



CERTIFICATE OF ANALYSIS

Product Description: MSD GOLD™ SULFO-TAG NHS-Ester 2 µmol
Catalog Number: R91AO-2
Lot Number: S0100124
Expiration Date 30 APR 2027

Functional Results:

	Specification	Measured Average
Purity (by HPLC):	≥ 90%	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
Review/Approval	Karen Hamilla	Quality		15 MAY 2024