

RayBio[®] G-Series Human RTK Phosphorylation Antibody Array 1

For Simultaneously Detecting the Relative Level of Tyrosine
Phosphorylation of Human Receptor Tyrosine Kinases (RTKs)

User Manual

(Revised Oct. 14th, 2020)

(Cat#: AAH-PRTK-G1-4 and AAH-PRTK-G1-8)



**We Provide You with Quality Products
and Excellent Service**

Tel: (770) 729-2992 Fax: (888) 547-0580
Web: www.raybiotech.com Email: info@raybiotech.com



RayBiotech Life, Inc.

**RayBio® G-Series Human RTK Phosphorylation Antibody Array 1
Protocol**

TABLE OF CONTENTS

I.	Introduction.....	2
	How It Works.....	3
II.	Materials Provided.....	4
III.	Additional Materials Required.....	5
IV.	Reagent Preparation.....	6
V.	Overview and General Considerations.....	7
	A. Preparation of Samples.....	7
	B. Handling Glass Slides.....	8
	C. Incubation.....	8
VI.	Protocol.....	9
	A. Dry the Array Slides.....	9
	B. Blocking and Incubation.....	9
	C. Fluorescence Detection.....	12
VII.	Interpretation of Results.....	13
VIII.	Troubleshooting Guide.....	16
IX.	Reference List.....	17

