

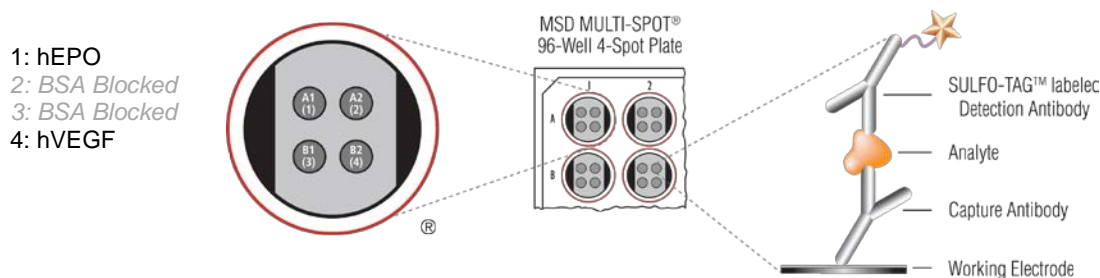
# MSD<sup>®</sup> 96-Well MULTI-SPOT<sup>®</sup> Human Hypoxia Serum/Plasma Assay

The following assay protocol has been optimized for quantifying erythropoietin (EPO) and 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-SPOT 96-well 4 Spot Human Hypoxia Plate(s)	2-8 °C
<input type="checkbox"/> SULFO-TAG <sup>™</sup> Anti-hEPO Antibody (100X) <sup>1</sup>	2-8 °C
<input type="checkbox"/> SULFO-TAG 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"/> Human VEGF Calibrator (1 µg/mL)	≤ -70 °C
<input type="checkbox"/> Human EPO Calibrator (20 IU/mL)	≤ -70 °C



The SECTOR<sup>®</sup> Imager data file will identify spots according to their well location, not by the coated capture antibody name.

## 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, which can be obtained from MSD Customer Service.

<sup>1</sup> SULFO-TAG conjugated detection antibodies should be stored in the dark.

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



## Other Materials & Equipment (not supplied)

**Notes:**

- ❑ Deionized water for diluting concentrated buffers
- ❑ Phosphate buffered saline with 0.05% Tween-20 (PBS-T) for plate washing
- ❑ Automatic plate washer or other efficient multi-channel pipetting equipment for washing 96-well plates
- ❑ Microtiter plate shaker
- ❑ Adhesive plate seals
- ❑ Pipettes that accurately dispense 25, 50, and 150  $\mu\text{L}$  into a 96-well microplate

*Read the entire detailed instructions before beginning work.*

## Protocol at a Glance

The following protocol describes a recommended assay format. 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 overnight.

- Step 1.** Add blocking solution, incubate 1 hour, wash.  
(Alternatively, block plates overnight at 4 °C.)
- Step 2.** Add 25  $\mu\text{L}$  of Diluent 7.  
Add 25  $\mu\text{L}$  of samples or calibrator, incubate 2 hours, wash.
- Step 3.** Add 25  $\mu\text{L}$  of detection antibody, incubate 2 hours, wash.
- Step 4.** Add 150  $\mu\text{L}$  of read buffer, read plate, and analyze data.

## Detailed Instructions

### Prepare Calibrator dilutions:

1. Determine the number of calibrator levels and replicates that will be run. Each well will require 25  $\mu\text{L}$  of calibrator. Thaw one vial of EPO and VEGF calibrator stock solutions and prepare the required calibrator dilution series using the stock solutions and Diluent 7.
  - Our 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 combined calibrator containing 100 ng/mL VEGF and 10 IU/mL of EPO by combining 20  $\mu\text{L}$  of VEGF stock solution at 1  $\mu\text{g/mL}$  and 100  $\mu\text{L}$  of the 20 IU/mL EPO stock solution with 80  $\mu\text{L}$  of Diluent 7.*
  - *Prepare 6 additional 1:4 serial dilutions, beginning with the high combined calibrator, by adding 50  $\mu\text{L}$  of the calibrator to 150  $\mu\text{L}$  Diluent 7.*



**Notes:**

- This will create 7 calibrators with 100000, 25000, 6250, 1563, 391, 98, 24 pg/mL of VEGF, and 10000, 2500, 625, 156, 39, 9.8, 2.4 mIU/mL of EPO.
  - The recommended 8<sup>th</sup> dilution is Diluent 7 alone (i.e., 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 following calibrators have been anchored and referenced to international standards when available. The table below summarizes the reference information.

Analyte	WHO Standard Reference Number	WHO Standard Units/ $\mu$ g	MSD Calibrator 1 $\mu$ g = WHO Units	WHO Units
hEPO	88/574	127	130	IU
hVEGF	01/424	n/a	0.5	$\mu$ g
hVEGF	02/286	1,000	360	U

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

**Prepare Detection Antibody Reagent:**

1. Each well will require 25  $\mu$ L of detection antibody reagent. Prepare 3 mL per plate.
2. In a 15 mL tube combine:
  - a. 2.94 mL Diluent 8
  - b. 30  $\mu$ L of 100X SULFO-TAG Anti-hEPO Antibody (final concentration: 1X)
  - c. 30  $\mu$ L of 100X SULFO-TAG Anti-hVEGF Antibody (final concentration: 1X)

**Prepare Diluted Read Buffer:**

- In a 50 mL tube combine (per plate):
- a. 5 mL 4X Read Buffer T
  - b. 15 mL deionized water

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



## Assay Protocol

Begin with a MULTI-SPOT 96-well 4 Spot Human Hypoxia plate.  
No pre-treatment is necessary.

1. Add 150  $\mu$ L/well of blocking solution C and incubate at room temperature for 1 hour or overnight at 2-8  $^{\circ}$ C.
2. Wash plates 3 times with phosphate buffered saline + 0.05% Tween-20 (PBS-T).
3. Add 25  $\mu$ L/well of Diluent 7.
4. Add 25  $\mu$ L/well calibrator or sample and incubate at room temperature with shaking for 2 hours.
5. Wash plates 3 times with PBS-T.
6. Add 25  $\mu$ L/well detection antibody reagent and incubate at room temperature with shaking for 2 hours.
7. Wash plates 3 times with PBS-T.
8. Prepare SECTOR Imager such that plate can be read immediately after read buffer addition.
9. Add 150  $\mu$ L/well 1X Read Buffer T.
10. Analyze immediately with SECTOR Imager.

*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.*

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

©2014 Meso Scale Diagnostics, LLC. All rights reserved.

