



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog #/Size:** R32AC-Series

**Lot Number:** W0017258S

**Expiration Date:** 31 DEC 2019

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

Total Volume: 100 µL

Concentration: 500 µg/mL

L/P: 8.0

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        |
|-----------------|--------------|-------------------|---------------------|-------------|
| Review/Approval | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 02 OCT 2017 |