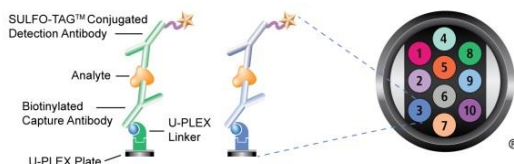


# Quick Guide

## U-PLEX® Biomarker Group 1 (Human) - Multiplex Assays

The MSD® U-PLEX platform provides a rapid and convenient method for creating your own multiplex assays. Using two simple tools – a 10-spot U-PLEX plate and unique Linkers, you can build custom multiplex panels for any combination of analytes. The U-PLEX assay menu is organized into groups which include a broad menu of analytes assembled together by species, abundance in serum and plasma samples, analytical compatibility, and expected use. As many as 10 U-PLEX assays from an assay group may be multiplexed on each plate for simultaneous quantification. For ultimate flexibility, custom panels can be created from a selection of MSD U-PLEX assays, your own antibodies, or a combination of both.



Read the U-PLEX Biomarker Group 1 (Human) Multiplex Assays product insert available at [www.mesoscale.com/U-PLEX-documents](http://www.mesoscale.com/U-PLEX-documents) for detailed instructions.

### Components

Reagent	Storage	Description
<b>10-spot 96-well U-PLEX Plate</b>	2–8°C	Foil sealed, with desiccant. The appropriate number of spots is activated to match the number of assays ordered.
<b>Linkers (1–10)</b>	2–8°C	The appropriate number of Linkers is provided to match the number of assays ordered.
<b>U-PLEX Antibody Sets</b>	2–8°C	Sets containing a biotinylated capture antibody and SULFO-TAG™ conjugated detection antibody.
<b>Calibrators</b>	2–8°C	Multi-analyte blends, each containing multiple recombinant human proteins lyophilized in a buffered diluent. Individual analyte concentration is provided in the lot-specific certificate of analysis.
<b>Diluent 43</b>	≤–10°C	Diluent for samples and Calibrators; contains serum, blockers, and preservatives.
<b>Diluent 3</b>	≤–10°C	Diluent for detection antibody; contains protein, blockers, and preservatives.
<b>Stop Solution</b>	2–8°C	Biotin-containing buffer to stop Linker-antibody coupling reaction.
<b>Read Buffer T (4X)</b>	RT	Buffer to catalyze the electrochemiluminescence reaction. Dilute to 2X and use at room temperature.

### Additional Materials

MSD Wash Buffer (20X, 100 mL, catalog # R61AA-1) or Phosphate-buffered saline (PBS) plus 0.05% Tween-20 (PBS-T) for plate washing. For one plate, combine 15 mL of MSD Wash Buffer (20X) with 285 mL of deionized water.

For Research Use Only. Not for use in diagnostic procedures.

# Reagent Preparation

**Note:** Bring all reagents to room temperature.

Please consult the U-PLEX Biomarker Group 1 (Human) Multiplex Assays product insert available at [www.mesoscale.com/U-PLEX-documents](http://www.mesoscale.com/U-PLEX-documents) for detailed instructions.

## Prepare the U-PLEX Plate

### Step 1 Create the Individual U-PLEX-Coupled Antibody Solutions

The biotinylated capture antibody is provided at a ready-to-use concentration. Couple an individual biotinylated antibody to a unique Linker and record the antibody identity next to the Linker number on the Spot Map below.

- ❑ Add 200  $\mu$ L of each biotinylated antibody to 300  $\mu$ L of the assigned Linker. For 1-plate assay packs, the biotinylated antibody may be added directly to the Linker vial. A different Linker must be used for each biotinylated antibody. Mix by vortexing. Incubate at room temperature for 30 minutes.
- ❑ Add 200  $\mu$ L of Stop Solution. Mix by vortexing. Incubate at room temperature for 30 minutes. Each individual U-PLEX-coupled antibody solution is at 10X the coating concentration and can be stored at 2-8°C for up to 7 days.

### Step 2 Prepare the Multiplex Coating Solution

- ❑ Combine 600  $\mu$ L of each U-PLEX-coupled antibody solution into a single tube and vortex. Up to 10 U-PLEX-coupled antibodies can be pooled.  
*Do not combine U-PLEX-coupled antibody solutions that share the same Linker.*
- ❑ When combining fewer than 10 antibodies, bring the solution up to 6 mL with Stop Solution to result in a final 1X concentration. Mix by vortexing. The 1X multiplex coating solution can be stored at 2-8°C for up to 7 days.



For Research Use Only. Not for use in diagnostic procedures.

