RayBio® C-Series Mouse RTK Phosphorylation Antibody Array C1

For the semi-quantitative detection of 71 phosphorylated Mouse receptor tyrosine kinases (RTKs) in cell and tissue lysates

Patent Pending Technology

User Manual (Revised September 23, 2021)

Cat# AAM-PRTK-1-2 (2 Sample Kit) Cat# AAM-PRTK-1-4 (4 Sample Kit) Cat# AAM-PRTK-1-8 (8 Sample Kit)

Please read manual carefully before starting experiment



Tel: (Toll Free) 1-888-494-8555 or +1-770-729-2992; Fax: +1-770-206-2393; Website: www.raybiotech.com Email: info@raybiotech.com



C-Series Antibody Arrays

TABLE OF CONTENTS

l.	Introduction	2
II.	How It Works	3
III.	Components and Storage	4
IV.	Additional Materials Required	4
V.	Sample Tips and General Considerations	5
	A. Sample Collection, Preparation, and Storage	5
	B. Sample Types and Recommended Dilutions/Amounts	5
	C. Handling Membranes	5
	D. Incubations and Washes	6
VI.	Chemiluminescence Detection Tips	6
VII.	Component Preparation	7
VIII.	Protocol	8
	A. Blocking	8
	B. Sample Incubation	8
	C. First Wash	8
	D. Biotinylated Antibody Cocktail Incubation	8
	E. Second Wash	9
	F. HRP-Streptavidin Incubation	9
	G. Third Wash	9
	H. Chemiluminescent Detection	9
	I. Storage	10
IX.	Typical Results	10
Χ.	Interpreting the Results	11
	A. Control Spots	11
	B. Data Extraction	11
	C. Data Analysis	11
XI.	Array Map	13
XII.	Troubleshooting Guide	14

I. INTRODUCTION

Protein phosphorylation plays an unusually prominent role in cell signaling, development and growth. The RayBio[®] Mouse RTK Phosphorylation Antibody Array C1 is a very rapid, convenient and sensitive assay to simultaneous detect multiple protein phosphorylations and can be used to monitor activation or function of important biological pathways.

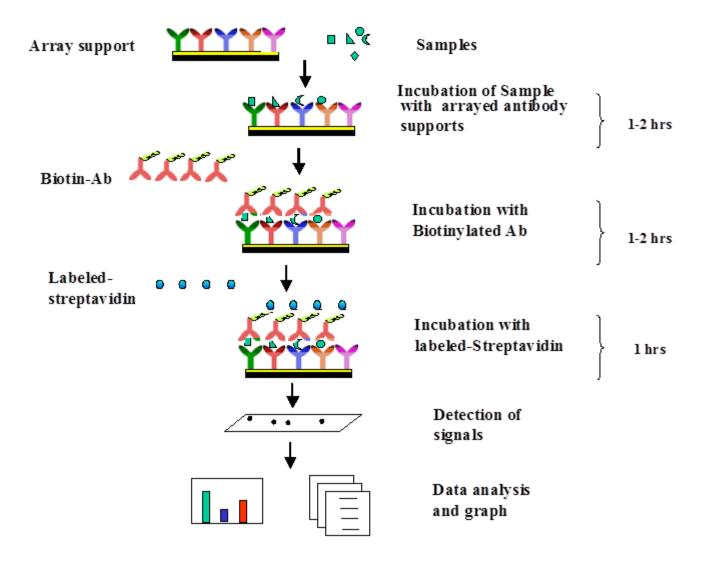
RayBiotech is committed to developing a series of phosphorylation antibody arrays. Our first product in this series is RayBio[®] Mouse RTK Phosphorylation Antibody Array C1, which is specifically designed for simultaneously identifying the relative levels of phosphorylation of 71 different Mouse receptor tyrosine kinases (RTKs) in cell lysates. By monitoring the changes in protein tyrosine phosphorylation in your experimental model system, you can verify pathway activation in your cell lines without spending excess time and effort in performing an analysis of immunoprecipitation and/or Western Blot.

By using RayBio[®] Mouse RTK Phosphorylation Antibody Array C1, treated or untreated cell lysate is added into antibody array membranes. The antibody array membranes are washed and biotinylated antiphosphotyrosine antibody is used to detect phosphorylated tyrosines on activated receptors. After incubation with HRP-streptavidin, the signals are visualized by chemiluminescence.

