



CERTIFICATE OF ANALYSIS: TITRATION PLATES

COA - 05560

Product Description: MSD GOLD™ 96-well Streptavidin SECTOR® PR Plate

Catalog #: L13SA-1 (1 plate); L13SA-2 (5 plate); L13SA-5 (30 plate); L13SA-6 (120 plate);
L13SA-7 (510 plate)

Lot Number: Z0020974

Expiration date: 30Jun2015

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

This product has an approximate binding capacity of 0.3 pmoles 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: R0010419

Concentration of SULFO-TAG™ Labeled Biotinylated IgG (BTI)	Mean ECL spec and tolerance range	Pass/Fail?
6 nM (0.3 pmole)	59,245 counts ± 15%	Pass
4 nM (0.2 pmole)	46,161 counts ± 15%	Pass
2 nM (0.1 pmole)	24,424 counts ± 15%	Pass
0 nM (0.0 pmole)	Not to exceed 200 counts	Pass

Concentration of SULFO-TAG™ Labeled Biotinylated IgG (BTI)	Intraplate CV or SD spec	Pass/Fail?
6 nM (0.3 pmole)	10%	Pass
4 nM (0.2 pmole)	10%	Pass
2 nM (0.1 pmole)	10%	Pass
0 nM (0.0 pmole)	35 count SD	Pass



CERTIFICATE OF ANALYSIS: UNIFORMITY PLATES

COA - 05560

This product has an approximate binding capacity of 0.3 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 CVs	$\leq 8\%$ for at least 92% 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, max to min range	$\leq 10\%$ for all plates	Pass
Median signal for columns, max to min range	$\leq 10\%$ for at least 92% of plates	Pass
Median signal for columns, max to min range	$\leq 15\%$ for all plates	Pass
Median signal for rows, max to min range	$\leq 10\%$ for at least 92% of plates	Pass
Median signal for rows, max to min 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	Heng Xu	Quality Control		15 Jun 2015
Review/Approval	Michael Grossman	Quality Assurance		17 Jun 2015



Meso Scale Diagnostics, L.L.C.

Memo: MSD GOLD 96-well Streptavidin SECTOR[®] PR Plate, Lot Z0020974
June 16, 2015

Based on additional testing, the date of expiration for this plate has been extended. The expiration date of Lot Z0020974 has been extended to 30 Jun 2016.

Function	Name	Signature	Date
Quality Assurance	Michael Grossman		16 Jun 2015

This memo supersedes all other statements of date of expiration for this plate, Lot Z0020974.