

**CERTIFICATE OF ANALYSIS**

**Product Description:** MSD GOLD™ 96-well Small Spot Streptavidin SECTOR® Plate  
**Catalog Number:** L45SA-Series  
**Lot Number:** Z0022157  
**Expiration Date:** 31 JAN 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	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 ≤ 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		21 AUG 2023