RayBio® C-Series Antibody Arrays have several advantages over detection of cytokines using single-target ELISA kits:

- 1. <u>More Data, Same or Less Sample</u>: Antibody arrays provide high-content screening using about the same sample volume as traditional ELISA.
- 2. <u>Global View of Cytokine Expression</u>: Antibody array screening improves the chances for discovering key factors, disease mechanisms, or biomarkers related to cytokine signaling.
- 3. <u>Similar (sometimes better) Sensitivity</u>: As little as 4 pg/ml of MCP-1 can be detected using the C-Series array format. In contrast, our similar MCP-1 ELISA assay has a sensitivity of 40 pg/ml of MCP-1.
- 4. <u>Increased Range of Detection</u>: ELISA assays typically detect a concentration range of 100- to 1000-fold, however, RayBiotech arrays can detect IL-2 at concentrations of 25 to 250,000 pg/ml, a range of 10,000-fold.
- 5. <u>Better Precision</u>: As determined by densitometry, the inter-array Coefficient of Variation (CV) of spot signal intensities is 5-10%, comparing favorably with ELISA testing (CV = 10-15%).

II. HOW IT WORKS



III. COMPONENTS AND STORAGE

Store kit at ≤ -20 °C immediately upon arrival. Kit must be used within the 6 months expiration

ПЕМ	COMPONENT	AAM-PRTK-1- 2	AAM-PRTK- 1-4	AAM-PRTK-1-	STORAGE TEMPERATURE AFTER THAWING**	
1	Antibody Arrays	2 membranes 4 membranes 8 membranes		≤-20°C		
2	Blocking Buffer	1 vial (25 ml)	2 vials (25 ml/ea)	≤-20 C	
3	Biotinylated Anti-Phosphotyrosine	1 vial	2 vials	4 vials	2-8°C	
3	Antibody	1 Viai 2 Viais		4 VIdiS	(for up to 3 days after dilution)	
4	1,000X HRP-Streptavidin	1 vial (50 μl)				
4	Concentrate					
5	20X Wash Buffer I Concentrate	1 vial (30 ml)		2 vial (30 ml/ea)	2 - 8 °C	
6	20X Wash Buffer II Concentrate	1 vial (30 ml)		1 vial (30 ml) 2 vial (30 ml/ea)		
7	2X Cell Lysis Buffer Concentrate	1 vial (10 ml)		1 vial (10 ml) 1 vial (16 ml)		
8	Detection Buffer C	1 vial (1.5 ml)		1 vial (1.5 ml) 1 vial (2.5 ml)		
9	Detection Buffer D	1 vial (1.5 ml)		1 vial (1.5 ml) 1 vial (2.5 ml)		
10	8-Well Incubation Tray w/ Lid	1 tray			Room Temperature	
11	Protease Inhibitor Cocktail	1 vial		1 vial 2 vials		
12	Phosphatase Inhibitor Cocktail II	1 vial		2 vials	≤-20°C	
Other	Other Kit Components: Plastic Sheets, Array Map Template, User Manual					

date.

IV. ADDITIONAL MATERIALS REQUIRED

- Pipettors, pipet tips and other common lab consumables
- Orbital shaker or oscillating rocker
- Tissue paper, blotting paper or chromatography paper
- Adhesive tape or plastic wrap
- Distilled or de-ionized water
- A chemiluminescent blot documentation system:
 - o CCD Camera
 - X-Ray Film and a suitable film processor
 - o Gel documentation system
 - Or another chemiluminescent detection system capable of imaging a western blot

^{*}Each package contains 2 or 4 membranes

^{**}For up to 3 months (unless stated otherwise) or until expiration date

V. SAMPLE TIPS AND GENERAL CONSIDERATIONS

A. Sample Collection, Preparation, and Storage

NOTE: Optimal methods will need to be determined by each experimenter empirically based on researched literature and knowledge of the samples.

