


**CERTIFICATE OF ANALYSIS**

Product Description: MSD GOLD™ SULFO-TAG NHS-Ester 2 µmol
Catalog Number: R91AO-2
Lot Number: S0100135
Expiration Date 30 JUN 2028

Functional Results:

	Specification	Measured Average
Purity (by HPLC)	≥ 95%	99%
Quantity	100% to 120% of nominal	109%

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		31 JUL 2025