

MSD[®] SULFO-TAG–Conjugated Anti-Hu/NHP Kappa 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

| | |
|-----------------------|--|
| Catalog Number | D20TF-6 |
| Contents | SULFO-TAG™–Conjugated Anti-Hu/NHP Kappa Light Chain Antibody |
| Size | 100 µg |
| Volume | 200 µL |
| Storage | 2–8°C, shielded from light |

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[®]

Summary and Intended Use

MSD SULFO-TAG–Conjugated Anti-Hu/NHP Kappa Light Chain Antibody 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 kappa light chains.

Specificity

MSD SULFO-TAG–Conjugated Anti-Hu/NHP Kappa Light Chain 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 Kappa Light Chain 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)

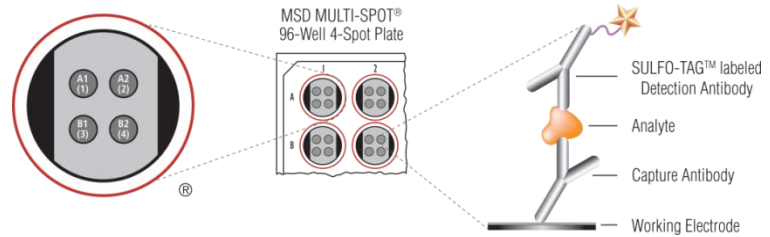
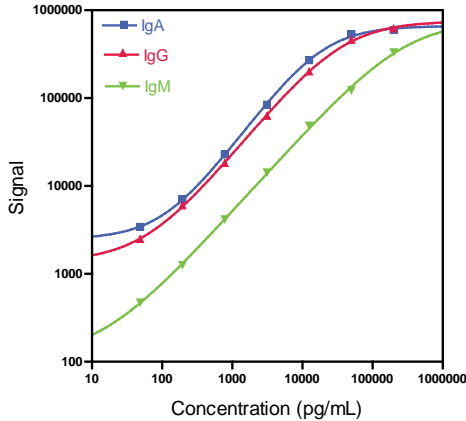
For initial assay development, the suggested concentrations of SULFO-TAG–Conjugated Anti-Hu/NHP Kappa Light Chain Antibody range from 0.2 to 2 µg/mL. As a starting point, MSD recommends diluting the antibody to 1 µg/mL and incubating for 1 hour (25 µL/well) following sample incubation. If background signals are higher than desired, reduce the concentration of MSD SULFO-TAG–Conjugated Anti-Hu/NHP Kappa Light Chain 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 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 Lambda Light Chain Antibody | D20QG |
| | 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.

