



## CERTIFICATE OF ANALYSIS

COA - 07607

**Product Description:** MSD GOLD Sulfo-TAG NHS Ester

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

**Lot Number:** S0080081

**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: 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.**

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Preparer</b>	Xiomara Talavera	Quality Control	<i>Xiomara Talavera</i>	07 DEC 2016
<b>Review/Approval</b>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	08 DEC 2016