- If not using fresh samples, freeze samples as soon as possible after collection.
- Avoid multiple freeze-thaw cycles. If possible, sub-aliquot samples prior to initial storage.
- It is strongly recommended to add a protease inhibitor cocktail to cell and tissue lysate samples.
- Avoid sonication of 1 ml or less as this can quickly heat and denature proteins
- Most samples will not need to be concentrated. If concentration is required, a spin column concentrator with a chilled centrifuge is recommended.
- Always centrifuge the samples hard after thawing (~10,000 RPM for 2-5 minutes) in order to remove any particulates that could interfere with detection.
- The Cell Lysate can be prepared as follows:
 - For attached cells, remove supernatant from cell culture, wash cells twice with cold 1X PBS (for suspension cells, pellet the cells by spinning down the cells at 1500 rpm for 10 min) making sure to remove any remaining PBS before adding Lysis Buffer. Solubilize the cells at 2x10⁷ cells/ml in 1X Lysis Buffer containing Protease Inhibitor Cocktail and Phosphatase Inhibitor Cocktail Set II (see preparation note shown on page 7 under Component Preparation Section). 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 14,000 x g for 10 min.

It is recommended that sample protein concentrations be determined using a total protein assay. For incubation with the Phosphorylation Antibody Array I, use at a protein concentration of 50-1000 μ g/ml for cell lysates.

Lysates should be used immediately or aliquot and stored at -70 °C. Thawed lysates should be kept on ice prior to use.

If you experience high background, you may further dilute your samples. If signals are too weak, the cell lysates can be pretreated by immunoprecipitations before incubation with array membranes. Immunoprecipitations can be done using anti-phosphotyrosine and protein A.

General tips for preparing lysate samples can be viewed on the online Resources page of the website.

B. Sample Types and Recommended Dilutions/Amounts

NOTE: Optimal sample dilutions and amounts will need to be determined by each experimenter empirically but the below recommendations may be used as a starting point. Blocking Buffer (ITEM 2) should be used to dilute samples. Normalize by loading equal amounts of protein per sample.

• Cell and Tissue Lysates: load 50 to 1000 μ g of total protein (after at least a 5-fold dilution to minimize the effect of any detergent(s)). Therefore the original lysate concentration should be 250 μ g to 5 mg/ml.

C. Handling Membranes

- The antibody printed side of each membrane is marked by a dash (-) or number (#) in the upper left corner.
- Do not allow membranes to dry out during the experiment or they may become fragile and break OR high and/or uneven background may occur.
- Grasp membranes by the corners or edges only using forceps. DO NOT touch printed antibody spots.

D. Incubations and Washes

- Perform <u>ALL</u> incubation and wash steps under gentle rotation or rocking motion (~0.5 to 1 cycle/sec) using an orbital shaker or oscillating rocker to ensure complete and even reagent/sample coverage. Rocking/rotating too vigorously may cause foaming or bubbles to appear on the membrane surface which should be avoided.
- All washes and incubations should be performed in the Incubation Tray (ITEM 10) provided in the kit.
- Cover the Incubation Tray with the lid provided during all incubation steps to avoid evaporation and outside debris contamination.
- Ensure the membranes are completely covered with sufficient sample or reagent volume during each incubation.
- Avoid forceful pipetting directly onto the membrane; instead, gently pipette samples and reagents into a corner of each well.
- Aspirate samples and reagents completely after each step by suctioning off excess liquid with a pipette. Tilting the tray so the liquid moves to a corner and then pipetting is an effective method.
- Optional overnight incubations may be performed for the following steps to increase overall spot signal intensities:
 - Sample Incubation
 - Biotinylated Antibody Cocktail Incubation
 - HRP-Streptavidin Incubation

NOTE: Overnight incubations should be performed at 4 °C (also with gentle rocking/shaking). Be aware that longer incubations can also increase the background response so complete liquid removal and washing is critical.

VI. CHEMILUMINESCENCE DETECTION TIPS

- Beginning with adding the detection buffers and ending with exposing the membranes should take no more than 10-15 minutes as the chemiluminescent signals may start to fade at this point.
- Trying multiple exposure times is recommended to obtain optimum results.
- A few seconds to a few minutes is the recommended exposure time range, with 30 seconds to 1 minute being suitable for most samples.

VII. COMPONENT PREPARATION

NOTE: Thaw all reagents to room temperature immediately before use. If wash buffers contain visible crystals, warm to room temperature and mix gently until dissolved.

