



CERTIFICATE OF ANALYSIS

Product Description: SULFO-TAG Goat Anti-Rat Antibody
Catalog Number: R32AH-5
Lot Number: W0022306S3
Expiration Date: 31 AUG 2029
Storage Temperature: 2-8°C

Size	Total Volume	Concentration	L/P	Final Buffer
50µg	100 µL	500 µg/mL	6.7	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