



CERTIFICATE OF ANALYSIS

COA - 03541

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

Catalog #: 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: Z0010553

Expiration date: Jun, 2014

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

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:

Concentration of SULFO-TAG™ Labeled Biotinylated IgG (BTI)	Mean ECL spec and tolerance range
18 nM (0.9 pmole)	Pass
12 nM (0.6 pmole)	Pass
6 nM (0.3 pmole)	Pass
0 nM (0.0 pmole)	Pass

Uniformity plates:

Metric	Pass/Fail?
Mean Intraplate CV	Pass
Intraplate CVs	Pass
Number of plates with intraplate CV >12%	Pass
Interplate CV	Pass
Plates where signal > 20% from plate mean occurs in same well on multiple plates	Pass
Wells with signal > 50% from plate mean	Pass
Median signal for concentric rings, max to min range	Pass
Median signal for columns, max to min range	Pass
Median signal for rows, max to min range	Pass
Number of plates sampled (N)	Pass

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

	Name	Function	Signature	Date
Preparer	Vernisha Greene	QC Analyst		15 Mar 2013
Review/Approval	Mike Eager	QC Scientist		15 MAR 2013