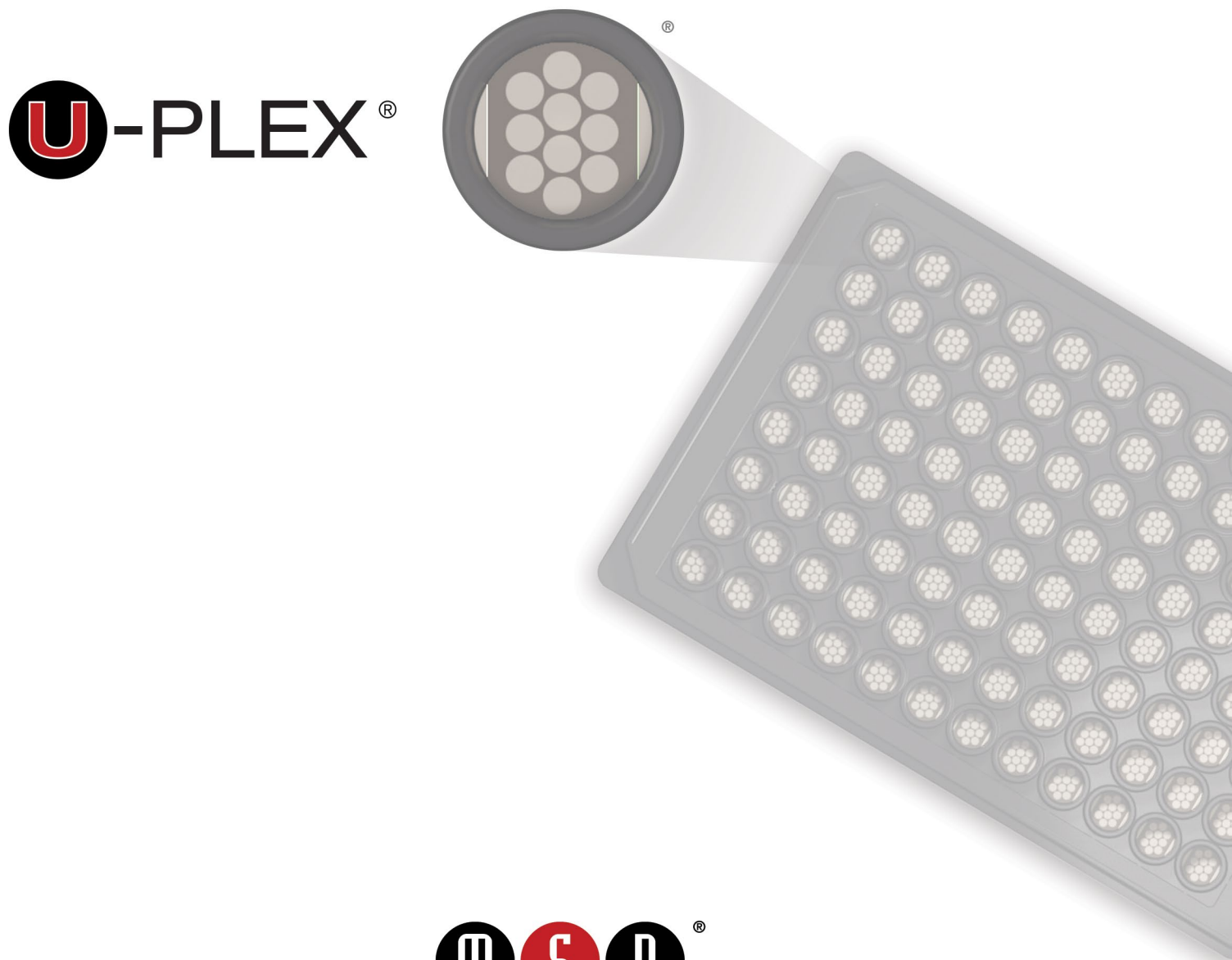


## U-PLEX<sup>®</sup> Metabolic Group 1 (mouse) Multiplex Assays



# MSD U-PLEX Platform

## U-PLEX Metabolic Group 1 (mouse) Multiplex Assays

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

For use with:

U-PLEX Metabolic Group 1 (mouse) Assays (catalog numbers K152ACM-1, K152ACM-2, K152ACM-4)

U-PLEX Metabolic Group 1 (mouse) 384-well Assays (catalog numbers K252ACM-2, K252ACM-4)

U-PLEX Metabolic Group 1 (mouse) Combos (catalog numbers are available at [https://www.mesoscale.com/en/products\\_and\\_services/assay\\_kits/u-plex\\_gateway/u-plex-combinations](https://www.mesoscale.com/en/products_and_services/assay_kits/u-plex_gateway/u-plex-combinations))

*This product 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](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, MesoSphere, Methodical Mind, Methodical Mind Enterprise, 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, 2020-2023 Meso Scale Diagnostics, LLC. All rights reserved.

# Table of Contents

Introduction .....	4
Principle of the Assay .....	4
Components.....	5
Additional Materials and Equipment .....	8
Safety .....	9
Assay Protocol (96-well plates) .....	15
Assay Performance .....	16
Appendix A .....	17
Appendix B .....	20
Assay Protocol (384-well plates) .....	22
Plate Diagrams.....	24

## 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

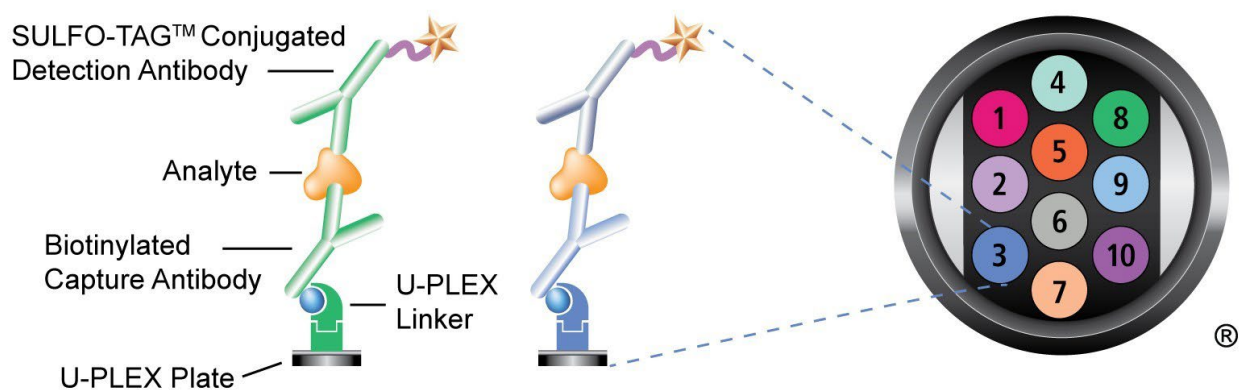
The U-PLEX platform allows the creation of custom multiplex assays for any combination of analytes in the same Group.

