



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD GOLD™ SULFO-TAG NHS-Ester
Catalog #/Size: R91AO-1
Lot Number: S0080089
Expiration Date: 31 MAY 2020

Purity specification (by HPLC): > 95%
Measured Average Purity: 98%

Quantity: 100% to 120% of nominal
Measured Average Quantity: 117%

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

	Name	Function	Signature	Date
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	31MAY2017