

MESO SCALE DISCOVERY® Labeled Reporters

MSD® SULFO-TAG Labeled Anti-Ubiquitinated Protein Antibody

Size	Part Number	Quantity
50 µg	R32AU-5	100 µL
200 µg	R32AU-1	400 µL

Description	MSD SULFO-TAG™ labeled Anti-Ubiquitinated Protein Antibody may be used as a primary detection reagent to generate electrochemiluminescence in MESOSCALE DISCOVERY assays. It allows the user to detect ubiquitinated proteins.
Clone	FK2
Isotype	IgG1
Specificity	MSD SULFO-TAG labeled Anti-Ubiquitinated Protein Antibody is specific for ubiquitin-protein conjugates and shows no reactivity with free ubiquitin. The antibody recognizes both mono- and poly-ubiquitinated species.
Host species	Mouse
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-Ubiquitinated Protein 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-Ubiquitinated Protein 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-Ubiquitinated Protein 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-Ubiquitinated Protein 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

