



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Sulfo-TAG Iodoacetamide, 500 nmol
Catalog #/Size: R90AA-1 / 500 nmol
Lot Number: S0150007
Expiration Date: 31 JUL 2019

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

Purity (by HPLC): > 90%
Quantity: 100% to 120% of nominal

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

	Name	Function	Signature	Date
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	29AUG2017