



## CERTIFICATE OF ANALYSIS

Page 1 of 1

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

**Catalog #/Size:** R91AO-1 / 150 nmol

**Lot Number:** S0080095

**Expiration Date:** 30 SEP 2020

**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: 109%

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

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	09 OCT 2017