



## 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):	≥ 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
<b>Review/Approval</b>	Karen Hamilla	Quality		28 FEB 2024