NOTE: The Biotinylated Antibody Cocktail (ITEM 3) and the HRP-Streptavidin Concentrate (ITEM 4) vials should be briefly centrifuged (~1000 g) before opening to ensure maximum recovery and mixed well as precipitates may form during storage.

^{*1} vial is enough to test 2 membranes

ПЕМ	COMPONENT	PREPARATION	EXAMPLE			
1	Antibody Arrays	No Preparation	N/A			
2	Blocking Buffer	NoPreparation				
3	Biotinylated Antibody Cocktail*	Pipette 2 ml of Blocking Buffer into each vial. Mix gently with a pipette.	N/A			
4	1,000X HRP-Streptavidin Concentrate	Dilute 1,000-fold with Blocking Buffer. Mix gently with a pipette.	10 μl of 1,000X concentrate + 9990 μl of Blocking Buffer = 10 ml of 1X working solution			
5	20X Wash Buffer I Concentrate	Dilute each 20-fold with distilled or deionized	10 ml of 20X concentrate + 190 ml of water			
6	20X Wash Buffer II Concentrate	water.	=200mlof1X working solution			
7	2X Cell Lysis Buffer Concentrate**	Dilute 2-fold with distilled or deionized water.	10 ml of 2X concentrate + 10 ml of water = 20 ml of 1X working solution			
8	Detection Buffer C					
9	Detection Buffer D	No Preparation	N/A			
10	8-Well Incubation Tray w/ Lid					
11	Protease Inhibitor Cocktail	Pipette 60 µl of 1X Cell Lysis Buffer into the vial to prepare 100X Protease Inhibitor Cocktail concentrate.				
12	Phosphatase Inhibitor Cocktail II	Add 180 µl of 1X Lysis Buffer into the vial to prepare 25X Phosphatase Inhibitor Cocktail Set II Concentrate. Dissolve the powder thoroughly by gentle mixing.				

^{**}Only for use for preparing cell or tissue lysates. General tips for preparing lysates and other common sample types can be found on the online Resources Page

Note: Prior to preparing cell or tissue lysates: Add 20 μ l Protease Inhibitor Cocktail Concentrate (100X) and 80 μ l Phosphatase Inhibitor Cocktail Set II Concentrate (25X) into 1.9 ml 1X Lysis Buffer immediately before use. Mix well.

VIII. PROTOCOL

NOTE: Prepare all reagents and samples immediately prior to use. See Sections V and VII. <u>ALL</u> incubations and washes must be performed under gentle rotation/rocking (~0.5-1 cycle/sec)

- 1) Remove the kit from storage and allow the components to equilibrate to room temperature (RT).
- 2) Carefully remove the Antibody Arrays (ITEM 1) from the plastic packaging and place each membrane (printed side up) into a well of the Incubation Tray (ITEM 10). One membrane per well.

NOTE: The antibody printed side is marked by a dash (-) or number (#) in the upper left corner.

A. **Blocking**

- 3) Pipette 2 ml of Blocking Buffer (ITEM 2) into each well and incubate for 1 hour at RT.
- 4) Aspirate blocking buffer from each well with a pipette.

B. Sample Incubation

5) Pipette 1 ml of diluted or undiluted sample into each well and incubate for 1.5 to 5 hours at RT OR overnight at 4 °C.

NOTE: Longer incubations can help maximize the spot signal intensities. However, doing so can also increase the background response so complete liquid removal and washing is critical.

6) Aspirate samples from each well with a pipette.

C. First Wash

NOTE: The 20X Wash Buffer Concentrates I and II (ITEM 5 and 6) must be diluted 20-fold before use. See Section VII for details.

- 7) Wash Buffer I Wash: Pipette 2 ml of **1X** Wash Buffer I into each well and incubate for 5 minutes at RT. Repeat this 2 more times for a total of 3 washes using fresh buffer and aspirating out the buffer completely each time.
- 8) Wash Buffer II Wash: Pipette 2 ml of **1X** Wash Buffer II into each well and incubate for 5 minutes at RT. Repeat this 1 more time for a total of 2 washes using fresh buffer and aspirating out the buffer completely each time.

