




EXTENSION MEMO

Page 1 of 1

Product Description(s): SULFO-TAG™ Amine
Catalog Number(s) / Size(s): R90AB-1 / 500 nmol
Lot Number: S0130004

Based on additional testing, the original date of expiration for this reagent has been extended by 29 months. The expiration date of S0130004 has been extended to 30 NOV 2024.

Function	Name	Signature	Date
Quality	Karen Hamilla		05 DEC 2023

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



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: SULFO-TAG™ Amine

Catalog Number: R90AB-1 / 500 nmol

Lot Number: S0130004


Expiration Date: 31 MAY 2022

Storage Temperature: ≤ -70°C

Statement: This product was manufactured according to MSD documents. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

Purity specification (by HPLC): > 90%

Quantity: 100% to 120% of nominal

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		03 JUN 2020