



RFE MEMO


Page 1 of 1

Product Description(s): SULFO-TAG Streptavidin

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

Lot Number: W0017855S

Based on additional testing, the original date of expiration for this detection antibody has been extended by 17 months. The expiration date of W0017855S has been extended to 30 SEP 2021

Function	Name	Signature	Date
Quality	Carissa B. Hubbard		07 APR 2021

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



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: SULFO-TAG™ Streptavidin

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

Lot Number: W0017855S

Expiration Date: 30 APR 2020

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

Total Volume: 400 µL; 100 µL

Concentration: 500 µg/mL

L/P: 10.2

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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	16 APR 2018