



CERTIFICATE OF ANALYSIS

COA - 07582

Product Description: MSD GOLD SULFO-TAG NHS-Ester 150 nmol

Catalog #/Size: R91AO-1

Lot Number: S0080080

Expiration Date: 31 OCT 2019



Statement: Purity specification (by HPLC): > 95%

Measured Average Purity: 97.6%

Quantity: 100% to 120% of nominal

Measured Average Quantity: 116%

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

	Name	Function	Signature	Date
Preparer	Carissa Hubbard	Quality Control		05 DEC 2016
Review/Approval	Dawn Stollar	Quality Assurance		06 DEC 2016