

Page 1 of 1

Product Description:	MSD GOLD [™] 96-well Small Spot Streptavidin SECTOR [®] Plate		
Catalog Number:	L45SA-Series		
Lot Number:	Z0022236		
Expiration Date:	31 MAY 2026		

The above plate lot was tested according to MSD's procedures and passed all specifications. Plates were tested using SULFO-TAGTM 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 $CV \le 8.0\%$	≥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	Likiam	02 JAN 2024