This product insert is for U-PLEX Combos and multiplex assays that contain any number of the 58 assays in the U-PLEX Metabolic Group 1 (mouse). Using open spots, custom combinations can include R-PLEX<sup>®</sup> Antibody Sets or your antibodies for other analytes.

Representative data for each U-PLEX assay is presented in the product-specific datasheets available at the [www.mesoscale.com](http://www.mesoscale.com)<sup>®</sup> website. The performance of MSD assays may vary when tested in combination. The data presented in the datasheets do not represent the product specifications.

## Principle of the Assay

Biotinylated capture antibodies are coupled to U-PLEX Linkers, which self-assemble onto unique spots on the U-PLEX plate. Analytes in the sample bind to the capture reagents. Detection antibodies conjugated with electrochemiluminescent labels (MSD GOLD<sup>™</sup> SULFO-TAG) bind to the analytes to complete the sandwich immunoassay (Figure 1). Once the sandwich immunoassay is complete, the U-PLEX 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 multiplex immunoassay on a U-PLEX 96-well 10-Assay Plate. U-PLEX 384-well 4-Assay plates are similar.*

# Components

Tables 1, 2, and 3 list the components provided with multiplex U-PLEX Metabolic Group 1 (mouse) Assays. You will only receive components relevant to the assays that you order.

**Table 1.** Reagents that are supplied with all U-PLEX Metabolic Group 1 (mouse) 96-well Assays

Reagent	Storage	Catalog No.	Size	Quantity Supplied			Description
				1 Plate	5 Plates	25 Plates	
Diluent 13	≤-10 °C	R56BB-4	10 mL	1 bottle	—	—	Diluent for samples and calibrators
		R56BB-3	50 mL	—	1 bottle	5 bottles	
Diluent 11	≤-10 °C	R55BA-5	10 mL	1 bottle	—	—	Diluent for detection antibody
		R55BA-3	50 mL	—	1 bottle	5 bottles	
Aprotinin (200,000 KIU/mL)	2-8 °C	R93AD-1	50 µL	1 vial	—	—	200X concentrated solution
		R93AD-2	250 µL	—	1 vial	5 vials	
Stop Solution	2-8 °C	R50AO-1	40 mL	1 bottle	1 bottle	5 bottles	Biotin-containing buffer to stop Linker-antibody coupling reaction
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	

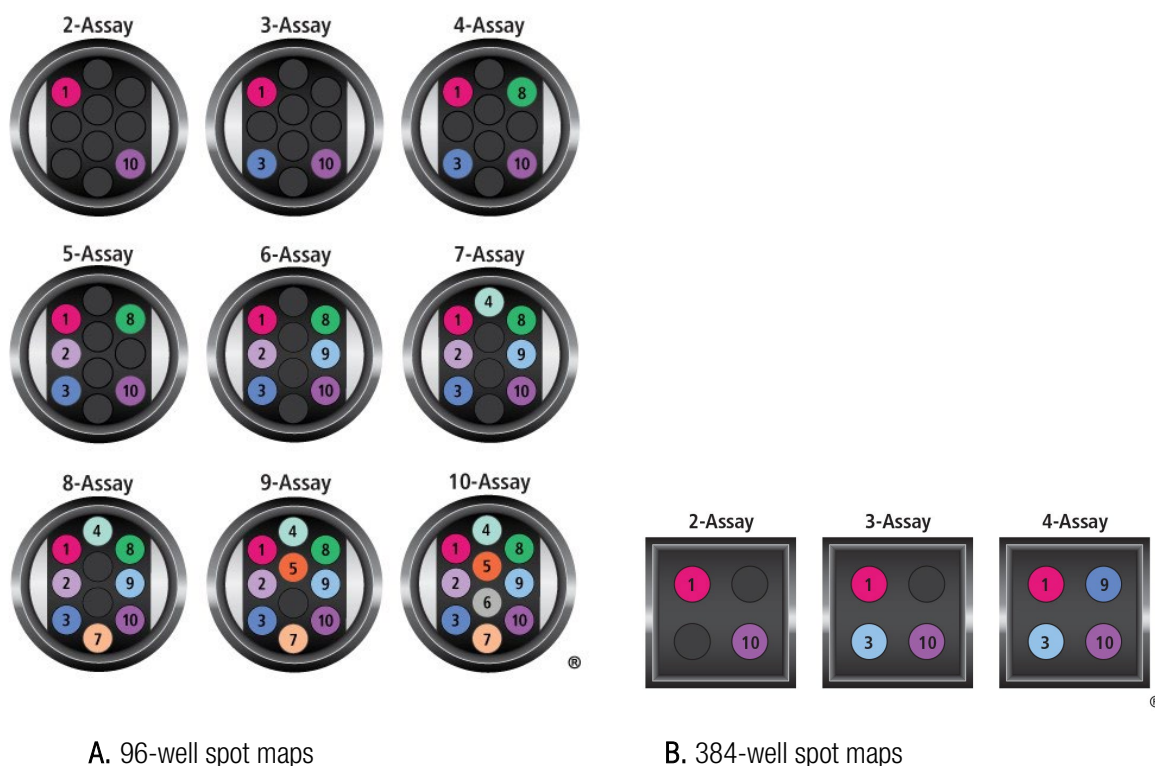
RT = room temperature  
Dash (—) = not applicable

## Assay-Specific Reagents

U-PLEX plates are provided in a sealed foil pouch with desiccant. The spots correspond to unique U-PLEX Linkers. The number and layout of the active spots on the plate depend on the plate well density (96 vs 384) and the number of assays to be multiplexed (Figure 2). For example, if 4 assays are to be multiplexed, either a U-PLEX 96-well or 384-well 4-Assay Plate will be provided.

### U-PLEX Plates

U-PLEX multiplex assays include plates that are specific to the assay with which they ship. Do not separate plates from the other refrigerated components.



