

MESO SCALE DISCOVERY® Labeled Reporters

MSD® SULFO-TAG Labeled Anti-Glutathione-S-Transferase Antibody (Goat)

Size	Part Number	Volume
50 µg	R32AA-5	100 µL
200 µg	R32AA-1	400 µL

Description	MSD SULFO-TAG™ labeled Anti-Glutathione-S-Transferase Antibody may be used as a primary detection reagent to generate electrochemiluminescence in MESOSCALE DISCOVERY assays. It allows the use of commercially available GST-tagged proteins as detection in MSD assays without directly labeling them.
Immunogen	Glutathione-S-Transferase
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-GST 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 GST 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-GST 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-GST 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-GST 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

