



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** SULFO-TAG™ Anti-PY20 Antibody

**Catalog #/Size:** R32AP-1

**Lot Number:** W0016439S2

**Expiration Date:** 31 AUG 2019

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

Total Volume: 400 µL

Concentration: 500 µg/mL

L/P: 7.1

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 JUN 2018