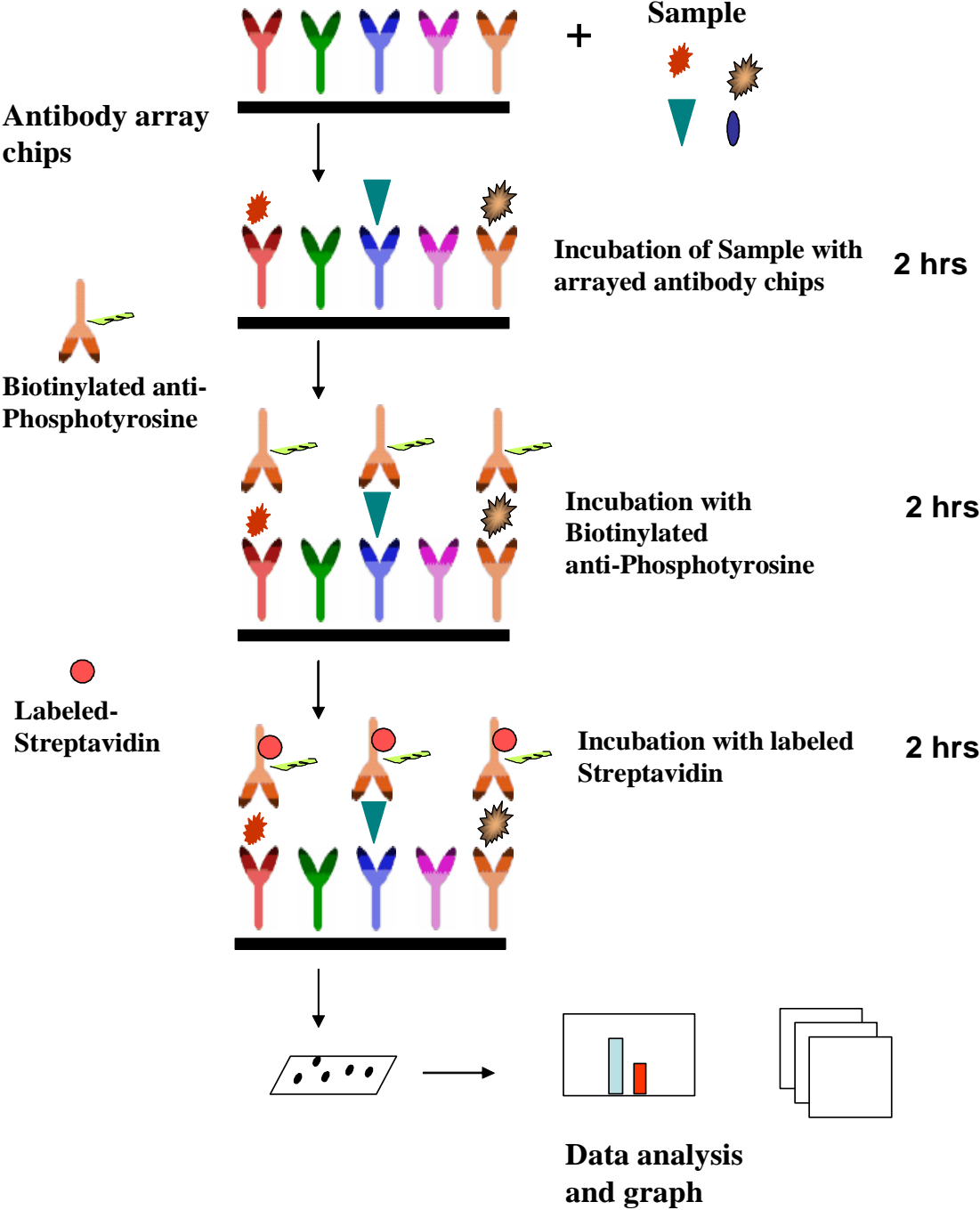
I. Introduction

Protein phosphorylation plays an unusually prominent role in cell signaling, development and growth. The RayBio® G-Series Human RTK Phosphorylation Antibody Array 1 is a very rapid, convenient, and sensitive assay that can simultaneously detect multiple protein phosphorylations and be used to monitor the 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 the RayBio® Human RTK Phosphorylation Antibody Array 1, which is specifically designed for simultaneous identification of the relative levels of phosphorylation of 71 different Human Receptor Tyrosine Kinases (RTKs) in cell lysate. 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 performing an analysis of immunoprecipitation and/or Western Blot.

To use the RayBio® G-Series Human RTK Phosphorylation Antibody Array 1, treated or untreated cell lysate is added into antibody array glass slide wells. The antibody array slide wells are washed, and biotinylated anti-phosphotyrosine antibodies are then used to detect the phosphorylated tyrosines on activated receptors. After incubation with a fluorescent dye-conjugated streptavidin (Cy3 equivalent), the slides can then be imaged using a laser scanner, such as the Axon GenePix, using the Cy3 channel.

Here's how it works



II. Materials Provided

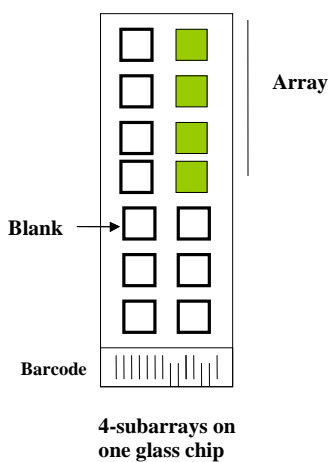
Upon receipt, the kit should be stored at $-20\text{ }^{\circ}\text{C}$ to $-80\text{ }^{\circ}\text{C}$. Please use within 6 months from the date of shipment. After initial use, the 2X Cell Lysis Buffer, Blocking Buffer, 20X Wash Buffer I, 20X Wash Buffer II, Biotin-Conjugated Anti-phosphotyrosine and Fluorescent dye-Conjugated Streptavidin (Cy3 equivalent) should be stored at $4\text{ }^{\circ}\text{C}$ to avoid repeated freeze-thaw cycles. The Array I Glass Slide, Protease Inhibitor Cocktail and Phosphatase Inhibitor Cocktail Set II should be kept at $-20\text{ }^{\circ}\text{C}$ to $-80\text{ }^{\circ}\text{C}$. Use within 3 months after initial use.

