



CERTIFICATE OF ANALYSIS

Product Description: SULFO-TAG™ Streptavidin (200 µg)
Catalog Number: R32AD-Series
Lot Number: W0022244S
Expiration Date: 31 MAR 2028
Storage Temperature: 2-8°C

Size	Total Volume	Concentration	L/P	Final Buffer
200 µg	400µL	500 µg/mL	10.2	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		14 NOV 2024