



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog #/Size: R32AG-Series

Lot Number: W0017540S

Expiration Date: 31 MAR 2020

Statement: This product was manufactured and tested according to approved MSD documents.

Total Volume: 400 µL

Concentration: 500 µg/mL

L/P: 10.7

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 |
|------------------------|--------------|-------------------|---------------------|-------------|
| Review/Approval | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 07 FEB 2018 |