

# RayBio<sup>®</sup> Label-Based (L-Series) Rat Antibody Array 90 (L-90)

## Patent Pending Technology User Manual (Revised Dec 9<sup>th</sup>, 2019)

For the simultaneous detection of the relative expression of 90 (L-90) rat proteins in serum, plasma, cell culture supernatants, cell/tissue lysates or other body fluids.

**L-Series Rat Antibody Array L-90**  
**Cat# AAR-BLG-1-4 (4 Sample Kit)**  
**Cat# AAR-BLG-1-8 (8 Sample Kit)**

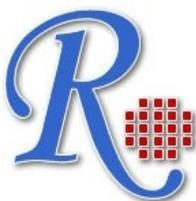
**Please read manual carefully  
before starting experiment**



**Your Provider for Excellent Protein Array Systems and Services**

**Tel: (Toll Free) 1-888-494-8555 or +1-770-729-2992; Fax: +1-770-206-2393;**

**Website: [www.raybiotech.com](http://www.raybiotech.com) Email: [info@raybiotech.com](mailto:info@raybiotech.com)**



## RayBiotech, Inc

---

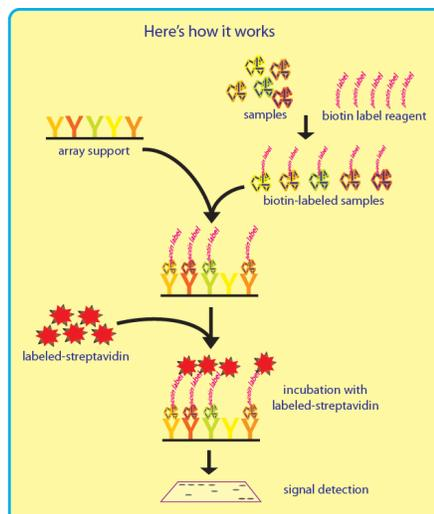
### **TABLE OF CONTENTS**

I.	Introduction and How It Works.....	2
II.	Materials Provided.....	3
	A. Storage Recommendations.....	3
	B. Additional Materials Required.....	4
III.	Overview and General Considerations.....	4
	A. Preparation and Storage of Samples.....	4
	B. Handling the Glass Slides.....	7
	C. Glass Slide Layout.....	8
	D. Incubation and Washes.....	8
IV.	Protocol.....	9
	A. Dialysis of Sample.....	9
	B. Biotin Labeling of Sample .....	9
	C. Drying of the Glass Chip.....	12
	D. Blocking and Incubations.....	13
	E. Fluorescence Detection.....	16
V.	Antibody Array Map.....	18
VI.	Interpretation of Results.....	20
VII.	Troubleshooting Guide.....	24
VIII.	Selected References.....	25

## I. Introduction

Recent technological advances by RayBiotech have enabled the largest commercially available antibody array to date. With the L-Series Antibody Array 90, researchers can now obtain a broad, panoramic view of cytokine expression. The expression levels of 90 rat target proteins can be simultaneously detected, including cytokines, chemokines, adipokine, growth factors, angiogenic factors, proteases, soluble receptors, soluble adhesion molecules and other proteins in cell culture supernatants, serum and plasma.

The first step in using the RayBio® L-Series Rat Antibody Array 90 is to biotinylate the primary amine of the proteins in serum or plasma samples, cell culture supernatant, cell lysate or tissue lysate. The glass slide arrays are then blocked, just like a Western blot, and the biotin-labeled sample is added onto the glass slide, which is pre-printed with capture antibodies, and incubated to allow for interaction of target proteins. Streptavidin-conjugated fluorescent dye (Cy3 equivalent) is then applied to the array. Finally, the glass slide is dried, and laser fluorescence scanning is used to visualize the signals.



## II. Materials Provided

### A. Storage Recommendations

Upon receipt, the kit should be stored at -20°C until needed. Please use within 6 months from the date of shipment. After initial use, remaining reagents should be stored at 4°C to avoid repeated freeze-thaw cycles (may be stored for up to 3 months, Labeling Reagent, Item B should be fresh preparation before use). Unused glass slides should be kept at -20 °C and avoid repeated freeze-thaw cycles (may be stored for up to 6 months).

### RayBio® L-Series Rat Antibody Array 90

ITEM	DESCRIPTION	Cat#: AAR-BLG-1-4	Cat#: AAR-BLG-1-8
A	Dialysis Vials	8 vials	16 vials
B	Labeling Reagent	2 vials	4 vials
D	Stop Solution	1 vial (50 ul)	
E	RayBio® L-series Rat Antibody Array L-90 Glass Slides*	1 L-90 Slide	2 L-90 Slides
F	Blocking Buffer	1 bottle (8 ml)	1 bottle (8 ml)
G	20X Wash Buffer I	1 bottle (30 ml)	1 bottle (30 ml)
H	20X Wash Buffer II	1 bottle (30 ml)	1 bottle (30 ml)
I	Cy3-Conjugated Streptavidin	1 vials	2 vials
J	Adhesive Plastic Strips		
K	Labeling Buffer	1 bottle (8 ml)	
L	Floating Dialysis Rack	1 rack	
N/A	2X Cell Lysis Buffer**	1 bottle (10 ml)	
M	30 ml Centrifuge Tube	1 tube	

