



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** Sulfo-TAG Iodoacetamide, 500 nmol

**Catalog Number:** R90AA-1

**Lot Number:** S0150010

**Expiration Date:** 30 SEP 2023

**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: 98%

Quantity: 100% to 120% of nominal

Measured Average Quantity: 106%

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		08 DEC 2021