**Figure 2.** Spot Maps of the different U-PLEX multiplex plates showing the placement of Linkers within a well. The colored spots represent the active U-PLEX binding spots. The numbering convention for the different spots is maintained in the software visualization tools and in the data files.

## Linkers

Each Linker has a biotin-binding domain that couples to the biotinylated capture antibody, as well as a domain that binds to its matching spot on the U-PLEX plate. The Linkers are color-coded and numbered with the spot to which they attach on the plate.

Record which antibody is coupled to each Linker when performing the coupling step (as described in the Reagent Preparation section).

## U-PLEX Antibody Sets

You will receive U-PLEX Antibody Sets containing the biotinylated capture antibody and the SULFO-TAG™ conjugated detection antibody (Table 2).

**Table 2.** Contents of U-PLEX Antibody Sets

Name	Storage	Size	Quantity Supplied			Description
			1 Plate	5 Plates	25 Plates	
U-PLEX Mouse 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

## U-PLEX Calibrators

Calibrators may be either lyophilized or frozen. Individual analyte concentrations are provided in lot-specific certificates of analysis (COA). Depending on the specific assays requested, one or more of the following Calibrators will be provided (Table 3).

**Table 3.** Analytes included in the Calibrators available for U-PLEX Metabolic Group 1 (mouse)

Name	Storage	Catalog No.	Analytes
Calibrator 5	2–8 °C	C0065-2	EPO, GM-CSF, IFN- $\gamma$ , IL-1 $\beta$ , IL-2, IL-4, IL-5, IL-6, IL-10, IL-12p70, IL-13, KC/GRO, TNF- $\alpha$ , VEGF-A
Calibrator 7	2–8 °C	C0073-2	IL-16, IL-17A, IL-17C, IL-17E/IL-25, IL-21, IL-22, IL-23
Calibrator 8	2–8 °C	C0074-2	IL-15, IL-17F, IL-31, IL-33, IL-27p28/IL-30
Calibrator 12	2–8 °C	C0092-2	IL-9, IL-17A/F, IP-10, MCP-1, MIP-1 $\alpha$ , MIP-1 $\beta$ , MIP-2, MIP-3 $\alpha$
Calibrator 16	$\leq$ -70 °C	C0295-2	6CKine/CCL21, BAFF, BCA-1/BLC, CD40, IFN- $\beta$ , MCP-5/CCL12, MDC
Calibrator 17	2–8 °C	C0296-2	Eotaxin, MMP-9 (total), NGAL/LCN2, RANTES, SDF-1 $\alpha$ , TARC, TNF-RI
Calibrator 18	2–8 °C	C0292-2	BDNF, C-Peptide, FGF-21, Ghrelin, GLP-1 (inactive), GLP-1 (total), PYY (total)
Calibrator 19	2–8 °C	C0293-2	Glucagon, Insulin, Leptin
Ghrelin (active)	2–8 °C	C016K-2	Ghrelin (active)
GLP-1 (active)	2–8 °C	C016L-2	GLP-1 (active)
Mouse IFN- $\alpha$ Calibrator	2–8 °C	C02W1-2	IFN- $\alpha$

# 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  $\mu\text{L}$ /well into a 96-well microtiter plate.
- Plate-washing equipment: automated plate washer or multichannel pipette.
- Microtiter plate shaker (rotary) capable of shaking at 500–1,000 rpm (1,000–1,500 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.
- A dipeptidyl peptidase IV (DPP-IV) inhibitor (such as diprotin A) is recommended for certain assays.
- MSD Blocker D–R (catalog number R93BR) is recommended for certain assays.
- Halt Protease Inhibitor Cocktail, EDTA-Free (Thermo Scientific, catalog number 87785). Recommended for Ghrelin (active) assay. Use as recommended by the manufacturer.
- Vortex mixer.

**Note:** If including Open Spots, you will also need:

- MSD GOLD SULFO-TAG NHS-Ester (catalog no. R91AO-1) for conjugating detection reagents or SULFO-TAG conjugated antispecies antibodies for use as reporters with unconjugated detection antibodies.
- Sulfo-NHS-LC-Biotin for biotinylating the capture reagents (e.g., EZ-Link Sulfo-NHS-LC-Biotin, Thermo Fisher Scientific, catalog no. 21327, or equivalent).
- Zeba Desalting Columns (Thermo Fisher Scientific, catalog numbers 87766-87773).
- Coating diluent such as 0.5% bovine serum albumin in PBS, or MSD Diluent 100 for diluting the capture antibody.

## Instrument Compatibility

MSD offers U-PLEX Assays designed for use on specific instrument platforms (Table 4).

**Table 4.** Instrument compatibility

Instrument	Assays on U-PLEX 96-well SECTOR™ Plate	Assays on U-PLEX 384-well SECTOR Plate
MESO® QuickPlex SQ 120	Y	—
MESO QuickPlex® SQ 120MM	Y	—
MESO SECTOR® S 600	Y	Y
MESO SECTOR S 600MM	Y	Y

Dash (—) = not applicable



# 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.

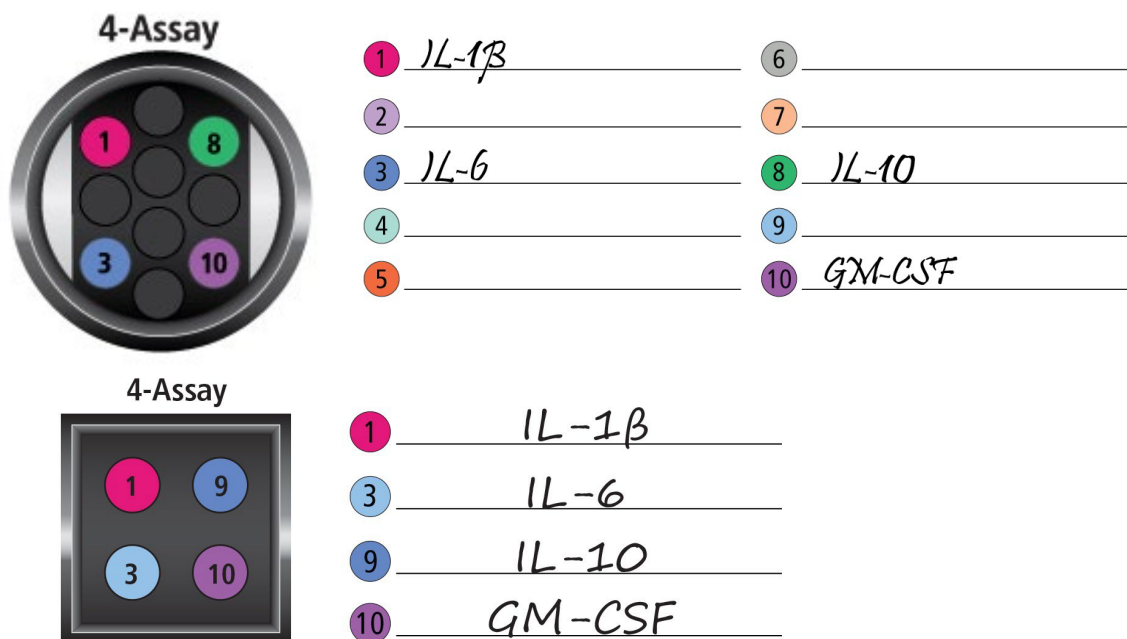
Additional product-specific safety information is available in the applicable safety data sheet(s) (SDS), which can be obtained from MSD Customer Service or at [www.mesoscale.com](http://www.mesoscale.com).