\*Each slide contains 4 identical subarrays

\*\*Only needed if testing cell or tissue lysates

## **B. Additional Materials Required**

- Distilled or de-ionized water
- KCl, NaCl,  $\text{KH}_2\text{PO}_4$  and  $\text{Na}_2\text{HPO}_4$
- Small plastic or glass containers
- Orbital shaker or oscillating rocker
- Beaker, stir plate and stir bar
- 1 ml tube
- Pipettors, pipette tips and other common lab consumables
- Laser scanner for fluorescence detection (list available online)
- Aluminum foil

## ***III. Overview and General Considerations***

### **A. Preparation and Storage of Samples**

#### 1) Preparation of Cell Culture Supernatants

- Seed cells at a density of  $1 \times 10^6$  cells in 100 mm tissue culture dishes (\*).
- Culture in complete culture medium for ~24–48 hours (\*\*).
- Replenish with serum-free or low-serum medium such as 0.2% FCS/FBS serum, and then incubate cells again for ~48 hours (\*\*, †). Recommended using membrane-based array if using high serum medium such as 10% FCS/FBS, the glass slide arrays tend to have extremely high background for high serum containing media samples.
- To collect supernatants, centrifuge at 1,000 g for 10 min and store as  $\leq 1$  ml aliquots at  $-80^\circ\text{C}$  until needed.

- Measure the total wet weight of cultured cells in the pellet and/or culture dish. You may then normalize between arrays by dividing fluorescent signals by total cell mass (i.e., express results as the relative amount of protein expressed/mg total cell mass). Or you can normalize between array by determining cell lysate concentration using a total protein assay (BCA Protein Assay Kit, Pierce, Prod# 23227).

*Note: \* The density of cells per dish used is dependent on the cell type. More or less cells may be required.*

*\*\* Optimal culture time may vary and will depend on the cell line, treatment conditions and other factors.*

*† Bovine serum proteins produce detectable signals on the RayBio® L-Series Rat Antibody Array 90 in media containing serum concentrations as low as 0.2%. When testing serum-containing media, we strongly recommend testing an uncultured media blank for comparison with sample results.*

## **2) Extracting Protein from Cells**

- For attached cells, remove supernatant from cell culture, wash cells twice with cold 1X PBS.  
For suspension cells, pellet the cells by centrifuging using a microcentrifuge at 1500 rpm for 10 min.
- Make sure to remove any remaining PBS before adding 1X

Cell Lysis Buffer (2X Cell Lysis Buffer should be diluted 2 fold with deionized or distilled water). Solubilize the cells at  $2 \times 10^7$  cells/ml in 1X Cell Lysis Buffer.

- Pipette up and down to resuspend cells and rock the lysates gently at 2–8 °C for 30 minutes. Transfer extracts to microfuge tubes and centrifuge at 13,000 rpm for 10 min at 2-8 °C \*.
- Transfer supernatant to a clean tube. Determining cell lysate concentration using a total protein assay (BCA Protein Assay Kit, Pierce, Prod# 23227). Aliquot the lysates and store at – 70°C.

Note \*: If the supernatant appears to be cloudy, transfer the supernatants to a clean tube, centrifuge again at 13,000 rpm for 20 minutes at 2-8°C. If the supernatant is still not clear, store the lysate at -70°C for 20 minutes. Remove from the freezer, immediately centrifuge at 13,000 rpm for 20 minutes at 2-8°C.

### **3) *Extracting Protein from Crude Tissue***

- Transfer approximate 100 mg crude tissue into a tube with 1 ml 1X Cell Lysis Buffer (2X Cell Lysis Buffer should be diluted 2 fold with deionized or distilled water).
- Homogenize the tissue according to homogenizer manufacturer instructions.
- Transfer extracts to microcentrifuge tubes and centrifuge for 20 min at 13,000 rpm (4°C).
- Transfer supernatant to a clean tube and store at – 70°C.

Note \*: If the supernatant appears to be cloudy, transfer the

supernatants to a clean tube, centrifuge again at 13,000 rpm for 20 minutes at 2-8°C. If the supernatant is still not clear, store the lysate at -70°C for 20 minutes. Remove from the freezer, immediately centrifuge at 13,000 rpm for 20 minutes at 2-8°C.

