



## CERTIFICATE OF ANALYSIS

COA - 07581

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

**Catalog #/Size:** R91AO-1

**Lot Number:** S0080079

**Expiration Date:** 31 OCT 2019



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

Measured Average Purity: 96.9%

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.**

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