## Reagent Preparation for 96-well Plates

**Important:** Upon the first thaw, aliquot diluents into suitably sized aliquots before refreezing.

### Prepare U-PLEX Plate

The preparation of a U-PLEX plate involves coating the provided plate with Linker-coupled capture antibodies. Assign each antibody to a unique Linker and record the antibody identity next to the assigned Linker, as shown in the examples below (Figure 3).



**Figure 3.** U-PLEX 4-Assay Plates and assigned Linkers: *(top)* 96-well plate, *(bottom)* 384-well plate.

This section describes the preparation of a multiplex coating solution for one plate. The volumes can be adjusted depending on the number of plates or wells, but the ratios of the reagents should remain the same (Table 5).

### STEP 1: Create Individual U-PLEX Linker-Coupled Antibody Solutions

A different Linker is used for each unique biotinylated antibody. Below are the steps to couple linkers with capture antibodies.

- Couple each biotinylated capture antibody to a unique Linker and record the antibody identity next to the Linker number on the Spot Map (blank Spot Maps are provided on page 23).
- Add 200  $\mu$ L of each biotinylated antibody to 300  $\mu$ L of the assigned Linker. Mix by vortexing. Incubate at room temperature for 30 minutes. Do not shake.

#### Notes:

- Each Linker vial has a matching colored cap and label.
  - To remove liquid from the cap, briefly centrifuge the Linker vial and open the cap gently.
  - Open one Linker at a time and close its cap as soon as you are done using it. Take precautions to avoid reagent contamination.
  - For studies using multiple plates of the same assay, it is recommended that the same Linker be coupled with the same antibody for the duration of the study.
- Add 200  $\mu$ L of Stop Solution, then mix by vortexing. Incubate at room temperature for 30 minutes.

**Note:** At the end of step 1, each individual U-PLEX Linker-coupled antibody solution is at 10X the coating concentration and can be stored at 2–8 °C. Do not store for more than 7 days.

Adjust the volumes for multiple plates (see Table 5). The volumetric ratio of Linker:antibody:Stop Solution is 3:2:2.

## STEP 2: Prepare the Multiplex Coating Solution for 96-well Plates

- ❑ Combine 600 µL of each U-PLEX Linker-coupled antibody solution into a 15 mL tube and mix by vortexing. Up to 10 U-PLEX Linker-coupled antibodies can be pooled. Do not combine U-PLEX Linker-coupled antibody solutions that share the same Linker.
- ❑ When combining fewer than 10 antibodies, bring the solution up to 6 mL with Stop Solution. This will result in a final 1X concentration. Mix by vortexing. For example, for a 4-assay coating solution, add 3.6 mL of Stop Solution to the 2.4 mL of combined antibodies.

**Note:** At the end of Step 2, the U-PLEX multiplex coating solution is at 1X and can be stored at 2–8 °C. Do not store for more than 7 days.

## STEP 3: Coat U-PLEX 96-well Plates

- ❑ Add 50 µL of multiplex coating solution to each well. Seal the plate with an adhesive plate seal and incubate at room temperature while shaking for 1 hour.
- ❑ Wash the plate 3 times with at least 150 µL/well of 1X MSD Wash Buffer.

**Note:** The plate is now coated and ready for use. Sealed plates may be stored in the original pouch for up to 7 days at 2–8 °C.

The recommended volumes of Linker, biotinylated capture antibody, and Stop Solution for coating one or multiple U-PLEX plates are provided below in Table 5. If using a partial plate, adjust the volumes used proportionally.

**Table 5.** Amount of each component required for U-PLEX coating solution per plate

No. of Plates	Individual Linker (µL)	Individual Biotinylated Antibody (µL)	Stop Solution (µL)
1	300	200	200
2	600	400	400
3	900	600	600
4	1,200	800	800
5	1,500	1,000	1,000
N	300 × N	200 × N	200 × N

## Prepare Metabolic Assay Working Solution

This procedure is used for preparing the calibrator, controls, and diluting the samples.

For one plate, combine the following in a 15 mL tube:

- 6,965  $\mu$ L of Diluent 13
- 35  $\mu$ L of aprotinin

### Notes:

- The addition of a DPP-IV inhibitor (such as diprotin A, not provided) to a final concentration of 0.1 mM in the Metabolic Assay Working Solution is strongly recommended. A DPP-IV inhibitor will aid in limiting the enzymatic action of DPP-IV present in serum/plasma and provides the most accurate measurement of some metabolic analytes.
- Ghrelin (active) assay only: Addition of Halt Protease Inhibitor Cocktail, EDTA-Free (Thermo Fisher Scientific, Catalog No. 87785) to a final concentration of 1X in the Metabolic Assay Working Solution is strongly recommended. The presence of Halt inhibitor may affect sample quantitation in other assays.

**Important:** Protease inhibitors should be added before use. The Metabolic Assay Working Solution should be kept on ice. Do not freeze the Metabolic Assay Working Solution for later use.

## Prepare Calibrator Standards

### For Lyophilized Calibrators

Bring the Calibrators to room temperature. Reconstitute by adding 250  $\mu$ L of Metabolic Assay Working Solution to the vial. This will result in a 10X concentrated stock of each Calibrator. 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.

