



CERTIFICATE OF ANALYSIS

Product Description: MSD GOLD™ SULFO-TAG NHS-Ester 150 nmol
Catalog Number: R91AO-1
Lot Number: S0080169
Expiration Date: 30 NOV 2026

Functional Results:

	Specification	Measured Average
Purity (by HPLC):	≥ 95%	99%
Quantity	100% to 120% of nominal	117%

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		02 JAN 2024