-3
U-PLEX Human IFN- γ Antibody Set	B21TT-2/-3	U-PLEX Human MCP-4 Antibody Set	B21UH-2/-3
U-PLEX NHP IL-1 α Antibody Set	B26UN-2/-3	U-PLEX Human M-CSF Antibody Set	B21XR-2/-3
U-PLEX Human IL-1 β Antibody Set	B21TU-2/-3	U-PLEX Human MDC Antibody Set	B21UI-2/-3
U-PLEX NHP IL-1RA Antibody Set	B26XP-2/-3	U-PLEX Human MIF Antibody Set	B21XJ-2/-3
U-PLEX NHP IL-2 Antibody Set	B26TV-2/-3	U-PLEX Human MIP-1 α Antibody Set	B21UJ-2/-3
U-PLEX NHP IL-4 Antibody Set	B26TW-2/-3	U-PLEX Human MIP-1 β Antibody Set	B21UK-2/-3
U-PLEX Human IL-5 Antibody Set	B21UO-2/-3	U-PLEX NHP MIP-3 α Antibody Set	B26UZ-2/-3
U-PLEX Human IL-6 Antibody Set	B21TX-2/-3	U-PLEX Human MIP-3 β Antibody Set	B21VA-2/-3
U-PLEX Human IL-7 Antibody Set	B21UP-2/-3	U-PLEX Human MIP-5 Antibody Set	B21XS-2/-3
U-PLEX Human IL-8 Antibody Set	B21TY-2/-3	U-PLEX NHP SDF-1 α Antibody Set	B26VB-2/-3
U-PLEX Human IL-9 Antibody Set	B21XK-2/-3	U-PLEX NHP TARC Antibody Set	B26UL-2/-3
U-PLEX Human IL-10 Antibody Set	B21TZ-2/-3	U-PLEX NHP TNF- α Antibody Set	B26UC-2/-3
U-PLEX Human IL-12/IL-23p40 Antibody Set	B21UQ-2/-3	U-PLEX Human TNF- β Antibody Set	B21UU-2/-3
U-PLEX NHP IL-12p70 Antibody Set	B26UA-2/-3	U-PLEX Human TPO Antibody Set	B21VK-2/-3
U-PLEX NHP IL-13 Antibody Set	B26UB-2/-3	U-PLEX Human TRAIL Antibody Set	B21XT-2/-3
U-PLEX Human IL-15 Antibody Set	B21UR-2/-3	U-PLEX Human VEGF-A Antibody Set	B21UV-2/-3
U-PLEX Human IL-16 Antibody Set	B21US-2/-3	U-PLEX Human YKL-40 Antibody Set	B21VL-2/-3
U-PLEX Human IL-17A Antibody Set	B21UT-2/-3		

Summary Protocols

Coat 96-well Plate

- ☐ Add 200 μ L of biotinylated capture antibody to 3.3 mL of Diluent 100. Mix by vortexing.
- ☐ Add 25 μ L of the above solution to each well of the provided MSD GOLD Small Spot Streptavidin Plate. Seal the plate with an adhesive plate seal and shake for 1 hour at room temperature.
- ☐ Wash the plate 3 times with at least 150 μ L/well of 1X MSD Wash Buffer. The plate is now coated and ready for use.

96-well Assay Protocol

STEP 1: Add Samples and Calibrators

- ☐ Add 25 μ L of Diluent 57 to each well. Tap the plate gently on all sides.
- ☐ Add 25 μ L of prepared Calibrator Standard or sample to each well. Seal the plate with an adhesive plate seal. Incubate at room temperature with shaking for 1 hour.

STEP 2: Wash and Add Detection Antibody Solution

- ☐ Wash the plate 3 times with at least 150 μ L/well of 1X Wash Buffer.
- ☐ Add 50 μ L of detection antibody solution to each well. Seal the plate with an adhesive plate seal and incubate at room temperature with shaking for 1 hour.

STEP 3: Wash and Read

- ☐ Wash the plate 3 times with at least 150 μ L/well of 1X Wash Buffer.
- ☐ Add 150 μ L of MSD GOLD Read Buffer B to each well. Analyze the plate on an MSD instrument. Incubation in Read Buffer is not required before reading the plate.

Coat 384-well Plate

- ☐ Add 240 μL of biotinylated capture antibody to 11.76 mL of Diluent 100. Mix by vortexing.
- ☐ Add 25 μL of the above solution to each well of the provided MSD 384-well Streptavidin Plate. Seal the plate with an adhesive plate seal and shake for 2 hours at room temperature.
- ☐ Wash the plate 3 times with 90 μL /well of 1X MSD Wash Buffer. The plate is now coated and ready for use and may be stored overnight at 4 °C.

