



RFE MEMO

Page 1 of 1

Product Description(s): SULFO-TAG™ Streptavidin

Catalog Number(s) / Size(s): R32AD-Series

Lot Number: W0016597S

Based on additional testing, the original date of expiration for this detection antibody has been extended by 12 months. The expiration date of W0016597S has been extended to 28 Feb 2020.

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar		18 APR 2019

This memo supersedes all other statements regarding date of expiration for this product.



CERTIFICATE OF ANALYSIS

COA - 06378

Product Description: SULFO-TAG™ Streptavidin

Catalog #/Size: R32AD-1/200 µg; R32AD-5/ 50 µg

Lot Number: W0016597S

Expiration Date: 28 FEB 2019

Statement: This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

Volume: 400 µL; 100 µL

Concentration: 500 µg/mL

L/P: 11.1

Diluent: 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
Preparer	Meredith Hutchinson	Production Control		09 DEC 2016
Review/Approval	Dawn Stollar	Quality Assurance		09 DEC 2016