



CERTIFICATE OF ANALYSIS

Product Description: SULFO-TAG™ Streptavidin
Catalog Number: R32AD-Series
Lot Number: W0022423S
Expiration Date: 31 JUL 2028
Storage Temperature: 2-8°C

Size	Total Volume	Concentration	L/P	Final Buffer
50 µg	100 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		02 APR 2025