



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD GOLD™ Read Buffer B

Catalog Number: R60AM-Series

Lot Number: Y0800026

Expiration Date: 31 JUL 2024

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

MSD GOLD™ Read Buffer B was tested using SULFO-TAG™ Conjugated Biotinylated IgG (BTI) lot # R0010769 and MSD GOLD™ 96-well Streptavidin SECTOR Plate Z0021725. 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) | 17042 - 23057 counts | 10 % | Pass |
| 0 nM BTI (0 pmol) | 24 - 44 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 | | 24 MAR 2022 |