



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog #/Size: R91AO-1

Lot Number: S0080101

Expiration Date: 28 FEB 2021

Purity specification (by HPLC): > 95%

Measured Average Purity: 98%

Quantity: 100% to 120% of nominal

Measured Average Quantity: 111%

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

| | Name | Function | Signature | Date |
|-----------------|-----------|-------------------|--------------|-----------|
| Review/Approval | Emily Lee | Quality Assurance | <i>E Lee</i> | 28FEB2018 |