



CERTIFICATE OF ANALYSIS

Product Description: MSD GOLD™ SULFO-TAG NHS-Ester 150 nmol
Catalog Number: R91AO-1
Lot Number: S0080183
Expiration Date 31 JUL 2028

Functional Results:

	Specification	Measured Average
Purity (by HPLC)	$\geq 95\%$	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
Review/Approval	Michelle Likiam	Quality		26 AUG 2025