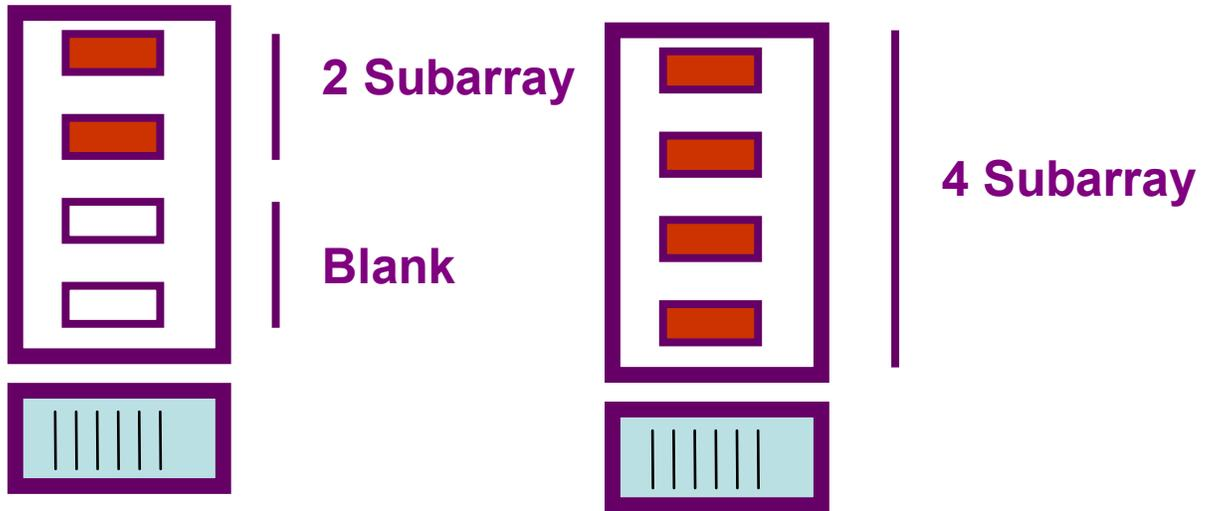
#### B. Handling the glass slides

- The microarray slides are delicate. Please do not touch the array surface with pipette tips, forceps or your fingers. Hold the slides by the edges only.
- Handle the slides with powder-free gloves and in a clean environment.
- Do not remove the glass slide from the chamber assembly until step 19, and take great care not to break the glass slide when doing so.
- Remove reagents/sample by gently applying suction with a pipette to corners of each chamber. Do not touch the printed area of the array, only the sides.



### C. Layout of Rat L-90 Glass Slide

Two or four identical sub-arrays on one slide



### D. Incubations and Washes

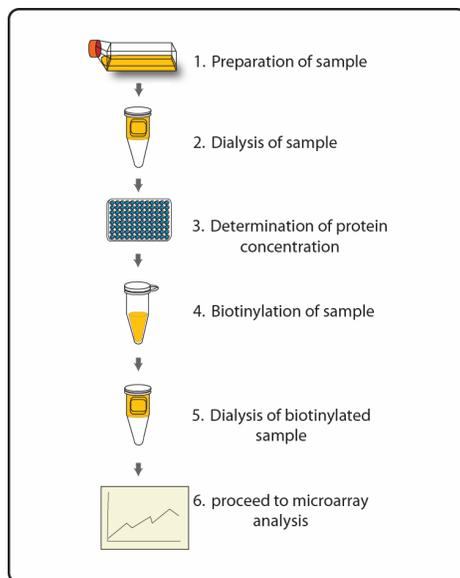
- Cover incubation chamber with a Plastic Adhesive Strip (Item J) to prevent evaporation during incubation or wash steps, particularly those lasting 2 hours or longer.
- During incubation and wash steps avoid foaming and be sure to remove all bubbles from the sub-array surface.
- Perform all incubation and wash steps under gentle rotation or rocking motion (~0.5 to 1 cycle/sec).
- Wash steps in Wash Buffer II and all incubation steps may be performed overnight at 4°C.
- Avoid cross-contamination of samples to neighboring wells. To remove Wash Buffers and other reagents from chamber wells, you may invert the Glass Slide Assembly to decant, and aspirate the remaining liquid.
- Unlike most Cy3 fluors, the streptavidin-conjugated fluor used in this kit is very stable at RT and resistant to

photobleaching on the hybridized glass slides. However, please protect glass slides from directly strong light and temperatures above RT.

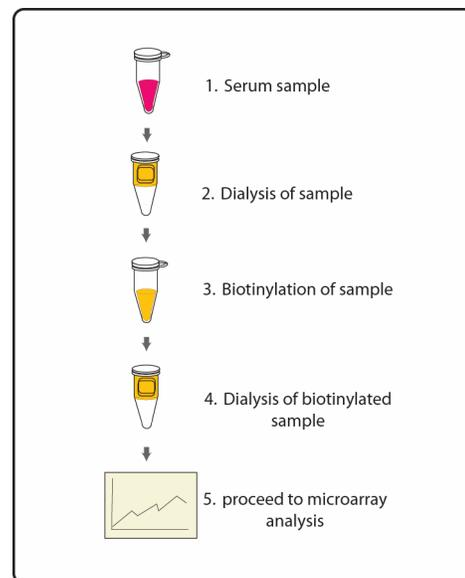
## IV. Protocol

### Assay Diagram

#### 1. Cell culture supernatants or cell/tissue lysates\*.



#### 2. Serum or plasma



\* If using cell or tissue lysates start at step 2. “Dialysis of sample”

#### A. Dialysis of Sample

*Note: Samples must be dialyzed prior to biotin-labeling (Steps 5–7).*

1. To prepare dialysis buffer (1X PBS, pH=8.0), dissolve 0.6 g KCl, 24 g NaCl, 0.6 g KH<sub>2</sub>PO<sub>4</sub> and 3.45 g Na<sub>2</sub>HPO<sub>4</sub> in 2500 ml de-ionized or distilled water. Adjust pH=8.0 with 1M NaOH and adjust final volume to 3000 ml with de-ionized or distilled water.
2. Add each sample into a separate Dialysis Tube (Item A). Load 200 µl cell culture supernatant\* or 100 µl cell lysates or tissue lysate (1~2 mg/ml total protein) or 20 µl serum or plasma + 80 µl dialysis buffer (5-fold dilution. Carefully place Dialysis Tubes into Floating Dialysis Rack (Item L).
3. Place Floating Dialysis Rack into ≥500 ml dialysis buffer in a large beaker. Place beaker on a stir plate and dialyze, for at least 3 hours at 4°C, stirring buffer gently. Then exchange the dialysis buffer and repeat dialysis for at least 3 h at 4°C. Transfer dialyzed sample to a clean eppendorf tube. Spin dialyzed samples for 5 min at 10,000 rpm to remove any particulates or precipitants, and then transfer the supernatants to a clean tube.

