



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: SULFO-TAG™ Anti-Human IgG (2A11) Antibody

Catalog #/Size: S52AC-1 / 1 vial

Lot Number: W0018415

Date of Manufacture: 11 DEC 2018

Storage Temperature: 2-8°C

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

Summary of Conjugation:

Raw Material Lot #	Final Conc. (µg/mL)	L/P	% Yield	Final Buffer	Final Volume (µL)
15/03-G1-A11	1917	6.3	62.1	PBS + 0.05% Azide	1620

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>	17 JAN 2019