



## CERTIFICATE OF ANALYSIS

Page 1 of 1

**Product Description:** MSD GOLD™ 96-well Streptavidin SECTOR® Plate

**Catalog Number:** L15SA-Series

**Lot Number:** Z0021722

**Expiration Date:** 30 APR 2023

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 |
|--|---------------|--------|
| Mean Intraplate CV                           | $\leq 6.0\%$  | Pass   |
| Intraplate CVs for 91.5% of plates           | $\leq 8.0\%$  | Pass   |
| Number of plates with Intraplate CV $> 12\%$ | 0 Plates      | Pass   |
| Interplate CV                                | $\leq 8\%$    | 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  |  | 12 NOV 2020 |