*Note: The sample volume may change during dialysis.*

*Note: Dialysis procedure may proceed overnight.*

*Note: Determine the total protein concentration for cell culture supernatants or cell/tissue lysate after dialysis procedure (Step 3). We recommended using a BCA total protein assay (eg, Pierce, Catalog # 23227).*

## B. Biotin-labeling Sample

*Note: Amines (e.g., Tris, glycine) and azides quench the biotinylation reaction. Avoid contaminating samples with these chemicals prior to biotinylation.*

4. Immediately before use, prepare 1X Labeling Reagent. Briefly spin down the Labeling Reagent tube (Item B). Add 100  $\mu$ l 1X PBS into the tube, pipette up and down or vortex slightly to dissolve the lyophilized reagent.
5. Add 1X Labeling Reagent to dialyzed samples.
  - a) **For labeling cell culture supernatants:** transfer 180  $\mu$ l dialyzed sample into a new tube. Add 36  $\mu$ l of 1X Labeling Reagent Solution per 1 mg total protein in dialyzed cell culture supernatant. Mix well. For example, if sample's total protein concentration is 0.5 mg/ml you need to add 3.24  $\mu$ l 1X Labeling Reagent to 180  $\mu$ l dialyzed sample.
  - b) **For labeling serum or plasma:** Add 22  $\mu$ l of 1X Labeling Reagent Solution into a new tube containing 35  $\mu$ l\* dialyzed serum or plasma sample and 155  $\mu$ l Labeling Buffer (Item K).

*Note \*:* To normalize serum/plasma concentrations during biotinylation, measure sample volume before and after dialysis. Then adjust the volumes of dialyze serum/plasma and Labeling Buffer to compensate (to keep same total protein amount and total volume). For example, if serum

*/plasma sample volume increased from 100  $\mu$ l to 200  $\mu$ l, add 70  $\mu$ l dialyzed serum and 120  $\mu$ l Labeling Buffer to keep same total volume, 212  $\mu$ l.*

- c) **For labeling cell or tissue lysates:** transfer 30  $\mu$ g (15  $\mu$ l of 2 mg/ml) cell or tissue lysates into a tube and add labeling buffer (Item K) for a total volume of 300  $\mu$ l. Then add 3.3  $\mu$ l of 1X Labeling Reagent Solution.
6. Incubate the reaction solution at room temperature with gentle rocking or shaking for 30 min. Mix the reaction solution by gently tapping the tube every 5 min.
7. 3  $\mu$ l Stop Solution (Item D) into each reaction tube. Collect and transfer each sample from reaction tube into a separate Dialysis Tube (Item A). Immediately dialyze samples as directed in Step 3 on pages 9.

*Note: Biotinylated samples can be stored at -20°C or -80°C until you are ready to proceed with the assay.*

### **C. Drying of the Glass Slide**

8. Remove the package containing the Assembled Glass Slide (Item E) from the freezer. Place unopened package on the bench top for approx. 15 min, and allow the Assembled Glass Slide to equilibrate to room temperature (RT).
9. Open package, and take the Assembled Glass Slide out of the sleeve (Do not disassemble the Glass Slide from the chamber assembly). Place glass slide assembly in laminar flow hood

or similar clean environment for 1-2 hours at RT.

*Note: Protect the slide from dust or others contaminants.*

#### **D. Blocking and Incubations**

***Note: Glass slide should be completely dry before adding Blocking Buffer to wells.***

10. Block sub-arrays by adding 400 µl of Blocking Buffer (Item F) into each well of Assembled Glass Slide and incubating at RT for 30 min. Ensure there are no bubbles on the array surfaces.

11. Immediately prior to sample incubation, spin biotin-labeled samples for 5 min at 10,000 rpm to remove any particulates or precipitants. Dilute samples with Blocking Buffer.\*

*\*Note: Recommended dilution of the biotin-labeled samples with Blocking Buffer prior to incubation is 2-10 fold for cell culture supernatants, 20-fold for serum/plasma or 30 fold cell/tissue lysate .*

*Note: Optimal sample dilution factor will depend on the abundance of target proteins. If the background or antigen-specific antibody signals are too strong, the sample can be diluted further in subsequent experiments. If the signal is too weak, more concentrated samples can be used.*

12. Completely remove Blocking Buffer from each well. Add 400  $\mu$ l of diluted samples into appropriate wells. Remove any bubbles on array surfaces. Incubate arrays with gentle rocking or shaking for 2 hours at RT or overnight at 4°C.

*Note: Avoid the flow of sample into neighboring wells.*

13. Dilute 20X Wash Buffer I Concentrate (Item G) 20-fold with de-ionized or distilled water. Decant the samples from each well, and wash 3 times with 800  $\mu$ l of 1X Wash Buffer I at RT with gentle rocking or shaking for 5 min per wash.

