



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: D20JJ-Series

Lot Number: W0020728S3

Expiration Date: 30 JUN 2026

Storage Temperature: 2-8°C

Size	Total Volume	Concentration	L/P	Final Buffer
100 µg	200 µL	500 µg/mL	2.8	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		31 JAN 2023