



CERTIFICATE OF ANALYSIS

COA - 07608

Product Description: MSD GOLD Sulfo-TAG NHS Ester

Catalog #/Size: R91AO-2

Lot Number: S0100039

Expiration Date: 30 NOV 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: 99%

Quantity: 100% to 120% of nominal

Measured Average Quantity: 113%

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

	Name	Function	Signature	Date
Preparer	Xiomara Talavera	Quality Control	<i>X Talavera</i>	09 DEC 2016
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	09 DEC 2016