



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: SULFO-TAG™ Human/NHP Lambda Detection Antibody
Catalog Number: D20QG-6
Lot Number: W0018775S
Expiration Date: 30 NOV 2022

Total Volume: 200 µL
Concentration: 500 µg/mL
L/P: 10.5
Diluent: Diluent 100 w/ Azide
Storage conditions: 2-8°C

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Carissa Hubbard	Quality		28 AUG 2019