



CERTIFICATE OF ANALYSIS

Product Description: MSD GOLD 96-Well 1-Spot Streptavidin SECTOR Plate
Catalog Number: L15SA-Series
Lot Number: Z0022529
Expiration Date: 31 JAN 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:

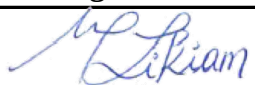
Metric	Specification	Result
0.300 pmol BTI, Ratio of signal to reference lot signal	85-115%	Pass
0.200 pmol BTI, Ratio of signal to reference lot signal	85-115%	Pass
0.100 pmol BTI, Ratio of signal to reference lot signal	85-115%	Pass
Max Background Signal (counts)	≤ 100	Pass
Standard Deviation of Background Signal (counts)	≤ 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
Review/Approval	Michelle Likiam	Quality		26 AUG 2025