



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: D20JP-6

Lot Number: W0019928S

Expiration Date: 31 DEC 2023

Total Volume: 200 µL


Concentration: 500 µg/mL

L/P: 12.1

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.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		11 SEP 2020