

**CERTIFICATE OF ANALYSIS**

Product Description: MSD GOLD 96-Well Small Spot Streptavidin SECTOR Plate
Catalog Number: L45SA-Series
Lot Number: Z0022708
Expiration Date: 31 DEC 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
0.075 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	$\leq 6.0\%$	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	Karen Hamilla	Quality		14 AUG 2026