



**CERTIFICATE OF ANALYSIS**

**Product Description:** MSD GOLD™ SULFO-TAG NHS-Ester 2 μmol  
**Catalog Number:** R91AO-2  
**Lot Number:** S0100140  
**Expiration Date** 31 MAR 2029

**Functional Results:**

	<b>Specification</b>	<b>Measured Average</b>
Purity (by HPLC)	≥ 95%	99%
Quantity	100% to 120% of nominal	107%

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		01 MAY 2026