

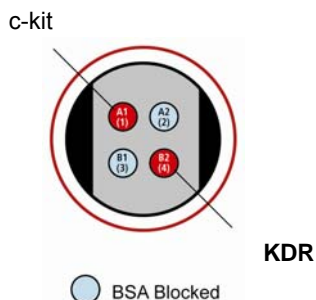
MSD[®] 96-Well MULTI-ARRAY[®] Human KDR Assay

The following assay protocol has been optimized for analysis of kinase domain insert (KDR) 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 [™] Anti-hKDR Antibody (100X) ¹ | 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[®] Imager data file will identify spots according to their well location, not by the coated capture antibody name.

¹ 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 μ 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**. 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 μ L of Calibrator or diluted Samples (diluted 50X in Diluent 10), incubate 2 hours, wash.
- Step 3.** Add 25 μ L of Detection Antibody, incubate 2 hours, wash.
- Step 4.** Add 150 μ L 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.



Notes:

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.
 - *Prepare 1:3 serial dilutions beginning with the High Calibrator, by adding 100 μL of High Calibrator to 200 μL of Diluent 10. Prepare 6 serial dilutions. The first Calibrator will be High Calibrator stock and the 8th 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 KDR.*
 - *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.*
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 μL of diluted sample.

Prepare Detection Antibody Reagent:

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

Detection Antibody Reagent is stable at RT for a few hours.

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.



Notes:

Assay Protocol

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

1. Add 150 μ L/well of Blocker A Solution and incubate at room temperature for 1-2 hours or overnight at 4 $^{\circ}$ C.
2. Wash plates 3 times with Phosphate Buffered Saline + 0.05% Tween-20 (PBS-T).
3. Add 50 μ 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 μ 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 μ 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.