14. Obtain a clean container (e.g., pipette tip box or slide-staining jar), place the Assembled Glass Slide into the container with enough volume of 1X Wash Buffer I to completely cover the entire assembly, and remove any bubbles in wells. Wash 2 times at RT with gentle rocking or shaking for 10 min per wash.

15. Dilute 20X Wash Buffer II Concentrate (Item H) 20-fold with de-ionized or distilled water. Decant the Wash Buffer I from each well, place the Assembled Glass Slide into the container with enough volume of 1X Wash Buffer II to completely cover the entire assembly, and remove any bubbles in wells. Wash 2 times at RT with gentle rocking or shaking for 5 min per wash.

16. Prepare 1X Cy3-Conjugated Streptavidin:

- a) Briefly spin down tube containing the Cy3-Conjugated Streptavidin (Item I) immediately before use.

- b) Add 1000  $\mu$ l of Blocking Buffer into the tube to prepare a concentrated Cy3-Conjugated Streptavidin stock solution. Pipette up and down to mix gently (do not store the stock solution for later use).
- c) Add 200  $\mu$ l of Cy3-Conjugated Streptavidin stock solution into a tube with 800  $\mu$ l of Blocking Buffer. Mix gently to prepare 1X Cy3-Conjugated Streptavidin.

17. Carefully remove Assembled Glass Slide from container. Remove all of Wash Buffer II from the wells. Add 400  $\mu$ l of 1X Cy3-Conjugated Streptavidin to each sub-array. Cover the incubation chamber with the plastic adhesive strips.

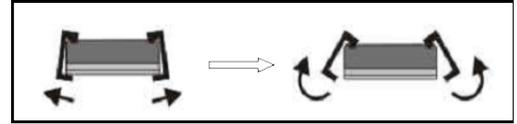
*Note: Avoid exposure to light in Steps 19–25 by covering the Glass Slide Assembly with aluminum foil or incubate in dark room.*

18. Incubate with Cy3-Conjugated Streptavidin at RT for 2 hours with gentle rocking or shaking.

*Note: Incubation may be done overnight at 4°C.*

19. Decant the solution and disassemble the glass slide from the incubation frame and chamber. Disassemble the device by pushing clips outward from the side, as shown below. Carefully remove the glass slide from the gasket.

**Note: Be careful not to touch the printed surface of the glass slide, which is on the same side as the barcode.**



20. Gently place the glass slide into 30 ml Centrifuge Tube (Item M). Add enough 1X Wash Buffer I to cover the entire glass slide. Wash with gentle rocking or shaking for 10 min. Remove the wash buffer. Repeat 2 times for a total of 3 washes.
21. Repeat step 20, this time with 1X Wash Buffer II. Repeat one time for a total of two washes for 5 min per wash.
22. Finally, wash the glass slide with 30 ml of de-ionized or distilled water for 5 min. Remove glass slide and decant water from Centrifuge Tube.
23. Remove excess liquid from Centrifuge Tube, and place glass slide into the tube. Centrifuge at 1,000 rpm for 3 minutes to remove water droplets. Make sure the finished glass slide is completely dry before scanning or storage.

**Note: Alternatively, you may gently dry the glass slide using a low-velocity Nitrogen gas stream or ambiently in a laminar flow hood or similar clean environment (Be sure to protect from light).**

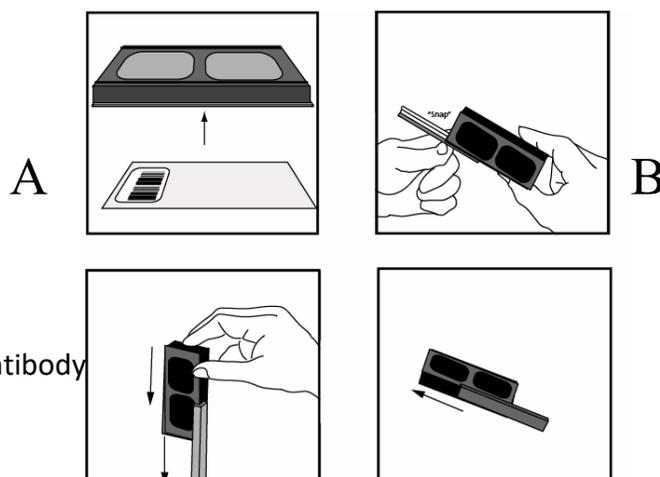
## **E. Fluorescence Detection**

24. You may proceed immediately to scanning or you may store the slide at -20 °C in the Centrifuge Tube provided or at RT and to scan at a later time.

