



**CERTIFICATE OF ANALYSIS**

**Product Description:** MSD GOLD™ SULFO-TAG NHS-Ester 150 nmol  
**Catalog Number:** R91AO-1  
**Lot Number:** S0080172  
**Expiration Date** 31 MAY 2027

**Functional Results:**

	<b>Specification</b>	<b>Measured Average</b>
Purity (by HPLC):	≥ 95%	98%
Quantity	100% to 120% of nominal	113%

This product was manufactured and tested according to MSD documents. This lot meets MSD’s specifications. This 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 2024