



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD GOLD SULFO-TAG NHS-Ester

Catalog #/Size: R91AO-1 / 150 nmol

Lot Number: S0080097

Expiration Date: 31 DEC 2020

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: 112%

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	11 DEC 2017