



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Sulfo-TAG Iodoacetamide, 500 nmol
Catalog #/Size: R90AA-1 / 500 nmol
Lot Number: S0150008
Expiration Date: 30 APR 2021

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: 114%

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

	Name	Function	Signature	Date
Review/Approval	Carissa Hubbard	Quality		17 JUN 2019