



CERTIFICATE OF ANALYSIS

Page 1 of 1

COA – 05338

Product Description: SULFO-TAG™ Anti-pSTAT-6 (Tyr641)

Catalog #/ Size: S52AC-1/1 vial

Lot Number: W0015625

Storage Condition: 2-8°C

Date of Manufacture: 27 MAR 2015



Expiration Date: 31 MAR 2017

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)
0079S0315	895	10.8	81.90%	PBS + 0.05% Azide	850

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

	Name	Function	Signature	Date
Preparer	Meredith Hutchinson	Inventory Technician		01 APR 2015
Review/Approval	Barbara Twum	Manufacturing Lead		01 APR 2015