**Note:** *Unlike most Cy3 fluors, the Streptavidin-Conjugated Fluorused in this kit is very stable at RT and resistant to photobleaching on completed glass slides. However, please protect glass slides from temperatures above RT and store them in the dark. Do not expose glass slide to strong light, such as sunlight or UV lamp.*

**Note:** *If you need to repeat any of the incubation after finishing the experiment, you must first re-assemble the glass slide into the incubation chamber by following step as shown in the figures below. To avoid breaking the printed glass slide, you may first want to practice assembling the device with a blank glass slide.*

1. Apply slide to incubation chamber barcode facing upward as in image A (below).
2. Gently snap one edge of a snap-on side as shown in image B.
3. Gently press other of side against lab bench and push in lengthwise direction (image C).
4. Repeat with the other side (image D)



# V. Antibody Array Map

C

D

## RayBio® Biotin Label-based Rat Antibody Array 1 Map

RayBio® L-Series Rat Antibod

	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15
1	Positive 1	Positive 1	Positive 2	Positive 2	Positive 3	Positive 3	Neg	Neg	Activin A	Activin A	ACTH	ACTH	ADFP	ADFP	Adiponectin/Acp30
2	CD106	CD106	CINC-2 alpha/beta	CINC-2 alpha/beta	CINC-3	CINC-3	CNTF	CNTF	CNTF R alpha	CNTF R alpha	CSK	CSK	CXCR4	CXCR4	EGFR
3	Fractalkine	Fractalkine	GFR alpha-1	GFR alpha-1	GFR alpha-2	GFR alpha-2	GM-CSF	GM-CSF	Growth Hormone	Growth Hormone	Growth Hormone R	Growth Hormone R	Hepassocin	Hepassocin	ICAM-1/CD54
4	IL-3	IL-3	IL-4	IL-4	IL-5	IL-5	IL-6	IL-6	IL-10	IL-10	IL-12/IL-23 p40	IL-12/IL-23 p40	IL-13	IL-13	Integrin alpha M beta 2
5	MIF	MIF	MIP-1 alpha	MIP-1 alpha	MIP-2	MIP-2	MIP-3 alpha	MIP-3 alpha	MIP-2	MIP-2	MMP-3	MMP-3	MMP-13	MMP-13	MusK
6	RA1/TMIG-6	RA1/TMIG-6	RELM beta	RELM beta	Resistin	Resistin	TLIA	TLIA	TGF-beta1	TGF-beta1	TGF-beta2	TGF-beta2	TGF-beta3	TGF-beta3	Thrombospondin
7	TROY	TROY	Ubiquitin	Ubiquitin	VEGF	VEGF	VEGF-C	VEGF-C	Neg	Neg	Neg	Neg	Neg	Neg	Neg

	16	17	18	19	20	21	22	23	24	25	26	27	28	29	30
Adiponectin/Acp30	AMPK alpha 1	AMPK alpha 1	AMPK alpha 1	B7-1/CD80	B7-1/CD80	BDNF	BDNF	beta-Casitin	beta-Casitin	beta-FCF	beta-FCF	beta-NGF	beta-NGF	CCR4	CCR4
EGFR	EG-VEGFPK1	EG-VEGFPK1	EG-VEGFPK1	E-Selectin	E-Selectin	FADD	FADD	Fas/TNFRSF6	Fas/TNFRSF6	Fas Ligand/TNRSF6	Fas Ligand/TNRSF6	FGF-BP	FGF-BP	Follistatin-like-1(FSL1)	Follistatin-like-1(FSL1)
ICAM-1/CD54	ICK	ICK	ICK	IDE (Insulin Degrading Enzyme)	IDE (Insulin Degrading Enzyme)	IFN-gamma	IFN-gamma	IL-1 alpha	IL-1 alpha	IL-1 beta	IL-1 beta	IL-1 R/IL-1 R p2	IL-1 R/IL-1 R p2	IL-2	IL-2
Integrin alpha M beta 2	Insulin	Insulin	Insulin	IP-10	IP-10	Leprotin (OB)	Leprotin (OB)	LIX	LIX	L-Selectin/CD62L	L-Selectin/CD62L	MCP-1	MCP-1	MDC	MDC
MusK	Neuropilin-2	Neuropilin-2	Neuropilin-2	NGR	NGR	Oxetin A	Oxetin A	Osteopontin/SPP1	Osteopontin/SPP1	PDGF-AA	PDGF-AA	Prolectin R	Prolectin R	RAGE	RAGE
Thrombospondin	THE-2	THE-2	THE-2	TIMP-1	TIMP-1	TIMP-2	TIMP-2	TIMP-3	TIMP-3	TLR4	TLR4	TNF-alpha	TNF-alpha	TRAIL	TRAIL
Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Positive 3	Positive 3	Positive 2	Positive 2	Positive 1	Positive 1

