



**CERTIFICATE OF ANALYSIS**

Page 1 of 1

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

**Catalog #/Size:** R32AB-Series

**Lot Number:** W0017282S

**Expiration Date:** 31 JUL 2019

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

Total Volume: 100 µL / 400 µL

Concentration: 500 µg/mL

L/P: 5.0

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.**

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	21 SEP 2017