



## CERTIFICATE OF ANALYSIS

**Product Description:** MSD GOLD™ 96-well Small Spot Streptavidin SECTOR® Plate  
**Catalog Number:** L45SA-Series  
**Lot Number:** Z0022245  
**Expiration Date:** 30 JUN 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 |
|---|---------------|--------|
| 0.075 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 |
|---|---------------|--------|
| 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        |
|------------------------|---------------|----------|-----------|-------------|
| <b>Review/Approval</b> | Karen Hamilla | Quality  |           | 22 JAN 2024 |