



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Sulfo-TAG Iodoacetamide, 500 nmol

Catalog Number: R90AA-1

Lot Number: S0150011

Expiration Date: 31 JUL 2024

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


Purity specification (by HPLC): > 90%

Measured Average Purity: 100%

Quantity: 100% to 120% of nominal

Measured Average Quantity: 112%

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		02 SEP 2022