



CERTIFICATE OF ANALYSIS

COA - 06694

Product Description: GOLD Sulfo-TAG NHS Ester

Catalog #/Size: R91AO-1

Lot Number: S0080070

Expiration Date: 31 MAR 2019

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

Purity specification (by HPLC): > 95%

Measured Average Purity: 98%

Quantity: 100% to 120% of nominal

Measured Average Quantity: 117%

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

	Name	Function	Signature	Date
Preparer	Morgan Manogue	Quality Control	<i>Morgan Manogue</i>	25 MAR 2016
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	25 MAR 2016