



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Quantity**

**Product Description:** MSD GOLD™ SULFO-TAG™ NHS-Ester

**Catalog #/Size:** R91AO-1 / 150 nmol

**Lot Number:** S0080109

**Expiration Date:** 31 JAN 2022

**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: 107%

**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>Review/Approval</b>	Dawn Stollar	Quality Assurance		30 APR 2019