



CERTIFICATE OF ANALYSIS

COA - 03578

Product Description: MULTI-ARRAY® 96-well High Bind Avidin Gold Plate

Catalog #/Size: L15AB-1/L11AB-1 (1 plate); L15AB-2/L11AB-2 (5 plates); L15AB-5/L11AB-5 (30 plates); L15AB-6/L11AB-6 (120 plates); L15AB-7/L11AB-7 (510 plates)

Lot Number: Z0010563

Expiration Date: 31 AUG 2014

Statement: This product was manufactured and tested according to approved MSD Standard Operating Procedures. This lot meets MSD's specifications.

This product has an approximate binding capacity of 0.9 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, tested using tagged BTI lot R0010381:

Concentration of SULFO-TAG™ Labeled Biotinylated IgG (BTI)	Mean ECL spec and tolerance range	Intraplate CV or SD spec	Pass/Fail?
18 nM (0.9 pmole)	76,752 counts \pm 15%	10%	PASS
12 nM (0.6 pmole)	62,123 counts \pm 15%	10%	PASS
6 nM (0.3 pmole)	35,647 counts \pm 15%	10%	PASS
0 nM (0.0 pmole)	Not to exceed 103 counts	15 count SD	PASS

Uniformity plates:

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; \leq 15% for all plates	PASS
Median signal for rows, max to min range	\leq 10% for at least 92% of plates; \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
Preparer	Sarah Robles	QC Technical Specialist		16 APR 2013
Review/Approval	Michael Eager	QC Scientist		08 MAY 2013