D. Biotinylated Antibody Cocktail Incubation

NOTE: The Biotinylated Antibody Cocktail (ITEM 3) must be prepared before use. See Section VII for details.

- 9) Pipette 1 ml of the **prepared** Biotinylated Antibody Cocktail into each well and incubate for 1.5 to 2 hours at RT OR overnight at 4°C.
- 10) Aspirate biotinylated antibody cocktail from each well.

E. Second Wash

11) Wash membranes as directed in Steps 7 and 8.

F. <u>HRP-Streptavidin Incubation</u>

NOTE: The 1,000X HRP-Streptavidin Concentrate (ITEM 4) must be diluted before use. See Section VII for details.

- 12) Pipette 2 ml of 1X HRP-Streptavidin into each well and incubate for 2 hours at RT OR overnight at 4°C.
- 13) Aspirate HRP-Streptavidin from each well.

G. Third Wash

14) Wash membranes as directed in Steps 7 and 8.

H. Chemiluminescence Detection

NOTE: Do not allow membranes to dry out during detection.

- 15) Transfer the membranes, printed side up, onto a sheet of chromatography paper, tissue paper, or blotting paper lying on a flat surface (such as a benchtop).
- 16) Remove any excess wash buffer by blotting the membrane edges with another piece of paper.
- 17) Transfer and place the membranes, printed side up, onto a plastic sheet (provided) lying on a flat surface.
- **NOTE:** Multiple membranes can be placed next to each other and fit onto a single plastic sheet. Use additional plastics sheets if necessary.
- 18) Into a single clean tube, pipette equal volumes (1:1) of Detection Buffer C (ITEM 8) and Detection Buffer D (ITEM 9). Mix well with a pipette.

EXAMPLE: 250 μ l of Detection Buffer C + 250 μ l of Detection Buffer D = 500 μ l (enough for 1 membrane)

- 19) Gently pipette 500 µl of the Detection Buffer mixture onto each membrane and incubate for 2 minutes at RT (DO NOT ROCK OR SHAKE). Immediately afterwards, proceed to Step 20.
- **NOTE:** Exposure should ideally start within 5 minutes after finishing Step 19 and completed within 10-15 minutes as chemiluminescence signals will fade over time. If necessary, the signals can usually be restored by repeating washing, HRP-Streptavidin and Detection Buffers incubations (Steps 11-19)

20) Place another plastic sheet on top of the membranes by starting at one end and gently "rolling" the flexible plastic sheet across the surface to the opposite end to smooth out any air bubbles. The membranes should now be "sandwiched" between two plastic sheets.

NOTE: Avoid "sliding" the top plastic sheet along the membranes' printed surface.

21) Transfer the sandwiched membranes to the chemiluminescence imaging system such as a CCD camera (recommended) and expose.

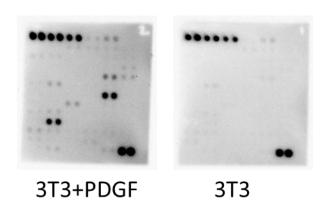
NOTE: Optimal exposure times will vary so performing multiple exposure times is strongly recommended. See Section VI for additional details.

I. Storage

22) To store, without direct pressure, gently sandwich the membranes between 2 plastic sheets (if not already), tape the sheets together or use plastic wrap to secure them, and store at \leq -20 °C for future reference.

IX. TYPICAL RESULTS

Typical results obtained with RayBio® C-Series Antibody Arrays



The preceding figures present typical images obtained with RayBio® C-Series Antibody Arrays. These membranes were probed with conditioned media from two different cell lines. Membranes were exposed with Kodak X-Omat® film at room temperature for 1 minute.

Note the strong signals of the Positive Control spots in the upper left and lower right corners. (See below for further details on the control spots.)

The signal intensity for each antigen-specific antibody spot is proportional to the relative concentration of the antigen in that sample. Comparison of signal intensities for individual antigen-specific antibody spots between and among array images can be used to determine relative differences in expression levels of each analyte sample-to-sample or group-to-group.

X. INTERPRETING THE RESULTS

A. Control Spots

<u>Positive Control Spots (POS)</u> – controlled amount of biotinylated antibody printed onto the array. Used for normalization and to orientate the arrays.

