



## CERTIFICATE OF ANALYSIS: TITRATION PLATES

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

**Catalog #:** L45SA - Series

**Lot Number:** Z0021591

**Date of Manufacture:** 16 Dec 2019

**Expiration Date:** 31 May 2022

**Statement:** This product was manufactured and tested according to approved MSD Standard Operating Procedures.

This product has an approximate binding capacity of 0.075 pmol of biotinylated IgG. For optimal assay results, the amount of biotinylated reagent used in the assay should not exceed this binding capacity.

Titration plates were tested using tagged BTI lot: R0010605

Concentration of SULFO-TAG™ Labeled Biotinylated IgG (BTI)	Referenced Signal	Pass/Fail?
1.5 nM (0.075 pmol)	45472 counts ± 15%	Pass
1 nM (0.05 pmol)	34498 counts ± 15%	Pass
0.5 nM (0.025 pmol)	18826 counts ± 15%	Pass
0 nM (0.0 pmol)	Not to Exceed 100 Counts	Pass

Concentration of SULFO-TAG™ Labeled Biotinylated IgG (BTI)	Intraplate CV or SD	Pass/Fail?
1.5 nM (0.075 pmol)	≤ 10%	Pass
1 nM (0.05 pmol)	≤ 10%	Pass
0.5 nM (0.025 pmol)	≤ 10%	Pass
0 nM (0.0 pmol)	≤ 15 Count SD	Pass



## CERTIFICATE OF ANALYSIS: UNIFORMITY PLATES

**Statement:** This product has an approximate binding capacity of 0.05 pmol of biotinylated IgG. For optimal assay results, the amount of biotinylated reagent used in the assay should not exceed this binding capacity.

Metric	Specification	Pass/Fail?
Mean Intraplate CV	$\leq 6\%$	Pass
Intraplate CVs	CV $\leq 8\%$ for 91.5% of Plates	Pass
Number of plates with intraplate CV $>12\%$	0 Plates	Pass
Interplate CV	$\leq 8\%$	Pass
Plates where signal $> 20\%$ from plate mean occurs in same well on multiple plates	0 Plates	Pass
Wells with signal $> 50\%$ from plate mean	0 Wells	Pass
Median signal for concentric rings, min to max range	$\leq 10.0\%$ for 100% of Plates	Pass
Median signal for columns, min to max range	$\leq 10.0\%$ for 91.5% of Plates	Pass
Median signal for columns, min to max range	$\leq 15.0\%$ for 100% of Plates	Pass
Median signal for rows, min to max range	$\leq 10.0\%$ for 91.5% of Plates	Pass
Median signal for rows, min to max range	$\leq 15.0\%$ for 100% of Plates	Pass

**The above product is intended for research use only. Not for use in diagnostic procedures.**

	Name	Function	Signature	Date
<b>Review/Approval</b>	Karen Hamilla	Quality		13 JAN 2020