

MSD[®] SULFO-TAG–Conjugated Anti-Hu/NHP Lambda Light Chain Antibody



Ordering Information

MSD Customer Service
Phone: 1-240-314-2795
Fax: 1-301-990-2776
Email: CustomerService@mesoscale.com
www.mesoscale.com/support

Scientific Support

Phone: 1-240-314-2798
Email: ScientificSupport@mesoscale.com
www.mesoscale.com/support

Company Address

MESO SCALE DISCOVERY[®]
A division of
Meso Scale Diagnostics, LLC.
1601 Research Boulevard
Rockville, MD 20850-3173 USA

www.mesoscale.com[®]

**For Research Use Only.
Not for use in
diagnostic procedures.**

Catalog Number	D20QG-6
Contents	SULFO-TAG [™] –Conjugated Anti-Hu/NHP Lambda Light Chain Antibody
Size	100 µg
Volume	200 µL
Storage	2–8°C, shielded from light

Summary and Intended Use

MSD SULFO-TAG–Conjugated Anti-Hu/NHP Lambda is a goat polyclonal antibody that reacts with all immunoglobulin isotypes and may be used as a primary or secondary detection reagent in MSD assays. This antibody has been tested with human and cynomolgus monkey samples, but may also cross-react with other non-human primate species. In addition to using this antibody for isotyping immunoglobulins in biological samples, this antibody can also be used for immunogenicity applications.

Immunogen

Human myeloma proteins with lambda light chains.

Specificity

MSD SULFO-TAG–Conjugated Anti-Hu/NHP Lambda antibody can recognize all immunoglobulin isotypes (IgA, IgG, IgM, and IgE). It is a goat polyclonal antibody.

Protein Concentration

The antibody is supplied as a 500 µg/mL solution in phosphate buffered saline (PBS) containing 0.05% azide and preservatives.

Purity

This product was purified with Zeba Spin Desalting Columns to remove any unconjugated SULFO-TAG.

Stability

Under recommended storage conditions, MSD SULFO-TAG–Conjugated Anti-Hu/NHP Lambda Antibody is stable until the expiration date on the label. Short term exposure of SULFO-TAG antibody to ambient light during the assay is generally not a concern.

Safety

Use safe laboratory practices and wear gloves, safety glasses, and lab coats when handling antibodies. Handle and dispose of all hazardous samples properly in accordance with local, state, and federal guidelines. Refer to the safety data sheet (SDS), included in the product package and available from MSD Customer Service, for additional product safety and handling information.

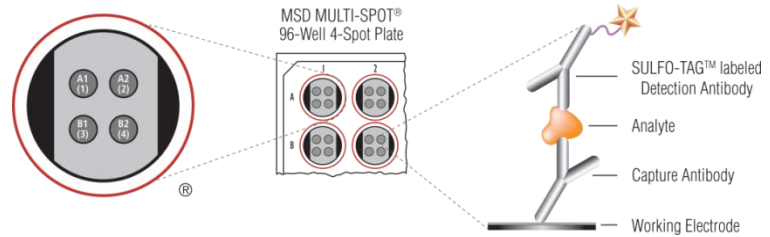
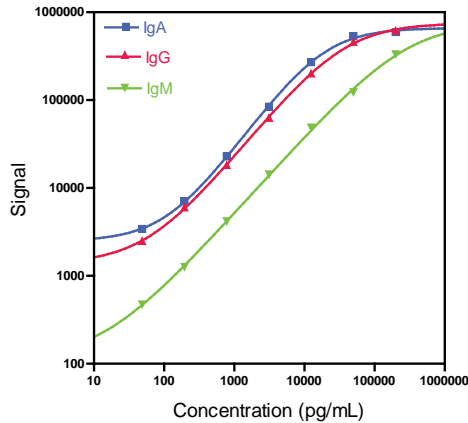
Recommended Dilution(s)

For initial assay development, the suggested concentrations of SULFO-TAG–Conjugated Anti-Hu/NHP Lambda antibody range from 0.2 to 2 µg/mL. As a starting point, MSD recommends diluting the antibody to 1 µg/mL and incubating for 1 hour (25 µL/well) following sample incubation. If background signals are higher than desired, reduce the concentration of MSD SULFO-TAG–Conjugated Anti-Hu/NHP Kappa antibody to 0.5 µg/mL or lower. Optimal concentrations for each application should be determined by the researcher.

MSD Labeled Reporters

Representative Data

The representative calibration curves for IgA, IgG, and IgM shown below were determined using MSD Isotyping Panel 1 (Human/NHP) MULTI-SPOT® Plate that is pre-coated with capture antibodies for IgA, IgG, and IgM on independent and well-defined spots in the layout shown below. A blend of SULFO-TAG–Conjugated Anti-Hu/NHP Kappa and Anti-Hu/NHP Lambda Light Chain antibodies was used as the detector in the assays. The reported data should be used only as a guide and may vary with between lots.



MULTI-SPOT 96-well Isotyping Panel 1 (Human/NHP) Plate

Spot Layout: A1: IgG, A2: IgA, B1: IgM, B2: BSA blocked

IgA		
Conc. (pg/mL)	Average Signal	%CV
0	2 054	6.2
49	3 431	1.0
195	7 076	5.1
781	22 890	1.0
3 125	84 314	0.6
12 500	270 399	3.8
50 000	536 227	9.8
200 000	593 623	14.5

IgG		
Conc. (pg/mL)	Average Signal	%CV
0	1 393	7.5
49	2 495	3.0
195	6 021	4.7
781	18 289	8.2
3 125	63 044	2.4
12 500	200 605	3.8
50 000	460 116	11.0
200 000	620 218	7.1

IgM		
Conc. (pg/mL)	Average Signal	%CV
0	188	5.3
49	473	0.6
195	1 261	2.1
781	4 186	4.7
3 125	14 261	6.8
12 500	48 785	0.3
50 000	122 944	8.9
200 000	332 323	4.3

Related Products

MSD offers a range of immunoassay kits and reagents for isotyping applications. The catalog numbers for some of the products are provided below. Visit www.mesoscale.com for a complete list of available products.

	Product Name	Catalog Number
Assay Kits	Isotyping Panel 1 (Human/NHP) Kit	K15203D
	Human/NHP IgA Kit	K150JJD
	Human/NHP IgG Kit	K150JLD
	Human/NHP IgM Kit	K150JPD
Labeled Reporters	SULFO-TAG–Conjugated Anti-Hu/NHP IgA Antibody	D20JJ
	SULFO-TAG–Conjugated Anti-Hu/NHP IgG Antibody	D20JL
	SULFO-TAG–Conjugated Anti-Hu/NHP IgM Antibody	D20JP
	SULFO-TAG–Conjugated Anti-Hu/NHP Kappa Light Chain Antibody	D20TF
	SULFO-TAG–conjugated Streptavidin	R32AD

MESO SCALE DISCOVERY, MSD, MULTI-SPOT, SULFO-TAG, www.mesoscale.com, 96 WELL 4-SPOT (design), MSD (design), and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC. ©2013 Meso Scale Diagnostics, LLC. All rights reserved.