### For Liquid Calibrators

Thaw the stock Calibrator(s) and gently mix. Keep on ice. Once thawed, the Calibrator is ready to use.

**Note:** We recommend that reconstituted or thawed Calibrators be used immediately. If storage is necessary, divide Calibrators into suitably sized aliquots (60  $\mu$ L aliquots are recommended) and store immediately at  $\leq -70$  °C.

Prepare Calibrator Standard 1 (top of the curve) in a polypropylene tube by mixing and diluting the reconstituted or thawed Calibrator as indicated in Table 6. Mix by vortexing.

**Table 6.** Combining Calibrators to generate the Calibrator Standard 1 (top of the curve) level

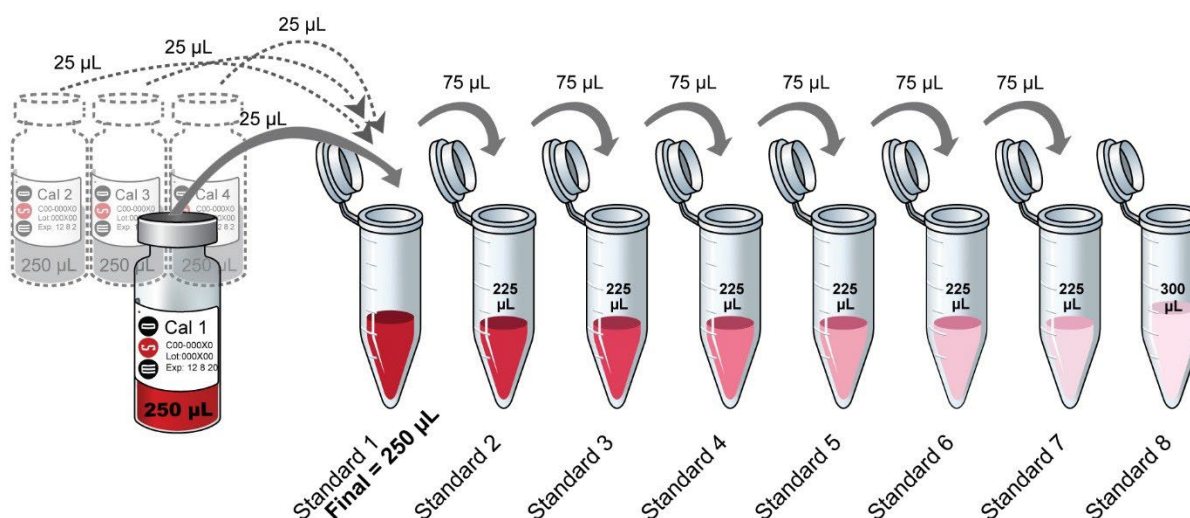
No. of Calibrators Provided	Volume of Reconstituted Calibrator ( $\mu$ L)	Metabolic Assay Working Solution ( $\mu$ L)	Total volume ( $\mu$ L)
2	25 each	200	250
3	25 each	175	250
4	25 each	150	250
N	25 each	250 - (25 x N)	250

Prepare the subsequent 6 dilutions for the curve (4-fold serial dilutions) in Metabolic Assay Working Solution (see Figure 4; Table 7). Use Metabolic Assay Working Solution for the Calibrator Standard 8 (zero Calibrator/blank). Mix by vortexing the tubes between each serial dilution.

**Table 7.** Serial dilution to generate the standard curve

Calibrator Standard No.	Tube No.	Source of Calibrator	Volume of Reconstituted Calibrator (µL)	Metabolic Assay Working Solution (µL)	Total volume (µL)
1	1	Calibrator Standard 1 (top of curve)	See Table 6		
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



**Figure 4.** Dilution schema for preparation of Calibrator Standards for U-PLEX Metabolic Group 1 (mouse) Assays.

## Sample Collection and Handling

Below are general guidelines for sample collection, storage, and handling for metabolic markers. We strongly suggest following these procedures if working with the active forms of protein analytes. If other methods are used, evaluate sample stability under the selected method as needed.

The assay requires 50 µL/well of the sample (25 µL for 384-well plates). Based on the number of replicates desired, prepare an adequate volume of the sample.

**Sample Collection (preferred method):** Samples should be collected using the BD P800 Collection and Preservation System, which contains a DPP-IV and other protease inhibitor cocktails (Product Number 366420 or 366421). The alternative collection method described below with K<sub>2</sub>EDTA tubes can also be used.

**Non-P800 collection method:** Collect blood in BD Vacutainer K<sub>2</sub>EDTA Tubes (Product Number 367841 or 366643). **Immediately** add a DPP-IV inhibitor (0.1 mM final concentration, not provided) and aprotinin (1,000 KIU/mL final concentration) and mix to avoid cleavage/degradation of metabolic peptides.

For BD tubes, process as follows:

- In a swing-out rotor centrifuge, spin the blood collection tubes as follows;
  - For 2 mL tubes —10 minutes at  $1,000 \times g$  (2–8 °C).
  - For 8.5 and 10 mL tubes—20 minutes at  $1,300 \times g$  (2–8 °C).
- Use the plasma immediately or the samples can be stored at 2–8 °C if used within 3 hours. For future use, aliquot the plasma and freeze in suitably sized aliquots at  $\leq -70$  °C.

For samples other than serum and plasma, add a DPP-IV inhibitor (0.1 mM final concentration, not provided) and aprotinin (1,000 KIU/mL final concentration) and use immediately or freeze at  $\leq -70$  °C.

Samples with hemolysis or significant lipemia may hinder accurate measurements.

Repeated freezing and thawing of samples is not recommended. After thawing, centrifuge samples at  $2,000 \times g$  for 3 minutes to remove particulates before use in the assay. If the samples are clear and no particulates are visible, you may not need to centrifuge. Hold on wet ice or 2–8 °C until processed and used in the assay.

## Dilute Samples

Depending on the sample set under investigation, dilution may be necessary. At least a two-fold dilution of serum and plasma is recommended to reduce matrix effects. Metabolic Assay Working Solution should be used for sample dilution. The dilution factor for the given sample type may need to be optimized.

**Note:** For BAFF, the concentrations in normal serum and EDTA plasma may exceed the standard working range of the assays. Refer to the product-specific datasheets for additional information.

