



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD GOLD™ Read Buffer A

**Catalog Number:** R92TG-Series

**Lot Number:** Y0600039

**Expiration Date:** 31 DEC 2025

**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](http://www.mesoscale.com).

MSD GOLD™ Read Buffer A was tested using SULFO-TAG™ Conjugated Biotinylated IgG (BTI) lot # R0010769 and MSD GOLD™ 96-Well Streptavidin Plate Z0021899. Specifications are BTI lot specific.

| Concentration of BTI    | Mean ECL Spec and Tolerance Range | Intraplate CV or SD spec | Pass/Fail |
|-------------------------|-----------------------------------|--------------------------|-----------|
| 2.00 nM BTI (0.10 pmol) | 17741 - 24002 counts              | 10 %                     | PASS      |
| 0 nM BTI (0 pmol)       | 48 - 68 counts                    | 15 counts                | PASS      |

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

|                 | Name          | Function | Signature | Date        |
|-----------------|---------------|----------|-----------|-------------|
| Review/Approval | Karen Hamilla | Quality  |           | 17 FEB 2023 |