



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: SULFO-TAG™ Anti-C-Peptide

Catalog #/Size: S52AC-1/1 vial

Lot Number: W0017460

Date of Manufacture: 18 SEP 2017

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)
N/A	1913	9.1	55.7%	PBS w/ 0.05% Azide	640

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>	27 SEP 2017