

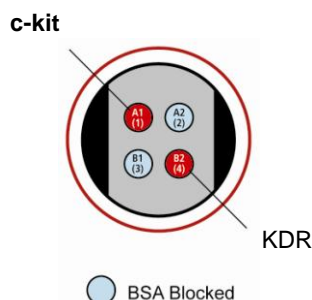
# MSD<sup>®</sup> 96-Well MULTI-ARRAY<sup>®</sup> Human c-Kit Assay

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

Storage

## MSD Materials

❑ Read Buffer T (4X) with surfactant	RT
❑ Blocker A Kit	RT
❑ MULTI-SPOT <sup>®</sup> 96-well 4 Spot Human Growth Factor II Plate(s)	2-8 °C
❑ SULFO-TAG <sup>™</sup> Anti-hc-kit Antibody (100X) <sup>1</sup>	2-8 °C
❑ Diluent 10	≤-10 °C
❑ Diluent 11	≤-10 °C
❑ 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.



## ***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 150µL into a 96-well micro plate

## ***Protocol at a Glance***

*Read the entire detailed instructions before beginning work.*

**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**. All reagents can be prepared hours ahead of time if desired.

- Step 1.** Add Blocking Solution, incubate 1-2 hours, wash. (alternatively, block plates overnight at 4 °C).
- Step 2.** Add 50 µL of Calibrator or diluted Samples (diluted 50X in Diluent 10), incubate 2 hours, wash.
- Step 3.** Add 25 µL of Detection Antibody, incubate 2 hours, wash.
- Step 4.** Add 150 µL of Read Buffer, read plate and analyze data.

## ***Preparation Instructions***

### **Prepare Blocker A:**

1. Prepare Blocker A Solution using the instructions provided with the Blocker A kit.
2. 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 µL of Calibrator or 50 µL of diluted sample per well. Thaw a vial of Diluent 10 and one vial of High Calibrator. Vortex briefly. Prepare the required Calibrator dilution series using Diluent 10.
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. *Note: The curve should be adjusted as necessary to provide the proper range for test samples.*

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



## Notes:

- *Prepare 1:3 serial dilutions beginning with the High Calibrator, by adding 100  $\mu$ L of High Calibrator to 200  $\mu$ L of Diluent 10. Prepare 6 serial dilutions. The first Calibrator will be High Calibrator stock and the 8<sup>th</sup> Calibrator should be Diluent 10 alone.*
  - *This will create seven Calibrators with 150 ng/mL, 50 ng/mL, 16.7 ng/mL, 5.56 ng/mL, 1.85 ng/mL, 0.62 ng/mL, 0.21 ng/mL, and 0 ng/mL c-Kit.*
  - *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 c-Kit Calibrator is 7500 ng/mL.*
3. Calibrators are stable at room temperature for a few hours. The High Calibrator 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 Detection Antibody Reagent:

1. Each well requires 25  $\mu$ L of Detection Antibody Reagent. Prepare 3 mL per plate.
2. In a 15 mL tube combine:
  - a. 2.97 mL Diluent 11
  - b. 30  $\mu$ L of 100X SULFO-TAG Anti-hc-Kit Antibody (final concentration: 1X)

### Dilute Read Buffer:

In a 50 mL tube combine (per plate):

1. 5 mL 4X Read Buffer T
2. 15 mL deionized water

*Diluted Read Buffer 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 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 + 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 Detection Antibody Reagent 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.*

