

## **CERTIFICATE OF ANALYSIS**

## Page 1 of 1

**Product Description:** MSD GOLD<sup>TM</sup> 96-Well Streptavidin SECTOR® Plate

Catalog Number: L15SA-Series

**Lot Number:** Z0022023

**Expiration Date:** 30 JUN 2025

The above plate lot was tested according to MSD's procedures and passed all specifications. Plates were tested using SULFO-TAG<sup>TM</sup> 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	Nithya Santhanam	Quality	Mit	09 Feb 2023

COA-17841-A QA-FM-266-A