



**CERTIFICATE OF ANALYSIS**

Page 1 of 1

**Product Description:** MSD GOLD™ SULFO-TAG™ NHS-Ester  
**Catalog Number:** R91AO-1 /150 nmol  
**Lot Number:** S0080141  
**Expiration Date:** 31 MAY 2024

**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: 112%

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>	Karen Hamilla	Quality		07 JUN 2021