# Reagent Preparation

## Step 3 Coat the U-PLEX Plate

- ❑ Add 50  $\mu\text{L}$  of 1X multiplex coating solution to each well. Seal the plate with an adhesive plate seal and incubate with shaking at room temperature for 1 hour or at 2–8°C overnight.
- ❑ Wash the plate 3 times with at least 150  $\mu\text{L}$ /well of 1X MSD Wash Buffer or PBS-T (PBS plus 0.05% Tween-20). The plate is now coated and ready for use. Plates may be stored in the original pouch with desiccant and sealed, for up to 7 days at 2–8°C.

## Prepare Calibrator Standards

Depending on the assays ordered, you may receive one or more multi-analyte Calibrator vials with your order. Individual analyte concentrations are provided in the lot-specific certificates of analysis (COA) that are shipped with the product.

To prepare 7 Calibrator Standard solutions plus a zero Calibrator Standard for up to 6 replicates:

- ❑ Reconstitute each vial of Calibrator by adding 250  $\mu\text{L}$  of Diluent 43 to the glass vial and inverting 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.

If you have more than one Calibrator, create a Calibrator Blend. Combine 50  $\mu\text{L}$  of each reconstituted Calibrator in a clean polypropylene tube and bring the total volume up to 250  $\mu\text{L}$  in Diluent 43.

**Note:** We recommend that reconstituted Calibrators are used immediately. If storage is necessary, divide into 60  $\mu\text{L}$  aliquots and store immediately at  $\leq -70^\circ\text{C}$ .

- ❑ Prepare Calibrator Standard 1 by adding 50  $\mu\text{L}$  of the Calibrator (or Calibrator Blend) to Diluent 43 to make a final volume of 250  $\mu\text{L}$ . Mix by vortexing.
- ❑ Prepare Calibrator Standard 2 by adding 75  $\mu\text{L}$  of Calibrator Standard 1 to 225  $\mu\text{L}$  of Diluent 43. 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 Diluent 43 as Calibrator Standard 8 (zero Calibrator).

## Dilute Samples

- ❑ Depending on the sample set under investigation, a dilution may be necessary. Diluent 43 (catalog # R50AG-1) may be used for sample dilution.

**Note:** TGF- $\beta$  samples require an acid treatment for activation prior to use:

1. Add 20  $\mu\text{L}$  of 1 M HCl per 100  $\mu\text{L}$  of sample volume. Vortex briefly.
2. Incubate sample for 10 min at room temperature.
3. Neutralize the sample by adding 14  $\mu\text{L}$  of 1.2M NaOH/0.5M HEPES per 100  $\mu\text{L}$  of sample volume. Vortex briefly. Samples are ready to use. Use immediately.

## Prepare Detection Antibody Solution

- ❑ The detection antibody is provided as a 100X stock solution. For one plate, combine 60  $\mu\text{L}$  of each supplied 100X detection antibody. Add Diluent 3 to bring the final volume to 6 mL.

For Research Use Only. Not for use in diagnostic procedures.

# Reagent Preparation/Assay Protocol

## Prepare 2X Read Buffer T

- ❑ For one plate, combine 10 mL of Read Buffer T (4X) and 10 mL of deionized water. You may keep excess diluted Read Buffer T in a tightly sealed container at room temperature for up to one month.

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

## Step 1 Add Samples and Calibrators

- ❑ Add 25  $\mu$ L of Diluent 43 to each well. Tap the plate gently on all sides.
- ❑ Add 25  $\mu$ 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  $\mu$ L/well of 1X MSD Wash Buffer or PBS-T.
- ❑ Add 50  $\mu$ 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  $\mu$ L/well of 1X MSD Wash Buffer or PBS-T.
- ❑ Add 150  $\mu$ L of 2X Read Buffer T to each well. Analyze the plate on an MSD instrument. Incubation in Read Buffer T is not required before reading the plate.

### 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)

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, MSD GOLD, DISCOVERY WORKBENCH, MULTI-ARRAY, MULTI-SPOT, QUICKPLEX, SECTOR, SECTOR PR, SECTOR HTS, SULFO-TAG, U-PLEX, S-PLEX, V-PLEX, STREPTAVIDIN GOLD, MESO, [www.mesoscale.com](http://www.mesoscale.com), SMALL SPOT (design), 96 WELL 1, 4, 7, 9, & 10-SPOT (designs), 384 WELL 1 & 4-SPOT (designs), MSD (design), U-PLEX (design), S-PLEX (design), V-PLEX (design), It's All About U, and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.

©2015-2016 Meso Scale Diagnostics, LLC. All rights reserved.

For Research Use Only. Not for use in diagnostic procedures.

MK-DS-232-v3-2016Nov