



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD GOLD™ SULFO-TAG™ NHS-Ester
Catalog Number: R91AO-1 / 150 nmol
Lot Number: S0080129
Expiration Date: 28 FEB 2023

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: 98%

Quantity: 100% to 120% of nominal
Measured Average Quantity: 112%

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		27 FEB 2020