



## CERTIFICATE OF ANALYSIS

**Product Description:** MSD GOLD 96-Well Small Spot Streptavidin SECTOR Plate  
**Catalog Number:** L45SA-Series  
**Lot Number:** Z0022521  
**Expiration Date:** 31 DEC 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)                          | $\leq 100$    | Pass   |
| Standard Deviation of Background Signal (counts)        | $\leq 15$     | 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 | Michelle Likiam | Quality  |  | 31 JUL 2025 |