- RayBio® G-Series Human RTK Phosphorylation Antibody Array 1 Glass Slide with Frame (each slide contains 4 Subarrays, with 1 slide included for 4-subarray slides, and 2 slides included for 8-subarray slides)
- 2X Cell Lysis Buffer (10 ml)
- Protease Inhibitor Cocktail (1/2 tubes, 1 tube included for the 4-subarray slides, and 2 for the 8-subarray slides)
- Phosphatase Inhibitor Cocktail Set II (1/2 tubes, 1 tube included for the 4-subarray slides, and 2 for the 8-subarray slides)
- Blocking Buffer (8 ml)
- 20X Wash Buffer I (30 ml)
- 20X Wash Buffer II (30 ml)
- Biotin-Conjugated Anti-phosphotyrosine (1/2 tubes, 1 tube included for the 4-subarray slides, and 2 for the 8-subarray slides)
- Fluorescent dye-Conjugated Streptavidin (Cy3 equivalent) (1/2 tubes, 1 tube included for the 4-subarray slides, and 2 for the 8-subarray slides)
- Wash Buffer III (20 ml)
- Adhesive film

III. Additional Materials Required

- Shaker
- Laser scanner for fluorescence detection
- Aluminum foil
- Distilled water
- Plastic box
- 50 ml Centrifuge tube
- Isopropanol (2-propanol)

Layout of Array Glass Slide



IV. Reagent Preparation

- 1. Protease Inhibitor Cocktail:** Briefly spin down the Protease Inhibitor Cocktail vial before use. Add 60 μ l of 1X Cell Lysis Buffer to the vial to prepare a 100X Protease Inhibitor Cocktail Concentrate.
- 2. Phosphatase Inhibitor Cocktail Set II:** Briefly spin down the Phosphatase Inhibitor Cocktail Set II vial before use. Add 180 μ l of 1X Cell Lysis Buffer to the vial to prepare a 25X Phosphatase Inhibitor Cocktail Set II Concentrate. **Dissolve the powder thoroughly by gentle mixing.**
- 3. 2X Cell Lysis Buffer:** The 2X Cell Lysis Buffer should be diluted 2-fold with deionized or distilled water to prepare a 1X Cell Lysis Buffer solution. Then, add 20 μ l of the Protease Inhibitor Cocktail Concentrate and 80 μ l of the Phosphatase Inhibitor Cocktail Set II Concentrate into 1.9 ml of the 1X Cell Lysis Buffer to prepare a 1X Cell Lysis Buffer with Protease and Phosphatase Inhibitor Cocktail solution. Mix well before use.
- 4. 20X Wash Buffer I or II:** If the 20X Wash Buffer Concentrate contains visible crystals, warm to room temperature and mix gently until dissolved. Dilute 25 ml of the 20X Wash Buffer Concentrate into deionized or distilled water to yield 500 ml of 1X Wash Buffer.
- 5. Biotinylated anti-Phosphotyrosine:** Briefly spin down the Detection Antibody vial before use. Add 65 μ l of Blocking Buffer to the vial to prepare a Biotinylated Anti-phosphotyrosine Concentrate. Pipette up and down to mix gently (the Concentrate can be stored at 4 $^{\circ}$ C for 5 days). Add 30 μ l of Detection Antibody Concentrate to a tube with 570 μ l of Blocking Buffer to prepare a 1X Biotinylated Anti-phosphotyrosine solution. Mix gently.
- 6. Fluorescent dye-Conjugated Streptavidin (Cy3 equivalent):** Briefly spin down the Fluorescent dye-Conjugated Streptavidin vial before use. Add 50 μ l of Blocking Buffer to the vial to prepare a Streptavidin

Concentrate. Pipette up and down to mix gently. Add 10 μ l of Streptavidin Concentrate to a tube with 1 ml of Blocking Buffer to prepare a 1X Fluorescent dye-Conjugated Streptavidin solution. Mix gently.

V. Overview and General Considerations

A. Preparation of Samples

Cells can be prepared using the following convention.

For attached cells, remove the supernatant from the cell culture, and wash the cells twice with cold 1X PBS (for cells in suspension, pellet the cells by spinning down at 1500 rpm for 10 min). Make sure to remove any remaining PBS. Then, solubilize the cells at 2×10^7 cells/ml in the 1X Cell Lysis Buffer with Protease and Phosphatase Inhibitor Cocktail solution. Pipette up and down to resuspend the cells, and rock the lysates gently at 2–8 °C for 30 min. Transfer the lysates to microcentrifuge tubes and centrifuge at 14,000 x g for 5 min.

