



CERTIFICATE OF ANALYSIS

Product Description: MSD GOLD SULFO-TAG NHS-Ester 150 nmol
Catalog Number: R91AO-1
Lot Number: S0080187
Expiration Date 31 MAR 2029

Functional Results:

	Specification	Measured Average
Purity (by HPLC)	≥ 95%	97%
Quantity	100% to 120% of nominal	113%

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		22 APR 2026