



## RFE MEMO

Page 1 of 1

**Product Description(s):** MSD Sulfo-TAG™ Amine 500 nmol

**Catalog Number(s) / Size(s):** R90AB-1 / 500 nmol

**Lot Number:** S0130002

Based on additional testing, the original date of expiration for this product has been extended by 58 months. The expiration date of S0130002 has been extended to 19 APR 2019.

Function	Name	Signature	Date
Quality Assurance	Dawn Stollar	<i>Dawn Stollar</i>	19 APR 2018

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



## CERTIFICATE OF ANALYSIS

COA - 11159

**Product Description:** MSD Sulfo-TAG Amine

**Catalog #/Size:** R90AB-1/500 nmol

**Lot Number:** S0130002

**Expiration Date:** 31 MAY 2014

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

Purity (by HPLC): > 90%

Quantity: 100% to 120% of nominal

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

	Name	Function	Signature	Date
Preparer	Elise Garrido	Reagent Manufacturing	<i>Elise Garrido</i>	07 FEB 2013
Review/Approval	Mike Eager	Quality Control	<i>Mike Eager</i>	07 FEB 2013