



**CERTIFICATE OF ANALYSIS**

Page 1 of 1

**Product Description:** MSD GOLD SULFO-TAG NHS-Ester  
**Catalog #/Size:** R91AO-1  
**Lot Number:** S0080090  
**Expiration Date:** 31 MAY 2020

Purity specification (by HPLC): > 95%  
Measured Average Purity: 97%

Quantity: 100% to 120% of nominal  
Measured Average Quantity: 116%

**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>	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	23 MAY 2017