

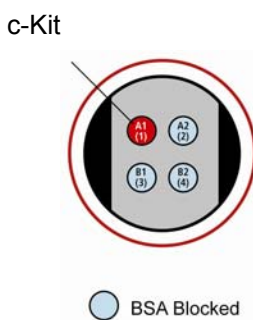
# MSD<sup>®</sup> 96-Well MULTI-ARRAY<sup>®</sup> Total c-Kit Assay

The following assay protocol has been optimized for quantifying Total c-Kit in whole cell lysate.

Storage

## MSD Materials

<input type="checkbox"/> Read Buffer T (with surfactant), 4X	RT
<input type="checkbox"/> Blocker A	RT
<input type="checkbox"/> MULTI-SPOT <sup>®</sup> 96-well 4 Spot Total c-Kit plates(s)	2-8°C
<input type="checkbox"/> SULFO-TAG <sup>™</sup> Anti-Total c-Kit Antibody (50X)	2-8°C
<input type="checkbox"/> Tris Wash Buffer (10X)	2-8°C
<input type="checkbox"/> Tris Lysis Buffer (1X)	2-8°C
<input type="checkbox"/> Phosphatase Inhibitor I (100X)	2-8°C
<input type="checkbox"/> Phosphatase Inhibitor II (100X)	2-8°C
<input type="checkbox"/> Protease Inhibitor Solution (50X)	≤-10°C



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

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



## Notes:

### *Other Materials & Equipment (not supplied)*

- Deionized water for diluting Wash Buffer and Read Buffer
- One 250 mL bottle
- Two 50 mL tube
- One 15 mL tube
- Adhesive plate seals
- Microtiter plate shaker
- Various microcentrifuge tubes for making serial dilutions of lysates (if desired)
- Automated 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  $\mu$ L and 150  $\mu$ L into a 96-well micro plate

*Read the entire detailed instructions before beginning work.*

### *Protocol at a Glance*

The following protocol describes the most conservative approach toward achieving highly sensitive results using MSD technology to quantify phosphoproteins. The protocol takes approximately 3 to 3½ hours to complete if each reagent is prepared during the preceding incubation. All reagents with the exception of diluted lysates can also be prepared ahead of time. This lengthens the overall time required for the assay but frees up time during incubation steps.

Once desired results are achieved, the protocol can be streamlined to eliminate multiple incubation and wash steps to increase throughput.

1. Add blocking solution, incubate 1 hour, wash.
2. Add sample or lysate, incubate 1 hour, wash.
3. Add detection antibody, incubate 1 hour, wash.
4. Add Read Buffer and analyze plate.

*A larger amount of Wash Buffer may be prepared and stored at room temperature for later use.*

### *Detailed Instructions*

#### ***Prepare a stock of 1X Tris Wash Buffer:***

- a) 1X Tris Wash Buffer will be used throughout the assay to make other reagents as well as wash plates. Approximately 250 mL per plate is required– more if using an automatic plate washer.
- b) In a 250 mL bottle combine:
  - 25 mL 10X Tris Wash Buffer
  - 225 mL deionized water



**Notes:**

Solutions containing Blocker A should be kept at 4°C and discarded after 14 days.

**Prepare Blocking Solution-A:**

- a) Prepare 20 mL per plate.
- b) In a 50 mL tube combine:
  - 20 mL 1X Tris Wash Buffer
  - 600 mg Blocker A (30 mg/mL or 3%)

**Prepare Antibody Dilution Buffer:**

- a) Prepare 3 mL per plate.
- b) In a 15 mL tube combine:
  - 1 mL Blocking Solution-A
  - 2 mL 1X Tris Wash Solution

Begin with a MULTI-SPOT Total c-Kit Plate.  
No pre-treatment is necessary.

**STEP 1**

Add 150 µL/well of Blocking Solution-A.

Incubate with shaking at room temperature for 1 hour. Prepare Complete Tris Lysis buffer, and prepare samples or dilute cell lysates during this time.

Plates may also be blocked overnight at 4°C.

Complete Tris Lysis Buffer should be made each day of experimentation.

**Prepare Complete Tris Lysis Buffer:**

- a) To 10 mL of Tris Lysis Buffer, add the following:
  - 200 µL Protease Inhibitor Solution (50X stock)
  - 100 µL Phosphatase Inhibitor I (100X stock)
  - 100 µL Phosphatase Inhibitor II (100X stock)
- b) Keep Complete Tris Lysis Buffer on ice until use.

The Complete Tris Lysis Buffer should be ice cold before use.

**Prepare samples or positive and negative cell lysates:**

*(Note: Recommendations for cell lysate handling are provided, however, the suggested concentrations listed below may need to be adjusted depending upon specific samples tested.)*

- a) Thaw cell lysate samples on ice and dilute immediately before use. Keep on ice during all manipulations and discard all remaining thawed unused material.
- b) Dilute cell lysate in Complete Tris Lysis Buffer to a final concentration of 0.4 µg/µL. This will deliver 10 µg/well in 25 µL. A dilution series may also be prepared if desired.

Wash plates four times with Wash Buffer.



**Notes:**

**STEP 2** Dispense 25  $\mu$ L/well of samples or diluted lysates.

Incubate with shaking at room temperature for 1 hour. Prepare Detection Antibody Solution during this time.

*Shaking a 96-well MSD MULTI-ARRAY or MULTI-SPOT plate accelerates capture at the working electrode.*

**Prepare Detection Antibody Solution:**

- a) Prepare 3.0 mL per plate.
- b) In a 15 mL tube combine:
  - a. 2.94 mL cold Antibody Dilution Buffer
  - b. 60  $\mu$ L 50X Anti-Total c-Kit Antibody  
(Final concentration: 1X)

Wash plates four times with Wash Buffer.

**STEP 3** Add 25  $\mu$ L/well of Detection Antibody.

Incubate with shaking at room temperature for 1 hour. Prepare Read Buffer during this time.

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

**Dilute Read Buffer:**

- a) Approximately 20 mL per plate is required.
- b) Dilute 4X Read Buffer T (with surfactant) to 1X with deionized water.

Wash plates four times with Wash Buffer.

*Bubbles in the Read Buffer will interfere with reliable imaging of the plate if carried into the wells.*

**STEP 4** Add 150  $\mu$ L of diluted Read Buffer T (with surfactant).

Analyze with SECTOR Imager instrument.

*Plates can be imaged immediately following the addition of read buffer. Most biological interactions tolerate incubation in Read Buffer however each unique assay should be tested for stability in read buffer before being left to sit for extended periods.*

