




## CERTIFICATE OF ANALYSIS

**Product Description:** MSD GOLD™ SULFO-TAG NHS-Ester 2 µmol  
**Catalog Number:** R91AO-2  
**Lot Number:** S0100136  
**Expiration Date:** 30 JUN 2028

**Functional Results:**

	Specification	Measured Average
Purity (by HPLC)	≥ 95%	99%
Quantity	100% to 120% of nominal	110%

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>	Karen Hamilla	Quality		06 AUG 2025