



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD GOLD™ SULFO-TAG™ NHS-Ester
Catalog #/Size: R91AO-2 / 2 μmol
Lot Number: S0100097
Expiration Date: 31 JAN 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: 111%

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		16 MAR 2022