It is recommended that sample protein concentrations be determined using a total protein assay. For incubation with the Phosphorylation Antibody Array G-series 1, use cell lysates at a concentration of 50–1000 μ g/ml (as a starting point, we recommend using 400 μ g/ml of cell lysate diluted at least 5-fold with the Blocking Buffer).

Lysates should be used immediately or aliquoted and stored at –80 °C. Thawed lysates should be kept on ice prior to use.

If you experience high background, you may further dilute your sample. If signals are too weak, cell lysate can be pretreated by immunoprecipitation before incubation with the array slides. Immunoprecipitation can be done using an anti-phosphotyrosine antibody and Protein A.

B. Handling glass slides

- The microarray slides are very sensitive. Do not touch the array surface with tips, forceps or hands. Hold the slides by the edges only.
- Handle all buffers and slides with latex free gloves.
- Avoid breaking the glass slide.
- Maintain a clean environment.

C. Incubation

- Completely cover the array area with sample or buffer during incubation, and cover the incubation chamber with the adhesive film or plastic sheet protector to avoid drying.
- Avoid foaming during incubation steps.
- Perform all incubation and wash steps under gentle rotation.
- Cover the incubation chamber with the adhesive film during incubation, particularly when the incubation is more than 2 hours or less than 50 μ l of sample or reagent is used.
- Avoid cross-contamination from overflowing solution to neighboring wells.
- Several incubation steps such as step 2 (sample incubation), step 6 (Biotin-conjugated Anti-phosphotyrosine incubation) or step 9 (Fluorescent dye-Conjugated Streptavidin incubation) may be done at 4 $^{\circ}$ C overnight. Please make sure to cover the incubation chamber tightly to prevent evaporation.
- Avoid exposing the array slide to light from step 9 in page 12 on.

VI. Protocol

A. Dry the Glass Slide

Open the box containing the Glass Slide with Frame and take it out. Then let it air dry for 1 hour in a clean environment before use.

Note: Protect the slide from dust or other contaminants.

B. Blocking and Incubation

1. Add 100 μ l of 1X Blocking Buffer to each well and incubate at room temperature with gentle shaking for 30 min to block the slides. Make sure no bubbles are in the wells.

Note: Only add reagents to wells printed with antibodies.

2. Decant the Blocking Buffer from each well (make sure to remove all of the buffer). Add 100 μ l of each sample into appropriate wells. Incubate the arrays with sample at room temperature with gentle shaking for 2 hours or at 4 °C overnight.

*Note: We recommend using 100 μ l of cell lysate at a concentration of 50–1000 μ g/ml (as a starting point, we recommend using 400 μ g/ml cell lysate). **Dilute the lysate at least 5-fold with the Blocking Buffer. Make sure there are no bubbles in the wells.***

Note: The amount of sample used depends on the abundance of target proteins. More sample can be used if signals are too weak. If signals are too strong, the sample can be diluted further. The optimal sample dilution must be determined empirically by the researcher.

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

Note: Cell lysate can be pretreated by immunoprecipitation before incubation with the array slides if signals are too weak. Immunoprecipitation can be done using an anti-phosphotyrosine antibody and Protein A. The elution samples from Protein A can then be diluted with Blocking Buffer and incubated with the array slides.

3. Decant the samples from each well, and wash 3 times, 5 min per wash, with 100 μ l of 1X Wash Buffer I at room temperature with gentle shaking.

Note: Avoid the solution overflowing into neighboring wells.

