


**CERTIFICATE OF ANALYSIS**

Product Description: MSD GOLD Read Buffer A
Catalog Number: R92TG-Series
Lot Number: Y0600068
Expiration Date 30 SEP 2028

Statement: MSD GOLD Read Buffer A is verified to ensure that consistent signals are obtained across lots. Details on the quality control testing of MSD GOLD Read Buffer A are provided in the product insert available at www.mesoscale.com.

| Concentration of BTI | Specification Metric | Allowable Range | Pass/Fail? |
|-------------------------|------------------------------------|------------------|------------|
| 2.00 nM BTI (0.10 pmol) | Test : Reference Signal Ratio | 90 - 110 % | Pass |
| | Coefficient of Variation | $\leq 10 \%$ | Pass |
| 0 nM BTI (0 pmol) | Test - Reference Signal Difference | ± 10 counts | Pass |
| | Standard Deviation | ≤ 15 counts | Pass |

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

| | Name | Function | Signature | Date |
|-----------------|---------------|----------|--|-------------|
| Review/Approval | Karen Hamilla | Quality |  | 14 NOV 2025 |