



MSD[®] Conjugation Services

High quality reagents accelerate the development of assays and help manage the variability that can occur from lot to lot. MESO SCALE DISCOVERY[®] offers custom Conjugation Services for linking antibodies, proteins, peptides, and oligonucleotides with either **biotin** or **SULFO-TAG[™]**, a label used with MSD's electrochemiluminescence platform. These services leverage MSD's expertise and knowledge in reagent manufacturing, providing researchers with a convenient and quick route to well-characterized and analytically robust assay development tools.

Service	MSD GOLD [™] Conjugation Services	MSD Conjugation Service
Antibody buffer exchanged to remove interfering materials	✓	✓
Add-on size exclusion chromatography to remove aggregates (optional)	✓	
Pre-conjugation analytical reagent characterization	✓	
Conjugation according to optimized procedures under MSD's Quality Management System	✓	✓
Post-conjugation purification to remove free label	✓	✓
Post-conjugation measurement of labels per protein	✓	✓
Post-conjugation characterization	✓	
Vialing service (optional)	✓	✓
Certificate of Analysis	Includes characterization data	✓

MSD GOLD Conjugation Service

- Critical for customers who need reproducibility and well-characterized assays
- Includes additional characterization and quality assessment steps
- Identifies common antibody issues such as degradation, aggregation, purity, and changes in source material
- Ensures superior antibody conjugation, compared to standard Conjugation Service, for better assay performance and longitudinal studies



Conjugation Optimization Packages

MSD offers the following service packages in conjunction with its Gold conjugation service to specifically support assay development programs.

Challenge Ratio Optimization:

- Provides ability to correlate functional performance to incorporated labels
- Antibody conjugated at three different challenge ratios

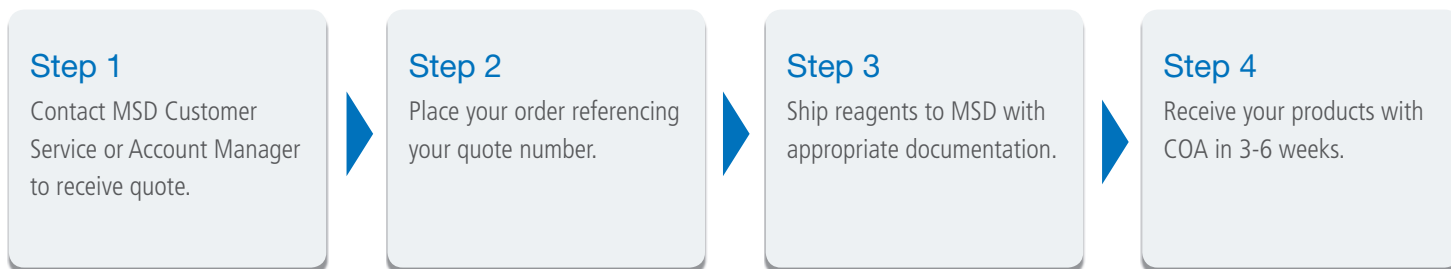
Multiple Conjugations:

- Measures antibody conjugation reproducibility with three separate conjugations of the same lot of material
- Conjugated on separate days for assessment of true lot-to-lot reproducibility

Multiple Reagent Lots:

- Identifies potential risks due to raw material variability
- Three separate lots of antibody raw materials are analytically characterized and conjugated

Contact us to discuss your needs.



Phone: 1-240-314-2795 . Fax: 1-301-990-2776
Email: customerservice@mesoscale.com

MESO SCALE DISCOVERY, MSD, SULFO-TAG, www.mesoscale.com, MSD (design), MSD GOLD, and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.
© 2015 Meso Scale Diagnostics, LLC. All rights reserved.



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

MK-DS-220-v1-2015Jan

Reproducibility of Conjugation

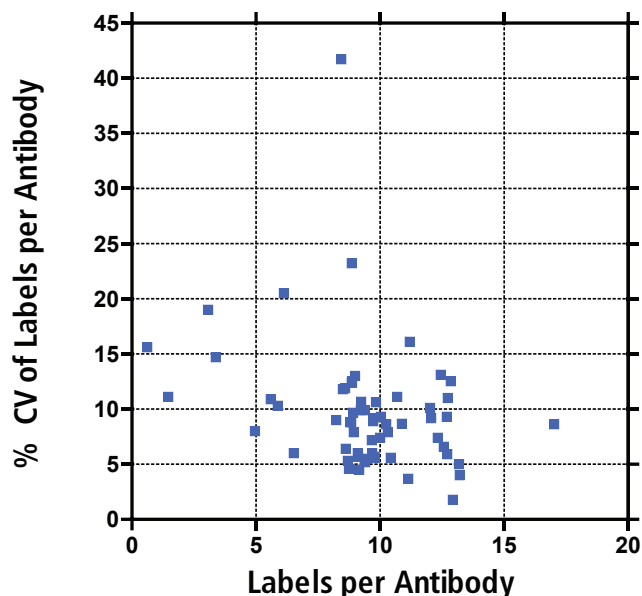


Figure 1. %CV of the labels per antibody for 61 antibodies conjugated more than five times each over a three-year period.