## RayBio® Biotin Label-based Rat Antibody Array 1 List

	Target protein		Target protein		Target protein		Target protein
1	Positive 1	28	Fas Ligand/TNFSF6	55	IP-10	82	TGF-beta3
2	Positive 2	29	FGF-BP	56	Leptin (OB)	83	Thrombospondin
3	Positive 3	30	Follistatin-like -1(FSL1)	57	LIX	84	TIE-2
4	Neg	31	Fractalkine	58	L-Selectin/CD62L	85	TIMP-1
5	Activin A	32	GFR alpha-1	59	MCP-1	86	TIMP-2
6	ACTH	33	GFR alpha-2	60	MDC	87	TIMP-3
7	ADFP	34	GM-CSF	61	MIF	88	TLR4
8	Adiponectin/Acrp30	35	Growth Hormone	62	MIP-1 alpha	89	TNF-alpha
9	AMPK alpha 1	36	Growth Hormone R	63	MIP-2	90	TRAIL
10	B7-1/CD80	37	Hepassocin	64	MIP-3 alpha	91	TROY
11	BDNF	38	ICAM-1/CD54	65	MMP-2	92	Ubiquitin
12	beta-Catenin	39	ICK	66	MMP-8	93	VEGF
13	basic-FGF	40	IDE (Insulin Degrading Enzyme)	67	MMP-13	94	VEGF-C
14	beta-NGF	41	IFN-gamma	68	MuSK	95	Neg
15	CCR4	42	IL-1 alpha	69	Neuropilin-2	96	Neg
16	CD106	43	IL-1 beta	70	NGFR	97	Neg
17	CINC-2 alpha/beta	44	IL-1 R6/IL-1 R rp2	71	Orexin A	98	Neg
18	CINC-3	45	IL-2	72	Osteopontin/SPP1	99	Neg
19	CNTF	46	IL-3	73	PDGF-AA	100	Neg
20	CNTF R alpha	47	IL-4	74	Prolactin R	101	Neg
21	CSK	48	IL-5	75	RAGE	102	Neg
22	CXCR4	49	IL-6	76	RALT/MIG-6	103	Positive 3
23	EGFR	50	IL-10	77	RELM beta	104	Positive 2
24	EG-VEGF/PK1	51	IL-12/IL-23 p40	78	Resistin	105	Positive 1
25	E-Selectin	52	IL-13	79	TL1A		
26	FADD	53	Integrin alpha M beta 2	80	TGF-beta1		
27	Fas/TNFRSF6	54	Insulin	81	TGF-beta2		

## VI. Interpretation of Results:

### A. Explanation of Controls Spots

- 1) **Positive Control spots (POS1, POS2, POS3)** are standardized amounts of biotinylated IgGs printed directly onto the array. All other variables being equal, the Positive Control intensities will be the same for each sub-array. This allows for normalization based upon the relative fluorescence signal responses to a known control, much as “housekeeping” genes or proteins are used to normalize results in PCR or Western blots, respectively.
- 2) **Negative Control (NEG) spots** contain a protein-containing buffer (used to dilute antibodies printed on the array). Their signal intensities represent non-specific binding of Biotin-conjugated anti-Cytokines and/or the Cy3-Conjugated Streptavidin. Negative control signal intensities are usually very close to background signals in each sub-array.

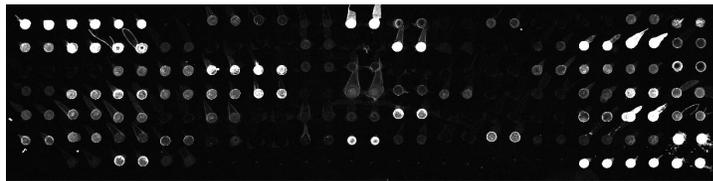
### B. Typical results obtained with RayBio® L-Series Rat Obesity Antibody Array L-90

The following figure shows the RayBio® L-Series Rat Antibody Array 90 probed with serum sample. The images were captured using a Axon GenePix laser scanner. The strong signals

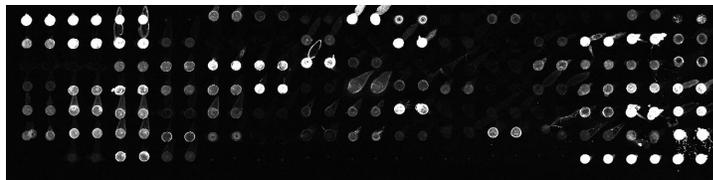
in row 20 and the upper left and lower right corners of each array are Positive Controls, which can be used to identify the orientation and help normalize the results between arrays.

## RayBio® L-Series Rat Antibody Array 90

### Sample-1



### Sample-2



If scanned using optimal settings, 3 distinct signal intensities will be seen: POS1>POS2>POS3. If all of these signals are of similar intensity, try increasing or decreasing laser power and/or signal gain settings.

Also, in the absence of an external standard curve for each protein detected, there is no means of assessing absolute or relative concentrations of different proteins in the same sample using immunoassays. If you wish to obtain quantitative data (ie,

concentrations of the various analytes in your samples), try using our Quantibody® Arrays instead.

### **C. Background Subtraction:**

Once you have obtained fluorescence intensity data, you should subtract the background and normalize to the Positive Control signals before proceeding to analysis.

Most laser fluorescence scanner software have an option to automatically measure the local background around each spot. For best results, we recommend comparing signal intensities representing the MEDIAN background signals minus local background. If your resulting fluorescence signal intensity reports do not include these values (e.g., a column labeled as “MED532-B532”), you may need to subtract the background manually or change the default settings on your scanner’s data report menu.

### **D. Normalization of Array Data:**

To normalize signal intensity data, one sub-array is defined as "reference" to which the other arrays are normalized. This choice is arbitrary. For example, in our Analysis Tool Software (described below), the array represented by data entered in the left-most column each worksheet is the default “reference array.”

**You can calculate the normalized values as follows:**

$$X(Ny) = X(y) * P1/P(y)$$

**Where:**

**P1 = mean signal intensity of POS spots on reference array**

**P(y) = mean signal intensity of POS spots on Array "y"**

**X(y) = mean signal intensity for spot "X" on Array "y"**

**X(Ny) = normalized signal intensity for spot "X" on Array "y"**

The RayBio® Analysis Tool software is available for use with data obtained using RayBio® Biotin Label-based Antibody Arrays. You can copy and paste your signal intensity data (with and without background) into the Analysis Tool, and it will automatically normalize signal intensities to the Positive Controls.

