



CERTIFICATE OF ANALYSIS

COA - 11355

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

Catalog #/Size: R32AJ-5/50 µg; R32AJ-1/200 µg

Lot Number: W0014709S

Expiration Date: 31 JAN 2016

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

Volume: 100 µL; 400 µL

Concentration: 500 µg/mL

L/P: 9.1

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 Yeager	Inventory Technician		27 Nov 2013
Review/Approval	Staci Leffel	Manufacturing Manager		27 Nov 2013