



CERTIFICATE OF ANALYSIS: TITRATION PLATES

Product Description: MSD GOLD™ 96-well High Bind Avidin SECTOR® Plate

Catalog #: L15AB - Series

Lot Number: Z0010670

Date of Manufacture: 08 Jul 2019

Expiration Date: 31 Dec 2021

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

This product has an approximate binding capacity of 0.9 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?
18 nM (0.9 pmol)	71164 counts ± 15%	Pass
12 nM (0.6 pmol)	58736 counts ± 15%	Pass
6 nM (0.3 pmol)	35280 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?
18 nM (0.9 pmol)	≤ 10%	Pass
12 nM (0.6 pmol)	≤ 10%	Pass
6 nM (0.3 pmol)	≤ 10%	Pass
0 nM (0.0 pmol)	≤ 15 Count SD	Pass



CERTIFICATE OF ANALYSIS: UNIFORMITY PLATES

Metric	Specification	Pass/Fail?
Mean Intraplate CV	≤ 6%	Pass
Intraplate CVs	CV ≤ 8% for 91.5% of Plates	Pass
Number of plates with intraplate CV >12%	0 Plates	Pass
Interplate CV	≤ 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	≤ 10.0% for 100% of Plates	Pass
Median signal for columns, min to max range	≤ 10.0% for 91.5% of Plates	Pass
Median signal for columns, min to max range	≤ 15.0% for 100% of Plates	Pass
Median signal for rows, min to max range	≤ 10.0% for 91.5% of Plates	Pass
Median signal for rows, min to max range	≤ 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	Karen Hamilla	Quality		02 OCT 2020