



CERTIFICATE OF ANALYSIS

Product Description: SULFO-TAG Goat Anti-Rabbit Antibody
Catalog Number: R32AB-1
Lot Number: W0022593S3
Expiration Date: 31 DEC 2029
Storage Temperature: 2-8°C

Size	Total Volume	Concentration	L/P	Final Buffer
200µg	400 µL	500 µg/mL	5.9	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		25 AUG 2026