



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog Number:** R91AO-1 /150 nmol

**Lot Number:** S0080158

**Expiration Date:** 30 NOV 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: 98%

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.

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Review/Approval</b>	Karen Hamilla	Quality		03 JAN 2023