



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD GOLD™ SULFO-TAG™ NHS-Ester (lyophilized)

Catalog Number: R91AO-2 / 2 µmol

Lot Number: S0100104

Expiration Date: 31 JUL 2025

Statement: This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.


Purity specification (by HPLC): > 95%

Measured Average Purity: 99%

Quantity: 100% to 120% of nominal

Measured Average Quantity: 110%

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		09 SEP 2022