



CERTIFICATE OF ANALYSIS

Product Description: MSD GOLD™ SULFO-TAG NHS-Ester 150 nmol
Catalog Number: R91AO-1
Lot Number: S0080178
Expiration Date 31 JAN 2028

Functional Results:

	Specification	Measured Average
Purity (by HPLC)	≥ 95%	96%
Quantity	100% to 120% of nominal	114%

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
Review/Approval	Karen Hamilla	Quality		20 FEB 2025