



CERTIFICATE OF ANALYSIS

COA - 07663

Product Description: MSD GOLD Sulfo-TAG NHS Ester

Catalog #/Size: R91AO-1

Lot Number: S0080082

Expiration Date: 31 DEC 2019

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: 98%

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.

| | Name | Function | Signature | Date |
|------------------------|------------------|-------------------|---------------------|-------------|
| Preparer | Xiomara Talavera | Quality Control | <i>Xiomara</i> | 20 DEC 2016 |
| Review/Approval | Dawn Stollar | Quality Assurance | <i>Dawn Stollar</i> | 20 DEC 2016 |