<u>Negative Control Spots (NEG)</u> – buffer printed (no antibodies) used to measure the baseline responses. Used for determining the level of non-specific binding of the samples.

Blank Spots (BLANK) – nothing is printed here. Used to measure the background response.

B. Data Extraction

Visual comparison of array images may be sufficient to see differences in relative protein expression. However, most researchers will want to perform numerical comparisons of the signal intensities (or more precisely, signal *densities*), using 2-D densitometry. Gel/Blot documentation systems and other chemiluminescent or phosphorescent detection systems are usually sold as a package with compatible densitometry software.

Any densitometry software should be sufficient to obtain spot signal densities from your scanned images. One such software program, ImageJ, is available for free from the NIH website along with an array plug-in.

We suggest using the following guidelines when extracting densitometry data from our array images:

- For each array membrane, identify a single exposure that the exhibits a high signal to noise ratio (strong spot signals and low background response). Strong Positive Control Spot signals but not too strong that that they are "bleeding" into one another is ideal. The exposure time does not need to be identical for each array, but Positive Control signals on each array image should have similar intensities.
- Measure the density of each spot using a circle that is roughly the size of one of the largest spots. Be sure to use the <u>same extraction circle</u> dimensions (area, size, and shape) for measuring the signal densities on every array for which you wish to compare the results.
- For each spot, use the <u>summed signal density</u> across the entire circle (ie, total signal density per unit area)

C. Data Analysis

NOTE: RayBiotech offers Microsoft® Excel-based Analysis Software Tools for each array kit for automatic analysis. Please visit the website at www.raybiotech.com or contact us for ordering information.

Once the raw numerical densitometry data is extracted, the background must be subtracted and the data normalized to the Positive Control signals to analyze.

<u>Background Subtraction</u>: Select values which you believe best represent the background. If the background is fairly even throughout the membrane, the Negative Control Spots (NEG) and/or Blank Spots (BLANK) should be similar and are accurate for this purpose.

<u>Positive Control Normalization:</u> The amount of biotinylated antibody printed for each Positive Control Spot is consistent from array to array. As such, the intensity of these Positive Control signals can be used to normalize signal responses for comparison of results across multiple arrays, much like housekeeping genes and proteins are used to normalize results of PCR gels and Western Blots, respectively.

To normalize array data, one array is defined as "Reference Array" to which the other arrays are normalized to. The choice of the Reference Array is arbitrary.

NOTE: The RayBio[®] Analysis Software Tools always designate Array 1/Sample 1 as the Reference Array.

Next, the simple algorithm below can be used to calculate and determine the signal fold expression between like analytes.

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

Where:

P1 = mean signal density of Positive Control spots on reference array P(y) = mean signal density of Positive Control spots on Array "y" X(y) = mean signal density for spot "X" on Array for sample "y" X(Ny)= normalized signal intensity for spot "X" on Array "y"

For example:

Let's determine the relative expression for IL-6 on two different arrays (Arrays 1 and 2). Let's assume that the duplicate signals for the IL-6 spots on each array are identical (or that the signal intensity used in the following calculation is the mean of the two duplicates spots). Also assume the following:

```
P1 = 2500
P2 = 2700
IL-6 (1) = 300
IL-6 (2) = 455
```

Then IL-6(N2) = 455 *2500/2700 = 421.30

The fold increase of IL-6(N2) vs IL-6(1) = 421.3/300 = 1.40-fold increase or a 40% increase in the signal intensity of IL-6 in Array 2 vs. Array 1.

