


**CERTIFICATE OF ANALYSIS**

Product Description: SULFO-TAG™ Iodoacetamide, 500 nmol
Catalog Number: R90AA-1
Lot Number: S0150014
Expiration Date 31 OCT 2027

Functional Results:

	Specification	Measured Average
Purity (by HPLC)	≥ 90%	96%
Quantity	100% to 120% of nominal	117%

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		03 DEC 2025