

# MSD<sup>®</sup> 96-Well MULTI-ARRAY<sup>®</sup> Human VEGF Assay

The following assay protocol has been optimized for analysis of Human Vascular Endothelial Growth Factor (VEGF) in human serum and plasma samples.

Storage

## MSD Materials

<input type="checkbox"/> Read Buffer T (4X), with surfactant	RT
<input type="checkbox"/> Blocker C	2-8 °C
<input type="checkbox"/> MULTI-ARRAY 96-well Small Spot VEGF Plate(s)	2-8 °C
<input type="checkbox"/> SULFO-TAG <sup>™</sup> Anti-hVEGF Antibody (100X) <sup>1</sup>	2-8 °C
<input type="checkbox"/> Diluent 7	≤-10 °C
<input type="checkbox"/> Diluent 8	≤-10 °C
<input type="checkbox"/> Diluent 9	≤-10 °C
<input type="checkbox"/> Human VEGF Calibrator (1µg/mL)	≤-70 °C

## Other Materials & Equipment (not supplied)

- Deionized water for diluting Read Buffer
- Phosphate Buffered Saline + 0.05% Tween-20 (PBS-T) for plate washing
- Adhesive plate seals
- Microtiter plate shaker
- Plate washer or other efficient multi-channel pipetting equipment for washing 96 well plates
- Appropriate liquid handling equipment for desired throughput that must accurately dispense 25 and 150 µL into a 96-well microplate

Meso Scale Discovery, Meso Scale Diagnostics, www.mesoscale.com, MSD, MSD (design), Discovery Workbench, Quickplex, Multi-Array, Multi-Spot, Sulfo-Tag and Sector are trademarks of Meso Scale Diagnostics, LLC.

© 2009 Meso Scale Discovery a division of Meso Scale Diagnostics, LLC. All rights reserved.

<sup>1</sup> SULFO-TAG labeled detection antibodies may be light-sensitive, so they should be stored in the dark.

FOR RESEARCH USE ONLY.  
NOT FOR USE IN DIAGNOSTIC OR THERAPEUTIC PROCEDURES.



## Notes:

Read the entire detailed instructions before beginning work.

The assay protocol was optimized for human serum samples. For significantly different sample matrices, it is recommended to use a Calibrator diluent that is similar to the sample matrix (e.g. lysis buffer + carrier protein).

## Protocol at a Glance

The protocol can be completed in approximately **5.5 hours** if each reagent is prepared during the preceding incubation. This time can be reduced to **4.5 hours** if the blocking reagent is added the night before.

1. Block plates for 1-2 hours at room temperature (alternatively block plates overnight at 4 °C).
2. Wash.
3. Add Diluent 7 and Calibrator and/or sample and incubate for 2 hours with shaking.
4. Wash.
5. Add Detection Antibody Reagent and incubate for 2 hours with shaking.
6. Wash.
7. Add Read Buffer and read immediately.

## Preparation Instructions

### Prepare Calibrator dilutions:

1. Determine how many Calibrator levels and replicates will be run. Each well will require 25  $\mu\text{L}$  of Calibrator. Thaw one vial of VEGF Calibrator stock solution and prepare the required Calibrator dilution series using the stock solution and Diluent 9.
  - A recommended Calibrator dilution procedure is listed below for up to 4 replicates of 7 Calibrator concentrations spanning a wide range, plus 1 zero-Calibrator point.
  - Prepare 200  $\mu\text{L}$  of a high Calibrator containing 100 ng/mL VEGF by combining 20  $\mu\text{L}$  of VEGF stock solution at 1  $\mu\text{g}/\text{mL}$  with 180  $\mu\text{L}$  of Diluent 9.
  - Prepare 6 additional 1:4 serial dilutions, beginning with the high Calibrator, by adding 50  $\mu\text{L}$  of the Calibrator to 150  $\mu\text{L}$  Diluent 9.
  - This will create 7 Calibrators with 100000, 25000, 6250, 1563, 391, 98, and 24 pg/mL of VEGF.
  - The recommended 8<sup>th</sup> dilution is Diluent 9 alone (e.g. zero Calibrator).
- ❖ Once the expected range of sample concentrations is known, the Calibrator concentrations can be adjusted appropriately.
2. Calibrators are stable at room temperature for a few hours.
3. The human VEGF Calibrator has been anchored and referenced to international standards. The table below summarizes the reference information.



**Notes:**

Analyte	WHO Standard Reference Number	WHO Standard Units / $\mu\text{g}$	MSD Calibrator $1\mu\text{g}$ = WHO Units	WHO Units
h VEGF	01/424	n/a	0.5	$\mu\text{g}$
h VEGF	02/286	1,000	360	U

\*\* MSD VEGF Calibrator previously used in Human VEGF Kits and Human Hypoxia Kits was anchored to WHO Standard Reference 01/424 with  $1\mu\text{g}$  of MSD Calibrator =  $1\mu\text{g}$  of WHO Standard

**Prepare Detection Antibody Reagent:**

1. Each well requires  $25\mu\text{L}$  of Detection Antibody Reagent. Prepare  $3\text{ mL}$  per plate.
2. In a  $15\text{ mL}$  tube combine:
  - a.  $2.97\text{ mL}$  Diluent 8
  - b.  $30\mu\text{L}$  of 100X SULFO-TAG Anti-hVEGF Antibody (final concentration: 1X)

**Dilute Read Buffer:**

1. Determine total number of wells in the experiment. Each well will receive  $150\mu\text{L}$  of Read Buffer T. Prepare an extra 20%.
2. Dilute 4X Read Buffer T to 1X with deionized water.
3. Diluted Read Buffer may be stored at room temperature for later use.

*Diluted Read Buffer may be kept in a tightly sealed container at room temperature for later use.*

**Assay Protocol**

Begin with a MULTI-ARRAY 96-well Small Spot VEGF plate.

No pre-treatment is necessary.

1. Add  $150\mu\text{L}$ /well of Blocker C and incubate at room temperature for 1-2 hours or overnight at  $4^\circ\text{C}$ .
2. Wash plates 3 times with Phosphate Buffered Saline + 0.05% Tween-20 (PBS-T).
3. Add  $25\mu\text{L}$ /well of Diluent 7.  
Add  $25\mu\text{L}$ /well Calibrator or sample and incubate at room temperature with shaking for 2 hours.
4. Wash plates 3 times with PBS-T.
5. Add  $25\mu\text{L}$ /well Detection Antibody Reagent and incubate at room temperature with shaking for 2 hours.
6. Wash plates 3 times with PBS-T.
7. Prepare SECTOR<sup>®</sup> Imager so that plate can be read immediately after Read Buffer addition.
8. Add  $150\mu\text{L}$ /well 1X Read Buffer T.
9. Analyze immediately with SECTOR Imager.

*Plates may also be blocked overnight at  $4^\circ\text{C}$  and stored for up to a week with blocker.*

*Shaking a 96-well MSD MULTI-ARRAY<sup>®</sup> or MULTI-SPOT plate typically accelerates capture at the working electrode.*

*Bubbles in the Read Buffer will interfere with reliable imaging of the plate.*

