

# MSD<sup>®</sup> SULFO-TAG–Conjugated Anti-Hu/NHP IgA Antibody



## Ordering Information

MSD Customer Service  
Phone: 1-240-314-2795  
Fax: 1-301-990-2776  
Email: [CustomerService@mesoscale.com](mailto:CustomerService@mesoscale.com)  
[www.mesoscale.com/support](http://www.mesoscale.com/support)

## Scientific Support

Phone: 1-240-314-2798  
Email: [ScientificSupport@mesoscale.com](mailto:ScientificSupport@mesoscale.com)  
[www.mesoscale.com/support](http://www.mesoscale.com/support)

## Company Address

MESO SCALE DISCOVERY<sup>®</sup>  
A division of  
Meso Scale Diagnostics, LLC.  
1601 Research Boulevard  
Rockville, MD 20850-3173 USA

[www.mesoscale.com](http://www.mesoscale.com)<sup>®</sup>

<b>Catalog Number</b>	D20JJ-6
<b>Contents</b>	SULFO-TAG <sup>™</sup> –Conjugated Anti-Hu/NHP IgA Antibody
<b>Size</b>	100 µg
<b>Volume</b>	200 µL
<b>Storage</b>	2–8°C, shielded from light

## Summary and Intended Use

MSD SULFO-TAG–Conjugated Anti-Hu/NHP IgA Antibody is a monoclonal antibody that reacts with immunoglobulin A (IgA) and may be used as a primary or secondary detection reagent in MSD assays. This antibody has been tested with human, cynomolgus monkey, and rhesus monkey samples, but may also cross-react with other non-human primate species. In addition to using this antibody for detection of IgA in biological samples, it can also be used for immunogenicity applications.

## Immunogen

Purified IgA from human colostrum.

## Specificity

MSD SULFO-TAG–Conjugated Anti-Hu/NHP IgA Antibody does not cross-react with other immunoglobulin isotypes (IgE, IgG, or IgM). This mouse monoclonal antibody has been shown to detect IgA in human and NHP sera at 250 000-fold dilution.

## 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 IgA 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)

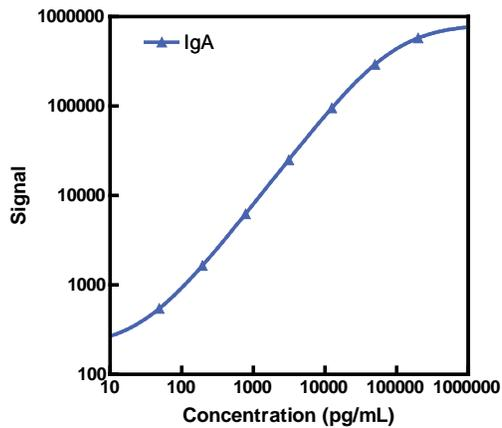
For initial assay development, the suggested concentrations of SULFO-TAG–Conjugated Anti-Hu/NHP IgA Antibody range from 0.2 to 2 µg/mL. As a starting point, MSD recommends diluting the antibody to 1 µg/mL and incubating for 1 hour (25 µL/well) following sample incubation. If background signals are higher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

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

# MSD Labeled Reporters

## Representative Data

The representative calibration curve for IgA shown below was determined using SULFO-TAG–Conjugated Anti-Hu/NHP IgA Antibody with MSD Isotyping Panel 1 (Human/NHP) Kit (catalog#K15203D). The reported data should be used only as a guide and may vary between lots.



IgA		
Conc. (pg/mL)	Average Signal	%CV
0	190	5.2
49	545	14.3
195	1 646	42.3
781	6 244	3.3
3 125	25 019	10.2
12 500	95 347	7.5
50 000	292 294	3.2
200 000	575 772	0.5

## 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](http://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 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 Anti-Hu/NHP Lambda Light Chain Antibody	D20QG
	SULFO-TAG–Conjugated Streptavidin	R32AD