



CERTIFICATE OF ANALYSIS: TITRATION PLATES

COA - 8959-B

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

Catalog #: L45SA- Series

Lot Number: Z0021357

Expiration date: 31 MAR 2020

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: R0010578

Concentration of SULFO-TAG™ Labeled Biotinylated IgG (BTI)	Reference Signal	Pass/Fail?
1.5 nM (0.075 pmol)	45,472 counts ± 15%	Pass
1 nM (0.05 pmol)	34,498 counts ± 15%	Pass
0.5 nM (0.025 pmol)	18,826 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)	Not to exceed 15 Counts	Pass



CERTIFICATE OF ANALYSIS: UNIFORMITY PLATES

COA - 8959-B

This product has an approximate binding capacity of 0.075 pmoles 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 CV	$\leq 8\%$ for at least 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 plates	Pass
Median signal for concentric rings, min to max range	$\leq 10\%$ for all plates	Pass
Median signal for columns, min to max range	$\leq 10\%$ for at least 91.5% of plates	Pass
Median signal for columns, min to max range	$\leq 15\%$ for all plates	Pass
Median signal for rows, min to max range	$\leq 10\%$ for at least 91.5% of plates	Pass
Median signal for rows, min to max range	$\leq 15\%$ for all plates	Pass
Number of plates sampled (N)	Lot size dependent; see SOP 3-2000-6P0002	Pass

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

	Name	Function	Signature	Date
Review/Approval	Carissa Hubbard	Quality Control		22 DEC 2017
Review/Approval	Dawn Stollar	Quality Assurance		22 DEC 2017