4. Put the Glass Slide with Frame into a box with Wash Buffer I (cover the whole glass slide and frame with Wash Buffer I), and wash at room temperature with gentle shaking for 20 min.
5. Decant the Wash Buffer I from each well. Put the Glass Slide with Frame into a box with Wash Buffer II (cover the whole glass slide and frame with Wash Buffer II), and wash 2 times, 5 min per wash, at room temperature with gentle shaking.
6. Remove all of Wash Buffer II from each well. Add 100 μ l of the 1X Biotin-conjugated Anti-phosphotyrosine solution to each corresponding well. Incubate at room temperature with gentle shaking for 2 hours.
7. Decant the antibody solution and wash as directed in step 4 three times (wash 3 times, 20 min per wash).
8. Wash as directed in step 5.
9. Remove all of Wash Buffer II from each well. Add 100 μ l of the 1X Fluorescent dye-Conjugated Streptavidin solution to each subarray. Cover the incubation chamber with the Adhesive film. Cover the plate

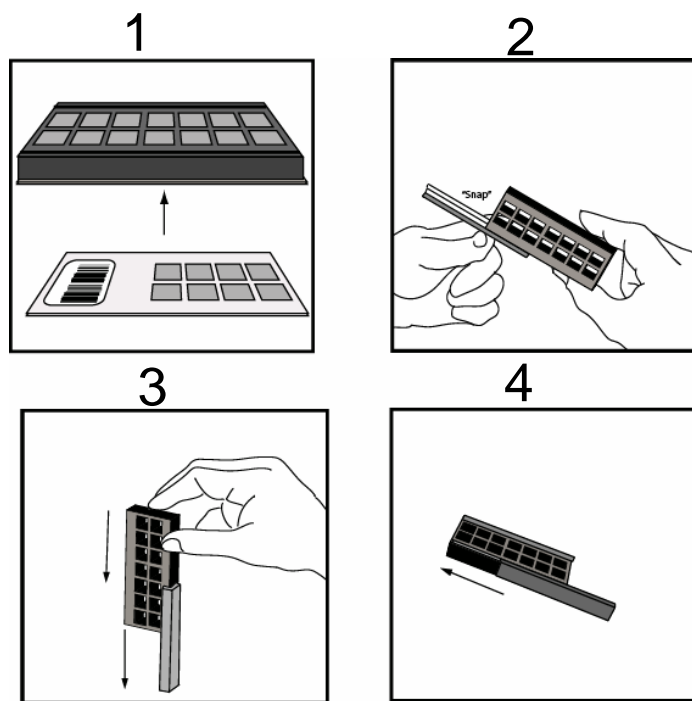
with aluminum foil to avoid exposure to light or incubate in a dark room.

Note: Avoid exposing the array slide to light from this step forward.

10. Incubate at room temperature with gentle shaking for 2 hours in the dark.

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

11. Decant the Fluorescent dye-Conjugated Streptavidin solution and disassemble the Glass Slide and Frame by removing the incubation frame and chamber from the slide as illustrated below.



Note: You may assemble and disassemble the glass slide into an incubation chamber and glass slide using the following steps.

1. To assemble, apply the incubation chamber to the slide with the printed side facing upward as illustrated in (1) above.

2. Gently snap one edge of a snap-on side as shown in (2).

3. *Adjust the position of the snap-on by gently pressing the edge of the snap-on side against a lab bench and pushing down as shown in (3).*

4. *Repeat steps 2 – 3 with a second snap-on as shown in (4).*

12. Gently put the glass slide into a 50 ml centrifuge tube or a plastic box with 40 ml of 1X Wash Buffer I as illustrated below. Gently roll or shake the tube for 5 min. Remove the Wash Buffer I. Repeat 2 more times for a total of 3 washes.



13. Wash the glass slide with 40 ml of Wash Buffer II for 5 min. Repeat one more time for a total of 2 washes.
14. Finally, wash the glass slide with 40 ml of deionized or distilled water.

C. Fluorescence Detection

1. To dry the glass slide, do one of the following:
- Put the glass slide into a 50 ml centrifuge tube and centrifuge at 1,000 rpm for 3 min
 - or*
 - Apply a compressed N₂ stream, or let glass slide air dry completely under clean air conditions (protected from light)

Make sure the slides are absolutely dry before scanning.

2. Image the slides using a laser scanner, such as the Axon GenePix, using the Cy3 channel.

Note: We recommend scanning the slides immediately after completing the experiment. Slides can also be stored at $-20\text{ }^{\circ}\text{C}$ in the dark for several days. If you do not have a laser scanner, we can scan and extract the data for free for you.

