



CERTIFICATE OF ANALYSIS

Product Description: SULFO-TAG™ Streptavidin (200 µg)
Catalog Number: R32AD-Series
Lot Number: W0022423S2
Expiration Date: 30 NOV 2028
Storage Temperature: 2-8°C

Size	Total Volume	Concentration	L/P	Final Buffer
200 µg	400 µL	500 µg/mL	12.4	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 JUL 2025