



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog #/Size:** R91AO-2 / 2  $\mu$ mol

**Lot Number:** S0100046

**Expiration Date:** 31 AUG 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: 116%

**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	<i>Dawn Stollar</i>	11 SEP 2017