



CERTIFICATE OF ANALYSIS

Product Description: SULFO-TAG Anti-Hu/NHP IgA Antibody
Catalog Number: D20JJ-6
Lot Number: W0023007S
Expiration Date: 30 SEP 2029
Storage Temperature: 2-8°C

Size	Total Volume	Concentration	L/P	Final Buffer
100 µg	200 µL	500 µg/mL	3.7	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		10 APR 2026