

# MESO SCALE DISCOVERY® Labeled Reporters

## MSD® SULFO-TAG Labeled Anti-Phospho-Tyrosine Antibody

Size	Part Number	Volume
50 µg	R32AP-5	100 µL
200 µg	R32AP-1	400 µL

<b>Description</b>	MSD SULFO-TAG™ labeled Anti-Phospho Tyrosine Antibody may be used as a primary detection reagent to generate electrochemiluminescence in MESOSCALE DISCOVERY assays. It allows the user to detect analytes containing phosphorylated tyrosine residues.
<b>Immunogen</b>	Phosphotyrosine conjugated to carrier protein
<b>Isotype</b>	IgG2b
<b>Molecular Weight</b>	The average molecular weight is of SULFO-TAG labeled Anti-Phospho Tyrosine Antibody is approximately 160,000 Da.
<b>Specificity</b>	MSD SULFO-TAG labeled Anti-Phospho Tyrosine Antibody detects phosphorylated tyrosine in native and denatured proteins. It is specific for phospho-tyrosine and does not react with either phospho-serine or phospho-threonine.
<b>Protein Concentration</b>	The Antibody is supplied as 500 µg/mL solution in Phosphate Buffered Saline (PBS) containing 2% Bovine Serum Albumin, 0.05% Azide and 0.02% Kathon CG.
<b>Storage</b>	MSD SULFO-TAG labeled Anti-Phospho Tyrosine Antibody should be stored at 2-8°C, shielded from light. It should be diluted to the working concentration only prior to immediate use.
<b>Stability</b>	Under recommended storage conditions, MSD SULFO-TAG labeled Anti-Phospho Tyrosine Antibody is stable until the expiration date on the label. Short term exposure of SULFO-TAG antibody to ambient light when carrying out assays is generally not a concern.
<b>Recommended Dilution(s)</b>	Suggested dilution concentrations range from 0.2 – 2 µg/mL for initial assay development. As a starting point, MSD recommends diluting MSD SULFO-TAG labeled Anti-Phospho Tyrosine Antibody to 1 µg/mL and incubating for 1 hour (25µL/well). If background signals are higher than desired, reduce the concentration of MSD SULFO-TAG labeled Anti-Phospho Tyrosine Antibody to 0.5 µg/mL. Optimal concentrations may be determined by the researcher for each application.

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

### Ordering Information

#### MSD Customer Service

Phone: 1-301-947-2085

Fax: 1-301-990-2776

Email: [CustomerService@mesoscale.com](mailto:CustomerService@mesoscale.com)

#### MSD Scientific Support

Phone: 1-301-947-2025

Email: [ScientificSupport@mesoscale.com](mailto:ScientificSupport@mesoscale.com)

### Company Address

#### Meso Scale Discovery

A division of Meso Scale Diagnostics, LLC.

9238 Gaither Road

Gaithersburg, MD 20877 USA

[www.mesoscale.com](http://www.mesoscale.com)

