

**CERTIFICATE OF ANALYSIS**

Product Description: MSD GOLD 96-Well Small Spot Streptavidin SECTOR Plate
Catalog Number: L45SA-Series
Lot Number: Z0022429
Expiration Date: 30 SEP 2027

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 | 80-120% | Pass |
| 0.050 pmol BTI, Ratio of signal to reference lot signal | 80-120% | Pass |
| 0.025 pmol BTI, Ratio of signal to reference lot signal | 80-120% | 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 |
|--|---------------|--------|
| Average Intraplate CV | $\leq 6.0\%$ | Pass |
| Percentage of plates with $CV \leq 8.0\%$ | $\geq 91.5\%$ | Pass |
| Number of plates with Intraplate $CV > 12\%$ | 0 Plates | Pass |
| Interplate CV | $\leq 8.0\%$ | 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 |  | 15 MAY 2025 |