



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** D20JJ-6

**Lot Number:** W0019824S

**Expiration Date:** 30 SEP 2023

Total Volume: 200 µL


Concentration: 500 µg/mL

L/P: 3.4

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>	Karen Hamilla	Quality		08 MAY 2020