

MESO SCALE DISCOVERY® Labeled Reporters

MSD® SULFO-TAG Labeled Anti-Phospho Tyrosine Antibody

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

Description	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.
Clone	PY20
Immunogen	Phosphotyrosine conjugated to carrier protein
Isotype	IgG2b
Molecular Weight	The average molecular weight is of SULFO-TAG labeled Anti-Phospho Tyrosine Antibody is approximately 160,000 Da.
Specificity	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.
Protein Concentration	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.
Storage	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.
Stability	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.
Recommended Dilution(s)	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

MSD Scientific Support

Phone: 1-301-947-2025

Email: 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

