



CERTIFICATE OF ANALYSIS

COA - 04758

Product Description: SULFO-TAG™ Goat Anti-Mouse Antibody

Catalog #/Size: R32AC-5/50 µg; R32AC-1/200 µg

Lot Number: W0014077S

Expiration Date: 30 JUN 2014

Statement: This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

Concentration: 500 µg/mL



Volume: 100 µL; 400 µL

L/P: 5.33

Diluent: Diluent 100 w/ Azide

Storage conditions: 2-8°C

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Preparer	Meredith Hutchinson	Inventory Technician		16 SEP 2014
Review/Approval	Staci Leffel	Manufacturing Manager		19 SEP 2014