




## CERTIFICATE OF ANALYSIS

**Product Description:** MSD GOLD™ SULFO-TAG NHS-Ester 2 µmol  
**Catalog Number:** R91AO-2  
**Lot Number:** S0100130  
**Expiration Date:** 31 OCT 2027

**Functional Results:**

	Specification	Measured Average
Purity (by HPLC):	≥ 95%	97%
Quantity	100% to 120% of nominal	115%

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.

	Name	Function	Signature	Date
<b>Review/Approval</b>	Hamilla, Karen	Quality		10 DEC 2024