



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: SULFO-TAG™ Anti-PA Antibody

Catalog #/Size: D20XX-1

Lot Number: W0016500-EM170630

Expiration Date: 30 NOV 2019

Statement: This product was manufactured and tested according to approved MSD documents.

Total Volume: 60 µL / 300 µL

Concentration: 50X

L/P: 13.1

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	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	15 SEP 2017