



**Product Description:** MSD GOLD<sup>TM</sup> 96-well Small Spot Streptavidin SECTOR Plate

Catalog Number:L45SA-SeriesLot Number:Z0022356Expiration Date:28 FEB 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                     | ≤ 6.0%        | Pass   |
| Percentage of plates with $CV \le 8.0\%$  | ≥ 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  | Haven Hamille | 26 SEP 2024 |

COA-21169-A QA-FM-266-E