



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog #/Size:** R91AO-2 / 2 µmol

**Lot Number:** S0100065

**Expiration Date:** 31 MAR 2022

**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: 114%

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality		18 APR 2019