## Prepare Detection Antibody Solution

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

For one plate, combine:

- 60  $\mu$ L of each 100X detection antibody
- Diluent 11 to bring the final volume to 6 mL

## Wash Buffer

Prepare a 1X working solution by diluting the 20X stock with deionized water.

## Read Buffer

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

# Assay Protocol (96-well plates)

**Note:** Follow Reagent Preparation before beginning this assay protocol.

## STEP 1: Add Sample or Calibrator Standard

- Add 50  $\mu\text{L}$  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 at least 150  $\mu\text{L}$ /well 1X MSD Wash Buffer.
- Add 50  $\mu\text{L}$  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  $\mu\text{L}$ /well of 1X MSD Wash Buffer.
- Add 150  $\mu\text{L}$  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-incubation:** 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, 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](http://www.mesoscale.com). 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 and with your specific multiplex, the assay may perform differently than the representative data shown.

## Specificity

To assess specificity, the Antibody Set for each analyte was tested individually against a larger panel of recombinant mouse analytes for nonspecific binding.

Nonspecific binding was less than 2.0% for all assays in U-PLEX Metabolic Group 1 (mouse) using the following calculation.

$$\% \text{ nonspecificity} = \frac{\text{nonspecific signal}}{\text{specific signal}} \times 100$$

Exceptions are noted below:

- The Desghrelin analyte cross-reacts 0.5% with the Ghrelin (active) assay. The Octanoylated Ghrelin analyte cross-reacts with the Ghrelin (total) assay as expected. We do not recommend multiplexing the Ghrelin (total) and Ghrelin (active) assays on the same plate. If measuring the Ghrelin (total) assay, the Ghrelin (active) Calibrator should not be combined with Calibrator 18.
- The GLP-1 (active) analyte cross-reacts with the GLP-1 (total) assay, and the GLP-1 (inactive) analyte cross-reacts with the GLP-1 (total) assay, as expected. We do not recommend multiplexing GLP-1 (active) or GLP-1 (inactive) assays with the GLP-1 (total) assay on the same plate. If measuring the GLP-1 (total) assay, GLP-1 (active) Calibrator should not be combined with Calibrator 18.
- IL-12p70 and IL-23 contain the IL-12p40 subunit. These analytes will cross-react with the IL-12/IL-23p40 assay as expected. We do not recommend multiplexing IL-12p70 or IL-23 assays on the same plate with the IL-12/IL-23p40 assay. If measuring IL-12/IL-23p40 assay, Calibrator 7 should not be combined with Calibrator 5.
- IL-17A and IL-17F analytes cross-react with the IL-17A/F assay as expected. The IL-17A/F analyte cross-reacts with the IL-17A and IL-17F assays as expected. IL-17A/F assay should not be multiplexed with IL-17A or IL-17F assays on the same plate. Any IL-17A/F in samples will be detected as both IL-17A and as IL-17F if these assays are multiplexed. Combining Calibrators 7 and 12 or Calibrators 8 and 12 may affect the quantification of IL-17A or IL-17F in samples.
- MIP-1 $\beta$  calibrator (Calibrator 12) nonspecifically binds to MIP-1 $\alpha$  capture antibody (~0.9%).
- The RANTES detection antibody nonspecifically binds (6%) to the MDC calibrator (Calibrator 16). We do not recommend multiplexing the RANTES assay with the MDC assay on the same plate.
- The Glucagon capture antibody also binds proglucagon which can bind the GLP-1 (total) detection antibody. We do not recommend multiplexing these assays on the same plate.



# Appendix A

## Best Practices

- Equilibrate all assay components to room temperature before use. Mix well before use. Bring plates to room temperature before opening the packet.
- Avoid bubbles at each stage of reagent addition because they can lead to variable results. This is very important when adding Read Buffer (just prior to reading the plate).
- 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. Keep the shaking speed and shaker model consistent for long-term studies.
- Tap the plate on a paper towel after washing to ensure the removal of residual fluid.
- 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 it 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.
- Dispense reagents and wash fluids at the side of the well towards the bottom corner away from the coated spots.
- Remove the plate seal before reading the plate in the instrument.
- Keep time intervals consistent between the addition of Read Buffer and reading the plate to improve interplate precision. Prepare an MSD instrument before 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.
- Avoid cross-contamination between Linkers and antibodies by following the techniques below:
  - Pulse centrifuge the vials to get all of the contents to the bottom of the vial.
  - Open one vial at a time. Close the cap after use.
  - Each Linker vial is color-coded; ensure that each cap and tube have matching colors when opening and closing vials.
  - Use a new filtered pipette for each reagent addition.
- Do not mix components between boxes of multiplex U-PLEX assays except for Stop Solution, Diluents, and Read Buffer. There will be a unique calibrator curve for each plate.

## Working with Partial Plates

A portion of a plate may be used when developing assays. Volumes should be adjusted proportionally when preparing reagents for partial plates.

When running a partial plate, seal the unused sectors to avoid contaminating unused wells. Remove all seals before reading. Partially used plates may be sealed and stored for up to 30 days at 2–8 °C in the original foil pouch with desiccant.

## Working with Multiplate Assays

U-PLEX assays can occupy more than one plate, depending on the number and compatibility of the selected assays. In such cases, each box contains components that are intended to be run together. Do not mix components between boxes, and do not separate the plate from the antibody sets. There will be a unique Calibrator curve for each box.

## Open Spots

### Prepare Conjugated Capture and Detection Antibodies

For assays that are being developed with your own antibody pairs, the capture antibodies (or other suitable capture reagents) must be biotinylated before starting the U-PLEX protocol. Similarly, the detection antibody must be conjugated with SULFO-TAG; however, you may choose to use a SULFO-TAG conjugated secondary detection antibody that is raised against the host of the detection antibody.

**Note:** Do not use a biotinylated detection antibody or SULFO-TAG streptavidin as a method for detection. SULFO-TAG streptavidin will cause high backgrounds, as it will bind to the biotin on the capture antibody.

### Prepare Biotinylated Capture Antibody

The working concentration of biotinylated capture antibody needed to prepare the multiplex coating solution for the U-PLEX Plate is 10 µg/mL. Prepare a stock solution of the biotinylated capture antibody by following the manufacturer's guidelines for the conjugation of an antibody to Sulfo-NHS-LC-Biotin (such as EZ-Link Sulfo-NHS-LC-Biotin [Thermo Fisher Scientific]) or an equivalent product. At least one biotin must be present on the capture antibody for it to be coupled to the U-PLEX Linker. We recommend starting with a biotin challenge ratio of 10 biotins to 1 capture antibody. This challenge ratio typically leads to the conjugation of an average of 2–4 biotins per antibody.

