



## CERTIFICATE OF ANALYSIS

Page 1 of 1

COA – 04762

**Product Description:** SULFO-TAG™ OSM Antibody

**Catalog #/ Size:** S52AC-1/1 vial

**Lot Number:** W0015408

**Storage Condition:** 2-8°C

**Date of Manufacture:** 16 SEP 2014



**Expiration Date:** 30 SEP 2016

**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	1344	5.6	86.00%	PBS + 0.05% Azide	5823

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

	Name	Function	Signature	Date
<b>Preparer</b>	Meredith Hutchinson	Inventory Technician		19 SEP 2014
<b>Review/Approval</b>	Barbara Twum	Manufacturing Lead		19 SEP 2014