



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MSD GOLD™ 96-Well Small Spot Streptavidin SECTOR® Plate

Catalog Number: L45SA-Series

Lot Number: Z0022062

Expiration Date: 31 AUG 2025

The above plate lot was tested according to MSD's procedures and passed all specifications. Plates were tested using SULFO-TAG™ labeled biotinylated IgG (BTI).

Titration Results:

| Metric | Specification | Result |
|---|---------------|--------|
| 0.075 pmol BTI, Ratio of signal to reference lot signal | 85-115% | Pass |
| 0.050 pmol BTI, Ratio of signal to reference lot signal | 85-115% | Pass |
| 0.025 pmol BTI, Ratio of signal to reference lot signal | 85-115% | Pass |
| Max Background Signal (counts) | ≤ 100 | Pass |
| Standard Deviation of Background Signal (counts) | ≤ 15 | Pass |

Uniformity Results:

Plates sampled from the above plate lot were tested by adding 0.050 pmol BTI to each well.

| Metric | Specification | Result |
|---|---------------|--------|
| Mean Intraplate CV | ≤ 6.0% | Pass |
| Intraplate CVs for 91.5% of plates | ≤ 8.0% | Pass |
| Number of plates with Intraplate CV > 12% | 0 Plates | Pass |
| Interplate CV | ≤ 8 % | 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 | | 06 APR 2023 |