**Note:** Free biotin will interfere with the U-PLEX assay signal. Therefore, after conjugation, it is recommended to purify the biotinylated antibody from the free biotin reagent by using Zeba Desalting Columns.

For long-term storage, perform a buffer exchange to store the final biotinylated antibody in the Conjugate Storage buffer.

### Prepare SULFO-TAG Conjugated Detection Antibody

The optimal concentration of the SULFO-TAG conjugated detection antibody concentration for use in the U-PLEX assay is typically in the range of 0.5–1 µg/mL. Prepare a concentrated stock solution of 100X for each SULFO-TAG conjugated detection antibody by following the guidelines for SULFO-TAG conjugation available at [www.mesoscale.com](http://www.mesoscale.com) (Please refer to the MSD GOLD SULFO-TAG Conjugation Quick Guide or the MSD GOLD SULFO-TAG NHS-Ester Technical Note). We recommend using a 20:1 challenge ratio for SULFO-TAG conjugation of antibodies. This leads to a typical conjugation ratio of 10 SULFO-TAG labels per antibody molecule. Optimization of the SULFO-TAG challenge ratio may be necessary to reduce backgrounds and increase assay signals.

To find out more details on optimizing the SULFO-TAG conjugation of the detection antibody, please refer to the MSD GOLD SULFO-TAG Conjugation Quick Guide or the MSD GOLD SULFO-TAG NHS-Ester technical note available at [www.mesoscale.com](http://www.mesoscale.com).

For long-term storage, purify the SULFO-TAG conjugated antibody to remove the unconjugated SULFO-TAG NHS-Ester. Antibody conjugates are typically stable for at least 1 year in conjugation storage buffer at 2–8 °C. Protect from direct exposure to light.

### **Prepare non-MSD Calibrator**

For assays that are being developed with your antibody pairs, a recombinant protein that is representative of the native protein can be used for the calibration curve. A good starting concentration is 10 ng/mL for the high calibrator and 0.001 ng/mL for the low calibrator. We recommend testing an 8-point titration curve and optimizing the calibrator diluent if required. Guidance on using recombinant protein calibrators can be found in the Development Pack Product Insert at [www.mesoscale.com/U-PLEX-documents](http://www.mesoscale.com/U-PLEX-documents).

# Appendix B

## Components for 384-well Assays

*Table 8. Reagents that are supplied with all U-PLEX Metabolic Group 1 (mouse) 384-well Assays*

Reagent	Storage	Catalog No.	Size	Quantity Supplied		Description
				5 Plates	25 Plates	
Diluent 13	≤-10 °C	R56BB-3	50 mL	2 bottles	10 bottles	Diluent for samples and calibrators
Diluent 11	≤-10 °C	R55BA-3	50 mL	2 bottles	10 bottles	Diluent for detection antibody
Aprotinin, 200,000 KIU/mL	2–8 °C	R93AD-2	250 µL	2 vials	10 vials	200X concentrated solution
Stop Solution	2–8 °C	R50AO-1	40 mL	2 bottles	10 bottles	Biotin-containing buffer to stop Linker-antibody coupling reaction
MSD GOLD Read Buffer B	RT	R60AM-2	90 mL	1 bottle	5 bottles	Buffer to catalyze the electrochemiluminescent reaction

## Reagent Preparation for 384-well Plates

**Important:** Upon the first thaw, aliquot diluents into suitably sized aliquots before refreezing.

### STEP 1: Create Individual U-PLEX Linker-Coupled Antibody Solutions

A different Linker is used for each unique biotinylated antibody. Below are the steps to couple linkers with capture antibodies.

- Couple each biotinylated capture antibody to a unique Linker and record the antibody identity next to the Linker number on the Spot Map (blank Spot Maps are provided on page 23).
- Add 200 µL of each biotinylated antibody to 300 µL of the assigned Linker. Mix by vortexing. Incubate at room temperature for 30 minutes. Do not shake.

#### Notes:

- Each Linker vial has a matching colored cap and label.
- To remove liquid from the cap, briefly centrifuge the Linker vial and open the cap gently.
- Open one Linker at a time and close its cap as soon as you are done using it. Take precautions to avoid reagent contamination.
- For studies using multiple plates of the same assay, it is recommended that the same Linker be coupled with the same antibody for the duration of the study.

- Add 200 µL of Stop Solution, then mix by vortexing. Incubate at room temperature for 30 minutes.

**Note:** At the end of step 1, each individual U-PLEX Linker-coupled antibody solution is at 10X the coating concentration and can be stored at 2–8 °C. Do not store for more than 7 days.

Adjust the volumes for multiple plates (see Table 5). The volumetric ratio of Linker:antibody:Stop Solution is 3:2:2.

### STEP 2: Prepare the Multiplex Coating Solution for 384-well Plates

- Combine 600 µL of each U-PLEX Linker-coupled antibody solution into a single tube and mix by vortexing. Up to 4 U-PLEX Linker-coupled antibodies can be pooled. Do not combine U-PLEX Linker-coupled antibody solutions that share the same Linker.
- Bring the solution up to 12 mL by mixing it with Stop Solution. Mix by vortexing.

**Note:** At the end of Step 2, the U-PLEX multiplex coating solution is at 0.5X and can be stored at 2–8 °C. Do not store for more than 7 days.

### STEP 3: Coat U-PLEX 384-well Plates

- Wash the plate 3 times with 80 µL/well of 1X Wash Buffer.
- Add 25 µL of the 0.5X multiplex coating solution to each well. Seal the plate with an adhesive plate seal and shake for 4 hours at room temperature.
- Wash the plate 3 times with 80 µL/well of 1X Wash Buffer.

**Note:** The plate is now coated and ready for use. Sealed plates may be stored in the original pouch with desiccant for up to 7 days at 2–8 °C.

## Prepare Metabolic Assay Working Solution

This procedure is used for preparing the calibrator, controls, and diluting the samples.