XI. ARRAY MAP

	Α	В	С	D	E	F	G	Н	I	J	K	L
1	POS1	POS1	POS2	POS2	POS3	POS3	ABL1	ABL1	ACK1	ACK1	ALK	ALK
2	NEG	NEG	NEG	NEG	Axl	Axl	Blk	Blk	вмх	вмх	Btk	Btk
3	Csk	Csk	Dtk	Dtk	EGFR	EGFR	EphA1	EphA1	EphA2	EphA2	EphA3	EphA3
4	EphA4	EphA4	EphA5	EphA5	EphA6	EphA6	EphA7	EphA7	EphA8	EphA8	EphB1	EphB1
5	EphB2	EphB2	EphB3	EphB3	EphB4	EphB4	EphB6	EphB6	ErbB2	ErbB2	ErbB3	ErbB3
6	ErbB4	ErbB4	FAK	FAK	FER	FER	FGFR1	FGFR1	FGFR2	FGFR2	FGFR2 (α isoform)	FGFR2 (α isoform)
7	Fgr	Fgr	FRK	FRK	Fyn	Fyn	Hck	Hck	HGFR	HGFR	IGF-IR	IGF-IR
8	Insulin R	Insulin R	Itk	Itk	JAK1	JAK1	JAK2	JAK2	JAK3	JAK3	LCK	LCK
9	LTK	LTK	Lyn	Lyn	MATK	MATK	M-CSFR	M-CSFR	MUSK	MUSK	NGFR	NGFR
10	PDGFR- α	PDGFR- α	PDGFR- ß	PDGFR- ß	PYK2	PYK2	RET	RET	ROR1	ROR1	ROR2	ROR2
11	ROS	ROS	RYK	RYK	SCFR	SCFR	SRMS	SRMS	SYK	SYK	Tec	Tec
12	Tie-1	Tie-1	Tie-2	Tie-2	TNK1	TNK1	TRKB	TRKB	TXK	TXK	NEG	NEG
13	Tyk2	Tyk2	TYRO10	TYRO10	VEGFR2	VEGFR2	VEGFR3	VEGFR3	ZAP70	ZAP70	POS4	POS4

POS = Positive Control Spot NEG = Negative Control Spot BLANK = Blank Spot

NOTE: Protein alternative names, accession numbers, and official symbols can be accessed on www.raybiotech.com via the Resources Page.

XII. TROUBLESHOOTING GUIDE

PROBLEM	CAUSE	RECOMMENDATION				
	Chemiluminescent imager is not working properly	Contact image manufacturer				
No signals	Too Short Exposure	Expose the membranes longer				
(not even the positive controls spots)	Degradation of components due to improper storage	Store entire kit at ≤ - 20°C. Do not use kit after expiration date. See storage guidelines.				
	Improper preparation or dilution	Centrifuge vial briefly before use, mix well, and do not				
	of the HRP-Streptavidin	dilute more than 1000-fold				
	Waiting too long before exposing	The entire detection process should be completed in 10-15 minutes				
Positive controls spots	Low sample protein levels	Decrease sample dilution, concentrate samples, or load more protein initially				
signals visible but no other	Skipped Sample Incubation Step	Samples must be loaded after the blocking step				
spots	Too Short of Incubations	Ensure the incubations are performed for the appropriate time or try the optional overnight incubation(s)				
	Bubbles present on or below membrane	Don't rock/rotate the tray too vigorously or pipette the sample or reagent with excessive force				
Uneven signals and/or	Insufficient sample or reagent volume	Load enough sample and reagent to completely cover the membrane				
background	Insufficient mixing of reagents	Gently mix all reagents before loading onto the membrane, especially the HRP-Streptavidin and Biotin Antibody Cockta				
	Rocking/Rotating on an uneven surface while incubating	Rock/rotate on a flat surface or the sample or reagent can "pool" to one side				
	Too much HRP-Streptavidin or Biotinylated Antibody Cocktail	Prepare these signal enhancing components precisely as instructed				
	Membranes dried out	Do not let the membranes dry out during the experiment. Cover the incubation tray with the lid to minimize evaporation				
High background signals or all spots visible	Too High of Sample Protein Concentration	Increase dilution of the sample or load less protein				
	Exposed Too Long	Decrease exposure time				
	Insufficient Washing	Ensure all the wash steps are carried out and the wash buffer is removed completely after each wash step				
	Non-specific binding	Ensure the blocking buffer is stored and used properly.				

RayBio® is the trademark of RayBiotech, Inc.

The RayBio[®] Cytokine Antibody Array C-Series is patent-pending technology developed by RayBiotech.

This product is intended *for research only* and is not to be used for clinical diagnosis. Our produces 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 6 months from the date of purchase 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.

Kodak X-Omat is a registered trademark of Eastman Kodak Company. Microsoft Excel is a registered trademark of Microsoft Corporation.

Note:

This product is for research use only.



©2010 RayBiotech, Inc.