

# MESO SCALE DISCOVERY® Labeled Reporters

## MSD® SULFO-TAG Labeled Anti-Ubiquitinylated Protein Antibody

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

<b>Description</b>	MSD SULFO-TAG™ labeled Anti-Ubiquitinylated Protein Antibody may be used as a primary detection reagent to generate electrochemiluminescence in MESOSCALE DISCOVERY assays. It allows the user to detect ubiquitinylated proteins.
<b>Isotype</b>	IgG1
<b>Specificity</b>	MSD SULFO-TAG labeled Anti-Ubiquitinylated Protein Antibody is specific for ubiquitin-protein conjugates and shows no reactivity with free ubiquitin. The antibody recognizes both mono- and poly-ubiquitinylated species.
<b>Host species</b>	Mouse
<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-Ubiquitinylated 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.
<b>Stability</b>	Under recommended storage conditions, MSD SULFO-TAG labeled Anti-Ubiquitinylated 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.
<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-Ubiquitinylated 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-Ubiquitinylated 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](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)

