



**CERTIFICATE OF ANALYSIS**

Page 1 of 1

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

**Catalog #/Size:** R91AO-2

**Lot Number:** S0100043

**Expiration Date:** 31 MAY 2020

Purity specification (by HPLC): > 95%

Measured Average Purity: 98%

Quantity: 100% to 120% of nominal

Measured Average Quantity: 108%

**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>	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	15JUN2017