

# MSD® SULFO-TAG Conjugated DIG 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)

<b>Catalog No.</b>	D20ARF-3
<b>Size</b>	5 plates
<b>Volume</b>	165 µL
<b>Storage</b>	2–8°C, shielded from light

## Technical Support

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

## Summary and Intended Use

MESO SCALE DISCOVERY® SULFO-TAG Conjugated DIG Antibody is a monoclonal antibody that reacts with digoxigenin and may be used as a detection reagent in MSD assays. This antibody can be used for analytes labeled with digoxigenin, including oligonucleotide probes, recombinant proteins, antibodies, or other biomolecules.

## Protein Concentration

The antibody is supplied as a 25 µg/mL (200X) solution in Diluent 100 containing 0.05% azide and preservatives.

## Purity

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

## Stability

Under recommended storage conditions, MSD SULFO-TAG Conjugated DIG 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 DIG Antibody range from 0.1 to 1.0 µg/mL. As a starting point, MSD recommends diluting the antibody to 0.125 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 DIG Antibody. Optimal concentrations for each application should be determined by the researcher.

## Recent Publications

Performance of a Digoxigenin Antibody-Based Electrochemiluminescent Hybridization Assay: A HELISA Alternative for Low Femtomolar Detection of an Antisense Oligonucleotide [Poster](#)

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

Please use the following link for a list of the trademarks and service marks owned by Meso Scale Diagnostics, LLC. and Methodical Mind, LLC. <https://www.mesoscale.com/trademarks>. All other trademarks and service marks are the property of their respective owners.

©2013-2026 Meso Scale Diagnostics, LLC. All rights reserved.