384-well Assay Protocol

STEP 1: Add Samples and Calibrators

- ☐ Wash the plate 3 times with 90 μL /well of 1X Wash Buffer.
- ☐ Add 25 μL of prepared Calibrator Standard or sample to each well. Seal the plate with an adhesive plate seal. Incubate at room temperature with shaking for 2 hours at room temperature.

STEP 2: Wash and Add Detection Antibody Solution

- ☐ Wash the plate 3 times with 90 μL /well of 1X Wash Buffer. Add 25 μL of detection antibody solution to each well. Seal the plate with an adhesive plate seal and incubate at room temperature with shaking for 2 hours at room temperature.

STEP 3: Wash and Read

- ☐ Wash the plate 3 times with 90 μL /well of 1X Wash Buffer.
Add 40 μL of MSD GOLD Read Buffer B to each well. Analyze the plate on an MSD instrument. Incubation in Read Buffer is not required before reading the plate.

Plate Diagrams

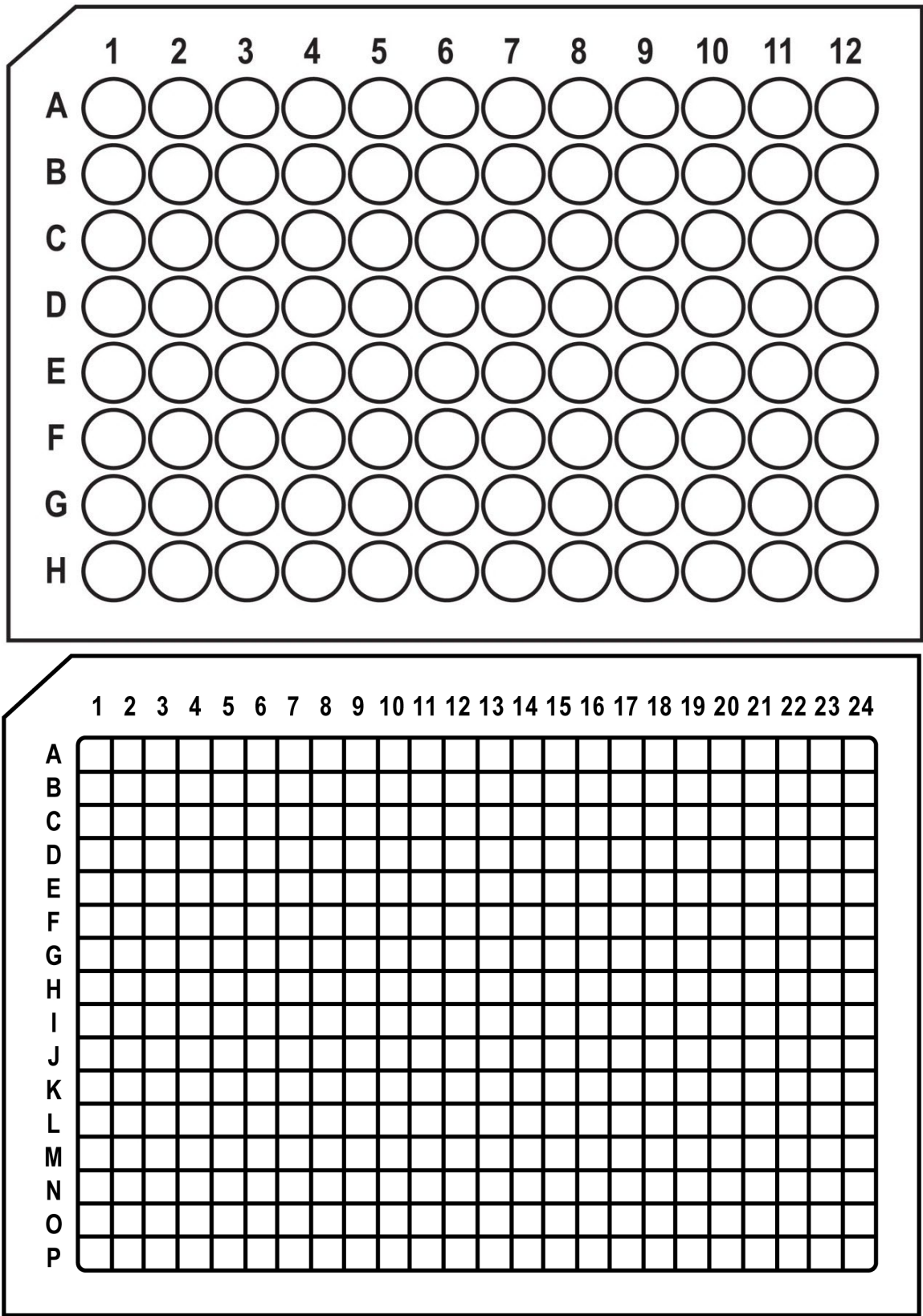


Figure 3: Plate diagrams. Similar plate layouts can be created in Excel and in the DISCOVERY WORKBENCH® software.