



## CERTIFICATE OF ANALYSIS

**Product Description:** MSD GOLD™ SULFO-TAG NHS-Ester 150 nmol  
**Catalog Number:** R91AO-1  
**Lot Number:** S0080171  
**Expiration Date** 31 MAR 2027

**Functional Results:**

	Specification	Measured Average
Purity (by HPLC):	≥ 95%	98%
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		23 MAY 2024