



CERTIFICATE OF ANALYSIS

COA - 05057

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

Catalog #/Size: R32AG-5/50 µg; R32AG-1/200 µg

Lot Number: W0015194S

Expiration Date: 30 NOV 2016

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

Concentration: 500 µg/mL

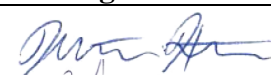

Volume: 100 µL; 400 µL

L/P: 12.6

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		09 JAN 2015
Review/Approval	Latrice Corsey	Manufacturing Lead		15 JAN 2015