

MESO SCALE DISCOVERY® Labeled Reporters

MSD® SULFO-TAG Labeled Anti-Rabbit Antibody (Goat)

Size	Part Number	Quantity
50 µg	R32AB-5	100 µL
200 µg	R32AB-1	400 µL

Description	MSD SULFO-TAG™ labeled Anti-Rabbit Antibody may be used as a secondary detection reagent to generate electrochemiluminescence in MESOSCALE DISCOVERY assays. It allows the use of commercially available rabbit antibodies as detection in MSD assays without directly labeling them.
Immunogen	Rabbit IgG whole molecule
Specificity	MSD SULFO-TAG labeled Anti-Rabbit Antibody is a goat polyclonal antibody and binds to heavy and light chains of rabbit IgG. No reaction has been observed against Bovine, Chicken, Goat, Guinea Pig, Hamster, Horse, Human, Mouse, Rat and Sheep antibodies.
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-Rabbit Antibody should be stored at 2-8°C, shielded from light. It should be diluted to the working concentration only prior to immediate use.
Purity	This product was prepared from monospecific antiserum by immunoaffinity chromatography using Rabbit IgG coupled to agarose beads followed by solid phase adsorption(s) to remove any unwanted reactivities.
Stability	Under recommended storage conditions, MSD SULFO-TAG labeled Anti-Rabbit 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-Rabbit Antibody to 1 µg/mL and incubating for 1 hour (25µL/well) following the primary antibody incubation. If background signals are higher than desired, reduce the concentration of MSD SULFO-TAG labeled Anti-Rabbit Antibody to 0.5 µg/mL. Alternatively, a single incubation step using both labeled and unlabeled antibodies at 1:1 molar ratio can be tested. 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

