


**CERTIFICATE OF ANALYSIS**

**Product Description:** MSD GOLD SULFO-TAG NHS-Ester 150 nmol  
**Catalog Number:** R91AO-1  
**Lot Number:** S0080185  
**Expiration Date** 30 NOV 2028

**Functional Results:**

|                  | Specification           | Measured<br>Average |
|------------------|-------------------------|---------------------|
| Purity (by HPLC) | $\geq 95\%$             | 99%                 |
| Quantity         | 100% to 120% of nominal | 106%                |

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        |
|------------------------|---------------|----------|---|-------------|
| <b>Review/Approval</b> | Karen Hamilla | Quality  |  | 05 JAN 2026 |