

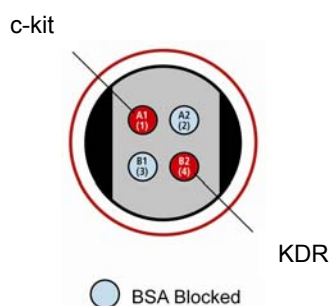
# MSD<sup>®</sup> 96-Well MULTI-SPOT<sup>®</sup> Human Growth Factor panel II Assay

The following assay protocol has been optimized for analysis of kinase domain insert (KDR) and the kit ligand receptor (c-Kit) in human serum and plasma samples.

Storage

## MSD Materials

<input type="checkbox"/> Read Buffer T (4X) with surfactant	RT
<input type="checkbox"/> Blocker A Kit	RT
<input type="checkbox"/> MULTI-SPOT 96-well 4 Spot Human Growth Factor II Plate(s)	2-8 °C
<input type="checkbox"/> SULFO-TAG <sup>™</sup> Anti-Human Growth Factor II Detection Antibody Blend (100X) <sup>1</sup>	2-8 °C
<input type="checkbox"/> Diluent 10	≤-10 °C
<input type="checkbox"/> Diluent 11	≤-10 °C
<input type="checkbox"/> Human Growth Factor II High Calibrator Blend (150 ng/mL c-Kit, 15 ng/mL KDR)	≤-70 °C



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

<sup>1</sup> Some 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:

### *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 washing equipment 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  $\mu$ L into a 96-well micro plate

*Read the entire detailed instructions before beginning work.*

### *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.

- Step 1.** Add Blocking Solution, incubate 1-2 hours, wash. (alternatively, block plates overnight at 4 °C).
- Step 2.** Add 50  $\mu$ L of Calibrator or Samples (diluted 50X), incubate 2 hours, wash.
- Step 3.** Add 25  $\mu$ L of Detection Antibody, incubate 2 hours, wash.
- Step 4.** Add 150  $\mu$ L of Read Buffer, read plate and analyze data.

### *Preparation Instructions*

#### **Prepare Blocker A:**

Prepare Blocker A Solution using the instructions provided with the Blocker A kit.

#### **Thaw Diluent:**

Thaw Diluent 10. Vortex briefly. Diluent is stable at 4°C for one week.

#### **Prepare Calibrator and Sample dilutions:**

1. Determine how many Calibrator levels and replicates will be tested in the experiment. Each well will require 50  $\mu$ L of Calibrator or 50  $\mu$ L of diluted sample per well. Thaw Diluent 10 and one vial of High Calibrator Blend. Vortex briefly. Prepare the required Calibrator dilution series using Diluent 10.



## Notes:

2. A recommended Calibrator dilution procedure is listed below for 3 replicates of 6 Calibrator concentrations spanning a wide range, plus 1 zero-Calibrator point.
  - *Prepare 1:3 serial dilutions beginning with the High Calibrator Blend, by adding 100  $\mu$ L of High Calibrator Blend to 200  $\mu$ L of Diluent 10. Prepare 6 serial dilutions. The first Calibrator will be High Calibrator Blend stock and the 8<sup>th</sup> Calibrator should be Diluent 10 alone.*
  - *This will create seven Calibrators with 15 ng/mL, 5 ng/mL, 1.67 ng/mL, 0.556 ng/mL, 0.185 ng/mL, 0.062 ng/mL, 0.021 ng/mL, and 0 ng/mL. The c-Kit concentrations will be 10X higher.*
  - *Since the sample will be diluted 1:50, the concentrations of the Calibrators need to be multiplied by 50 if samples are read directly from the calibration curve. Thus, the dilution-corrected High KDR Calibrator is 750 ng/mL, and the dilution-corrected High c-Kit Calibrator is 7500 ng/mL.*
3. Calibrators are stable at room temperature for a few hours. The High Calibrator Blend stock solution is stable for one day at 4 °C or one additional freeze-thaw. Diluent 10 is stable for one week at 4 °C.
4. Dilute samples 1:50 in Diluent 10. Each well will require 50  $\mu$ L of diluted sample.

### Prepare the 1X Detection Antibody Solution:

1. In a 15 mL tube combine:
  - ☐ 30  $\mu$ L of 100X SULFO-TAG Detection Antibody Blend
  - ☐ 2.97 mL of Diluent 11
2. This will yield 3 mL of diluted Detection Antibody Solution at the working concentration with sufficient volume for one plate.

*Detection Antibody Solution is stable at RT for a few hours.*

### Dilute Read Buffer:

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

*Diluted Read Buffer may be stored at room temperature for later use.*



## ***Assay Protocol***

Begin with a MULTI-SPOT 96-well 4 Spot Human Growth Factor II Plate.  
No pre-treatment is necessary.

1. Add 150  $\mu$ L/well of Blocker A Solution and incubate at room temperature for 1-2 hours or overnight at 4 °C.
2. Wash plates 3 times with phosphate buffered saline with 0.05% Tween-20 (PBS-T).
3. Add 50  $\mu$ L/well of Calibrator or diluted sample and incubate at room temperature with shaking for 2 hours.
4. Wash plates 3 times with PBS-T.
5. Add 25 $\mu$ L/well of 1X Detection Antibody Solution and incubate at room temperature with shaking for 2 hours.
6. Wash plates 3 times with PBS-T.
7. Prepare SECTOR Imager so that plate can be read immediately after Read Buffer addition.
8. Add 150  $\mu$ L/well 1X Read Buffer T.
9. Analyze immediately with SECTOR Imager.

*Avoid bubbles while adding the Read Buffer; it will interfere with accurate reading of the plate.*