For one plate, combine the following in a 15 mL tube:

- 13.93 mL of Diluent 13
- 70 µL of aprotinin

### Notes:

- The addition of a DPP-IV inhibitor (such as diprotin A, not provided) to a final concentration of 0.1 mM in the Metabolic Assay Working Solution is strongly recommended. A DPP-IV inhibitor will aid in limiting the enzymatic action of DPP-IV present in serum/plasma and provides the most accurate measurement of some metabolic analytes.
- Ghrelin (active) assay only: Addition of Halt Protease Inhibitor Cocktail, EDTA-Free (Thermo Fisher Scientific, Catalog No. 87785) to a final concentration of 1X in the Metabolic Assay Working Solution is strongly recommended. The presence of Halt inhibitor may affect sample quantitation in other assays.

**Important:** Protease inhibitors should be added before use. The Metabolic Assay Working Solution should be kept on ice. Do not freeze the Metabolic Assay Working Solution for later use.

## Prepare Detection Antibody Solution

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

For one plate, combine:

- 60 µL of each 100X detection antibody
- Diluent 45 to bring the final volume to 12 mL

# Assay Protocol (384-well plates)

## STEP 1: Add Samples and Calibrators

- Add 25  $\mu\text{L}$  of the prepared Calibrator Standard or sample to each well of a prepared plate. Seal the plate with an adhesive plate seal. Incubate at room temperature with shaking for 4 hours.

## STEP 2: Wash and Add Detection Antibody Solution

- Wash the plate 3 times with 80  $\mu\text{L}$ /well of 1X MSD Wash Buffer.
- Add 25  $\mu\text{L}$  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 80  $\mu\text{L}$ /well of 1X MSD Wash Buffer.
- Add 40  $\mu\text{L}$  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 suggestion below may be useful for simplifying the protocol.

- Alternate Protocol, Shortened Incubations:** Some assays in 384-well plates may achieve acceptable performance with shorter incubations. Consider incubating samples in the plate for 2 hours.

# Spot Maps



- 1 \_\_\_\_\_
- 2 \_\_\_\_\_
- 3 \_\_\_\_\_
- 4 \_\_\_\_\_
- 5 \_\_\_\_\_
- 6 \_\_\_\_\_
- 7 \_\_\_\_\_
- 8 \_\_\_\_\_
- 9 \_\_\_\_\_
- 10 \_\_\_\_\_



- 1 \_\_\_\_\_
- 3 \_\_\_\_\_
- 9 \_\_\_\_\_
- 10 \_\_\_\_\_

Figure 5. Spot maps. (top) 96-well plate, (bottom) 384-well plate.

# Plate Diagrams

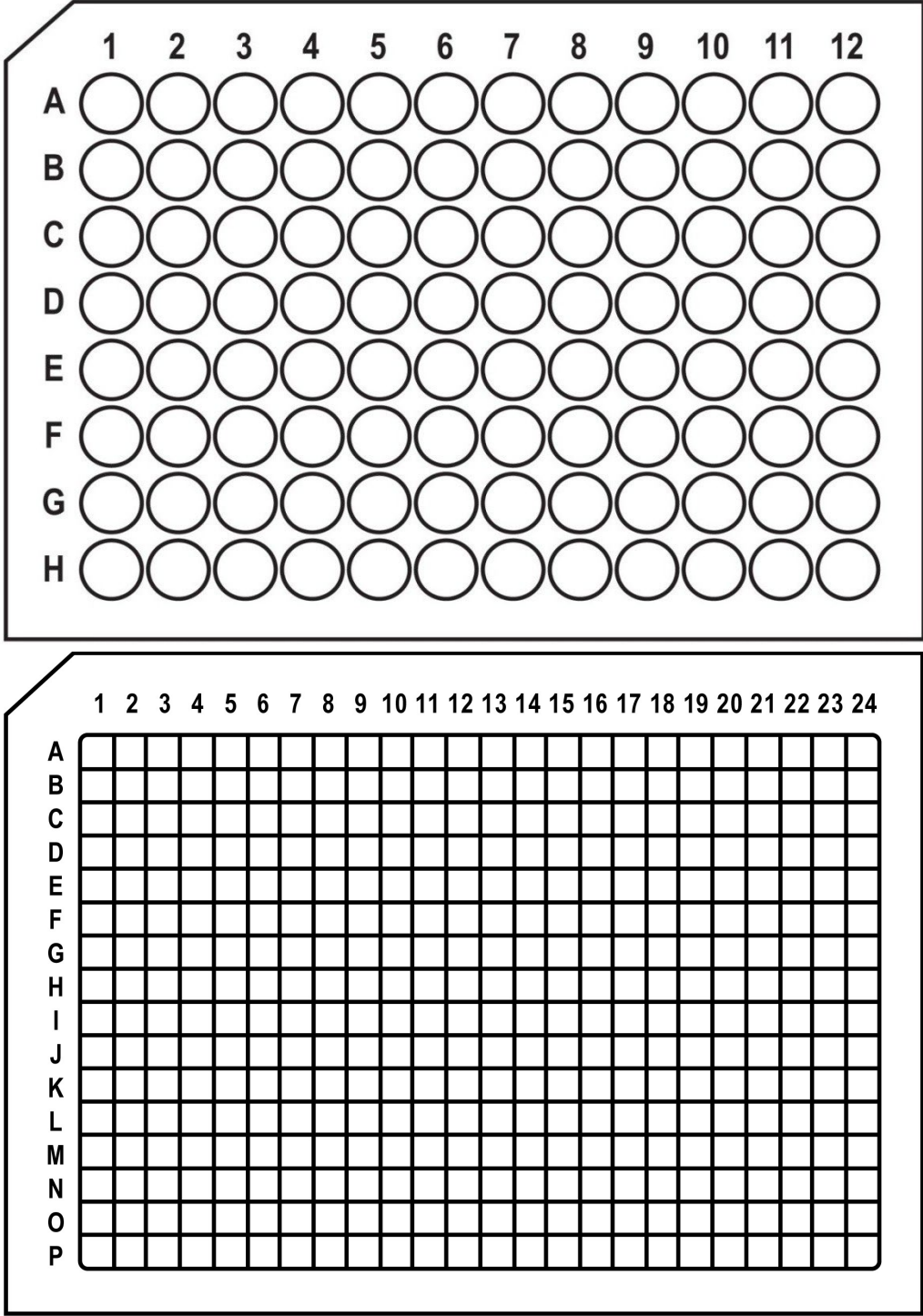


Figure 6. Plate diagrams; similar plate layouts can be created in Excel and in the DISCOVERY WORKBENCH® software.