

Meso Scale Discovery Labeled Reporter Reagents

MSD [®] Label	Reporter	Host Species	Immunogen	Quantity	Catalog Number
SULFO-TAG [™]	Anti-Rabbit Antibody	Goat	Rabbit IgG Whole Molecule	50 µg	R32AB-5
SULFO-TAG	Anti-Rabbit Antibody	Goat	Rabbit IgG Whole Molecule	200 µg	R32AB-1
SULFO-TAG	Anti-Mouse Antibody	Goat	Mouse IgG Whole Molecule	50 µg	R32AC-5
SULFO-TAG	Anti-Mouse Antibody	Goat	Mouse IgG Whole Molecule	200 µg	R32AC-1
SULFO-TAG	Anti-Goat Antibody	Donkey	Goat IgG Whole Molecule	50 µg	R32AG-5
SULFO-TAG	Anti-Goat Antibody	Donkey	Goat IgG Whole Molecule	200 µg	R32AG-1
SULFO-TAG	Anti-Rat Antibody	Goat	Rat IgG Whole Molecule	50 µg	R32AH-5
SULFO-TAG	Anti-Rat Antibody	Goat	Rat IgG Whole Molecule	200 µg	R32AH-1
SULFO-TAG	Anti-Human Antibody	Goat	Human IgG Whole Molecule	50 µg	R32AJ-5
SULFO-TAG	Anti-Human Antibody	Goat	Human IgG Whole Molecule	200 µg	R32AJ-1
SULFO-TAG	Anti-Sheep Antibody	Donkey	Sheep IgG Whole Molecule	50 µg	R32AI-5
SULFO-TAG	Anti-Sheep Antibody	Donkey	Sheep IgG Whole Molecule	200 µg	R32AI-1
SULFO-TAG	Streptavidin	N/A	N/A	50 µg	R32AD-5
SULFO-TAG	Streptavidin	N/A	N/A	200 µg	R32AD-1
SULFO-TAG	Anti-GST Antibody	Goat	Glutathione-S-transferase	50 µg	R32AA-5
SULFO-TAG	Anti-GST Antibody	Goat	Glutathione-S-transferase	200 µg	R32AA-1
SULFO-TAG	Anti-Phosphotyrosine Antibody	Mouse	KLH-pTyr, Gly & Ala (equal parts) (PY20)	50 µg	R32AP-5
SULFO-TAG	Anti-Phosphotyrosine Antibody	Mouse	KLH-pTyr, Gly & Ala (equal parts) (PY20)	200 µg	R32AP-1
SULFO-TAG	Anti-Ubiquitylated Protein Antibody	Mouse	Poly-ubiquitylated lysozyme	50 µg	R32AU-5
SULFO-TAG	Anti-Ubiquitylated Protein Antibody	Mouse	Poly-ubiquitylated lysozyme	200 µg	R32AU-1
SULFO-TAG	Anti Bio-Block A Antibody	Mouse	QuickPlex [™] Bio-Block A	50 µg	R32DA-5
SULFO-TAG	Anti Bio-Block A Antibody	Mouse	QuickPlex Bio-Block A	200 µg	R32DA-1
SULFO-TAG	Anti Bio-Block B Antibody	Mouse	QuickPlex Bio-Block B	50 µg	R32DB-5
SULFO-TAG	Anti Bio-Block B Antibody	Mouse	QuickPlex Bio-Block B	200 µg	R32DB-1
SULFO-TAG	Anti Bio-Block C Antibody	Rat	QuickPlex Bio-Block C	50 µg	R32DC-5
SULFO-TAG	Anti Bio-Block C Antibody	Rat	QuickPlex Bio-Block C	200 µg	R32DC-1
SULFO-TAG	Anti Bio-Block D Antibody	Rabbit	QuickPlex Bio-Block D	50 µg	R32DD-5
SULFO-TAG	Anti Bio-Block D Antibody	Rabbit	QuickPlex Bio-Block D	200 µg	R32DD-1
TAG	Anti-Rabbit Antibody	Goat	Rabbit IgG Whole Molecule	200 µg	R31AB-1
TAG	Anti-Mouse Antibody	Goat	Mouse IgG Whole Molecule	200 µg	R31AC-1
TAG	Streptavidin	N/A	N/A	500 µg	R31AD-2
TAG	Anti-GST Antibody	Goat	Glutathione-S-transferase	200 µg	R31AA-1

General Guidelines for Secondary Antibody Conjugates:

- Suggested dilution concentrations range from 0.2 - 2 µg/mL for initial assay development.
- As a starting point, MSD recommends diluting the labeled secondary antibody to 1 µg/mL, and incubating for 1 hour (25 µL/well) following the unlabeled antibody incubation.
- If background signals are higher than desired, reduce the labeled secondary antibody concentration to 0.5 µg/mL.
- Ultimately, the protocol should evolve toward a single incubation step using both labeled and unlabeled antibodies at a 1:1 molar ratio.

General Guidelines for SULFO-TAG[™] Streptavidin Reagent:

- Suggested dilution concentrations range from 0.2 - 2 µg/mL for initial assay development.
- SULFO-TAG Streptavidin can be used to convert ELISA-based assays to MSD[®] technology by directly replacing streptavidin detection systems (e.g. SA-HRP).
- In these cases, the capture antibody used in the ELISA can be directly immobilized on the surface of an MSD[®] plate.
- As a starting point, MSD recommends diluting the SULFO-TAG Streptavidin to 1 µg/mL, and incubating for 1 hour (25 µL/well).

For research use only; not for use in diagnostic or therapeutic procedures.



©2005 Meso Scale Discovery, LLC. All Rights Reserved.