

**CERTIFICATE OF ANALYSIS**

Product Description: MSD GOLD™ 96-well Streptavidin SECTOR Plate
Catalog Number: L15SA-Series
Lot Number: Z0022309
Expiration Date: 30 NOV 2026

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 |
|---|---------------|--------|
| 0.300 pmol BTI, Ratio of signal to reference lot signal | 85-115% | Pass |
| 0.200 pmol BTI, Ratio of signal to reference lot signal | 85-115% | Pass |
| 0.100 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 0.200 pmol BTI to each well.

| Metric | Specification | Result |
|---|---------------|--------|
| Average Intraplate CV | ≤ 6.0% | Pass |
| Percentage of plates with CV ≤ 8.0% | ≥ 91.5% | Pass |
| Number of plates with Intraplate CV > 12% | 0 Plates | Pass |
| Interplate CV | ≤ 8.0% | 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 | | 09 AUG 2024 |