




Extension Memo

Product Description: Sulfo-TAG Iodoacetamide, 500 nmol
Lot Number: S0150012
New Expiration Date: 31 JAN 2027
Catalog Number(s): R90AA-1

Based on additional testing, the original date of expiration for this reagent has been extended by 12 months.

Function	Name	Signature	Date
Quality	Karen Hamilla		22 Dec 2025




CERTIFICATE OF ANALYSIS

Product Description: Sulfo-TAG Iodoacetamide, 500 nmol
Catalog Number: R90AA-1
Lot Number: S0150012
Expiration Date 31 JAN 2026

Functional Results:

	Specification	Measured Average
Purity (by HPLC):	$\geq 90\%$	99%
Quantity	100% to 120% of nominal	113%

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		28 FEB 2024