

MESO SCALE DISCOVERY® Labeled Reporters

MSD® SULFO-TAG Labeled Streptavidin

Size	Part Number	Quantity
50 µg	R32AD-5	100 µL
200 µg	R32AD-1	400 µL

Description	MSD SULFO-TAG™ labeled Streptavidin may be used as a secondary detection reagent to generate electrochemiluminescence in MESOSCALE DISCOVERY assays. SULFO-TAG Streptavidin is used to report biotin-labeled molecules. It allows the use of commercially available biotinylated detection antibodies in MSD assays. The average molecular weight of SULFO-TAG Streptavidin is approximately 60 kDa.
Specificity	Streptavidin binds to biotin and biotin analogs with high specificity and high affinity. The dissociation constant for streptavidin and biotin is approximately 10^{-15} M.
Protein Concentration	MSD SULFO-TAG labeled Streptavidin 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 Streptavidin 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 Streptavidin is stable until the expiration date on the label. Short term exposure of SULFO-TAG Streptavidin 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 Streptavidin 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 Streptavidin 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

