



CERTIFICATE OF ANALYSIS

COA - 07529

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

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

Lot Number: W0016553S

Expiration Date: 31 JAN 2019

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

Volume: 400 µL; 100 µL



Concentration: 500 µg/mL

L/P: 10.0

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 | Production Control |  | 13 JAN 2017 |
| Review/Approval | Dawn Stollar | Quality Assurance |  | 17 JAN 2017 |