



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD GOLD Sulfo-TAG NHS Ester
Catalog #/Size: R91AO-1
Lot Number: S0080088
Expiration Date: 29 FEB 2020

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

Quantity: 100% to 120% of nominal
Measured Average Quantity: 118%

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>	30 MAR 2017