



CERTIFICATE OF ANALYSIS

Page 1 of 1

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

Catalog Number: L15SA-Series

Lot Number: Z0021932

Expiration Date: 31 JAN 2025

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
Mean Intraplate CV	≤ 6.0%	Pass
Intraplate CVs for 91.5% of plates	≤ 8.0%	Pass
Number of plates with Intraplate CV > 12%	0 Plates	Pass
Interplate CV	≤ 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		29 AUG 2022