To order the Analysis Tool, please contact us at +1-770-729-2992 or [info@raybiotech.com](mailto:info@raybiotech.com) for more information.

#### **E. Threshold of significant difference in expression:**

After subtracting background signals and normalization to Positive Controls, comparison of signal intensities between and among array images can be used to determine relative differences in expression levels of each protein between samples or groups.

Any  $\geq 1.5$ -fold increase or  $\leq 0.65$ -fold decrease in signal intensity for a single analyte between samples or groups may be considered a measurable and significant difference in expression, provided that both sets of signals are well above background (Mean background + 2 standard deviations, accuracy  $\approx 95\%$ ).

## VII. Troubleshooting Guide

<b>Problem</b>	<b>Cause</b>	<b>Recommendation</b>
Weak signal	Inadequate detection	Check laser power and PMT parameters
	Inadequate reagent volumes or improper dilution	Check pipettors and ensure correct preparation
	Short incubation times	Ensure sufficient incubation time and change sample incubation step to overnight
	Too low protein concentration in sample	Don't make too low dilution Or concentrate sample
	Improper storage of kit	Store kit at suggested temperature
High background	Sample is too concentrated	Use more diluted sample
	Excess of streptavidin	Make sure to use the correct amount of streptavidin
	Inadequate detection	Check laser power and PMT parameters
	Inadequate wash	Increase the volume of wash buffer and incubation time
Uneven signal	Bubbles formed during incubation	Avoid bubble formation during incubation
	Arrays are not completely covered by reagent	Completely cover arrays with solution

## VIII. Selected References

1. Christina Scheel et al. Paracrine and Autocrine Signals Induce and Maintain Mesenchymal and Stem Cell States in the Breast. *Cell*. 2011;145, 926–940
2. Lin Y, Huang R, Chen L, et al. Profiling of cytokine expression by biotin-labeled-based protein arrays. *Proteomics*. 2003, 3: 1750–1757.
3. Huang R, Jiang W, Yang J, et al. A Biotin Label-based Antibody Array for High-content Profiling of Protein Expression. *Cancer Genomics Proteomics*. 2010; 7(3):129–141.
4. Liu T, Xue R, Dong L, et al. Rapid determination of serological cytokine biomarkers for hepatitis B–virus-related hepatocellulare carcinoma using antibody arrays. *Acta Biochim Biophys Sin*. 2011; 43(1):45–51.
5. Cui J, Chen Y, Chou W-C, et al. An integrated transcriptomic and computational analysis for biomarker identification in gastric cancer. *Nucl Acids Res*. 2011; 39(4):1197–1207.
6. Jun Zhong et al. Temporal Profiling of the Secretome during Adipogenesis in Humans. *Journal of Proteome Research*. 2010, 9, 5228–5238

7. Chowdury UR, Madden BJ, Charlesworth MC, Fautsch MP. Proteomic Analysis of Human Aqueous Humor. *Invest Ophthalmol Visual Sci.* 2010; 51(10):4921–4931.
8. Wei Y, Cui C, Lainscak M, et al. Type-specific dysregulation of matrix metalloproteinases and their tissue inhibitors in end-stage heart failure patients: relationship between MMP-10 and LV remodeling. *J Cell Mol Med.* 2011; 15(4):773–782.
9. Kuranda K, Berthon C, Lepêtre F, et al. Expression of CD34 in hematopoietic cancer cell lines reflects tightly regulated stem/progenitor-like state. *J Cell Biochem.* 2011; 112(5):1277–1285.
10. Toh HC, Wang W-W, Chia WK, et al. Clinical Benefit of Allogenic Melanoma Cell Lysate-Pulsed Autologous Dendritic Cell Vaccine in MAGE-Positive Colorectal Cancer Patients. *Clin Chem Res.* 2009; 15:7726–7736.
11. Zhen Hou. Cytokine array analysis of peritoneal fluid between women with endometriosis of different stages and those without endometriosis Biomarkers. 2009;14(8): 604-618.
12. Yao Liang Tang, et al. Hypoxic Preconditioning Enhances the Benefit of Cardiac Progenitor Cell Therapy for Treatment of Myocardial Infarction by Inducing CXCR4. *Circ Res.* 2009;109:197723

RayBio<sup>®</sup> L-series Antibody Arrays are patent-pending technology developed by RayBiotech.

This product is intended for research only and is not to be used for clinical diagnosis. Our products may not be resold, modified for resale, or used to manufacture commercial products without written approval by RayBiotech, Inc.

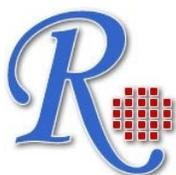
Under no circumstances shall RayBiotech be liable for any damages arising out of the use of the materials.

Products are guaranteed for six months from the date of shipment when handled and stored properly. In the event of any defect in quality or merchantability, RayBiotech's liability to buyer for any claim relating to products shall be limited to replacement or refund of the purchase price.

RayBio<sup>®</sup> is a registered trademark of RayBiotech, Inc.

GenePix<sup>®</sup> is a registered trademark of Molecular Devices, Inc.

**This product is for research use only.**



*©2011 RayBiotech, Inc.*