




CERTIFICATE OF ANALYSIS

Product Description: MSD GOLD™ SULFO-TAG NHS-Ester 2 μmol
Catalog Number: R91AO-2
Lot Number: S0100133
Expiration Date 31 MAR 2028

Functional Results:

| | Specification | Measured Average |
|------------------|-------------------------|------------------|
| Purity (by HPLC) | ≥ 95% | 99% |
| Quantity | 100% to 120% of nominal | 108% |

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

| | Name | Function | Signature | Date |
|-----------------|---------------|----------|---|-------------|
| Review/Approval | Karen Hamilla | Quality |  | 22 MAY 2025 |