



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: SULFO-TAG™ Anti-PY20 Antibody

Catalog Number: R32AP-Series

Lot Number: W0019093S4

Expiration Date: 31 DEC 2025

Storage Temperature: 2-8°C

Size	Total Volume	Concentration	L/P	Final Buffer
200 µg	400 µL	500 µg/mL	7.0	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		31 AUG 2022