



CERTIFICATE OF ANALYSIS: TITRATION PLATES

Product Description: MSD GOLD™ 96-Well Small Spot Streptavidin SECTOR Plate
Catalog #: L45SA - Series
Lot Number: Z0021397
Date of Manufacture: 05 APR 2018
Expiration date: 31 AUG 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)	Referenced 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)	≤15 Count SD	Pass



CERTIFICATE OF ANALYSIS: UNIFORMITY PLATES

Statement: 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.

Metric	Specification	Pass/Fail?
Mean Intraplate CV	$\leq 6\%$	Pass
Intraplate CVs	$\leq 8\%$ 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	$\leq 10.0\%$ 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
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	25 APR 2018