



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

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

**Lot Number:** S0080145

**Expiration Date:** 30 NOV 2024

**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: 114%

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

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		03 JAN 2022