



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** SULFO-TAG™ Anti-Hu/NHP IgG Antibody

**Catalog #/Size:** D20JL - Series

**Lot Number:** W0016994S

**Expiration Date:** 30 SEP 2019

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

Total Volume: 200 µL

Concentration: 500 µg/mL

L/P: 2.8

Diluent: Diluent 100

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>	22 MAY 2017