

**CERTIFICATE OF ANALYSIS**

**Product Description:** MSD GOLD 96-Well 1-Spot Streptavidin SECTOR Plate  
**Catalog Number:** L15SA-Series  
**Lot Number:** Z0022580  
**Expiration Date:** 31 MAR 2028

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)                          | $\leq 100$    | Pass   |
| Standard Deviation of Background Signal (counts)        | $\leq 15$     | Pass   |

**Uniformity Results:**

Plates sampled from the above plate lot were tested by adding 0.200 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        |
|------------------------|---------------|----------|--------------------------------------------------------------------------------------|-------------|
| <b>Review/Approval</b> | Karen Hamilla | Quality  |  | 20 NOV 2025 |