



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog #/Size: D20JP - Series

Lot Number: W0017441S

Expiration Date: 29 FEB 2020

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

Total Volume: 200 µL

Concentration: 500 µg/mL

L/P: 10.5

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>	23 MAY 2018