



CERTIFICATE OF ANALYSIS

Product Description: MSD GOLD™ SULFO-TAG NHS-Ester
Catalog Number: R91AO-1
Lot Number: S0080189
Expiration Date: 30 JUN 2029

Functional Results:

	Specification	Measured Average
Purity (by HPLC)	≥ 95%	99%
Quantity	100% to 120% of nominal	116%

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		17 AUG 2026