



CERTIFICATE OF ANALYSIS

Product Description: SULFO-TAG™ Anti-Hu/NHP IgM Antibody
Catalog Number: D20JP-Series
Lot Number: W0022130S
Expiration Date: 31 DEC 2027
Storage Temperature: 2-8°C

Size	Total Volume	Concentration	L/P	Final Buffer
100µg	200µL	500µg/mL	13.3	Diluent 100 w/ Azide

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		26 SEP 2024