Note: Put the glass slide into a tube with 40 ml of 30% Wash Buffer III in isopropanol (add 15 ml of Wash Buffer III to a tube with 35 ml of isopropanol and mix well) and incubate for 10 min at room temperature if the background is not even or too high (cover the tube with aluminum foil to avoid exposure to light or incubate in a dark room). Dry the slide completely and re-scan the slide.

VII. Interpretation of Results

The following figure shows the RayBio[®] G-Series Human RTK Phosphorylation Antibody Array 1 probed with different cell lysates. The images were captured using a laser scanner. A biotinylated protein produces positive control signals, which can be used to identify the orientation of the slide and to normalize the results for comparison of different wells.

The antibody affinity to its target varies significantly between different antibodies. The fluorescence intensity detected on the array with each antibody depends on this affinity; therefore, the signal intensity comparison can only be performed within the same antibody/antigen system and not between different antibodies on the same slide. Certain proteins containing phosphorylated tyrosine may not be recognized by biotinylated anti-phosphotyrosine because of steric hindrance of the recognition site.

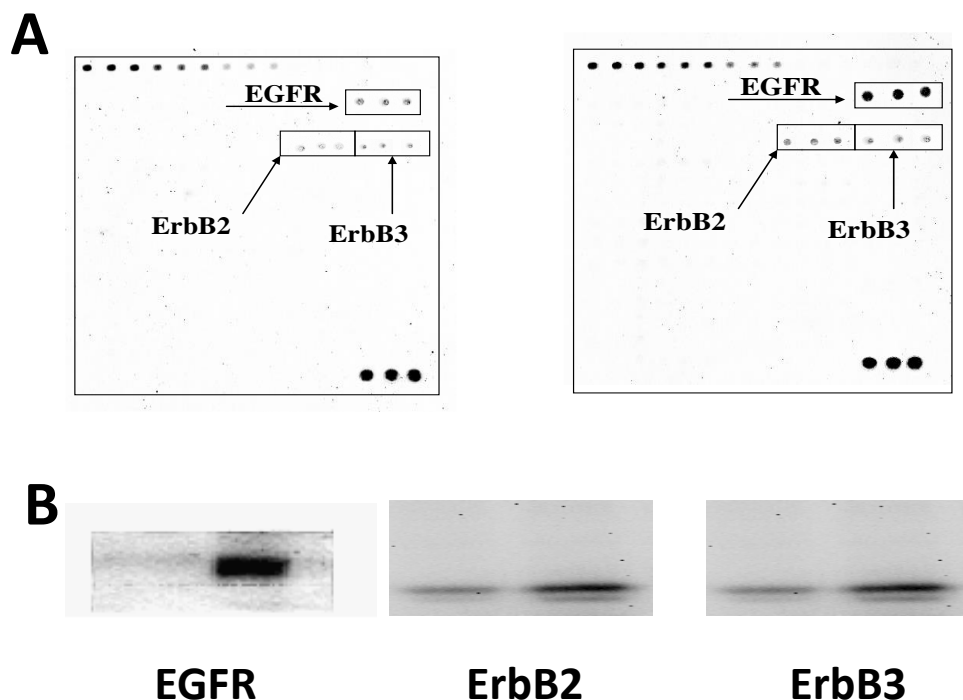


Fig 1. **(A)** Cells from a human epidermoid carcinoma cell line, A431, at 80-90% confluence were serum starved overnight, then exposed to 100 ng/ml EGF for 10 minutes at 37 °C. Control cells were serum starved without the subsequent stimulation with EGF. Cell lysates were prepared following the "Sample Preparation" portion IV of our protocol. To use the RayBio® G-Series Human RTK Phosphorylation Antibody Array 1, treated or untreated cell lysate was added to wells on the antibody array glass slides. The antibody array slides were washed and biotinylated anti-phosphotyrosine antibody was used to detect phosphorylated tyrosines on activated receptors. After incubation with Fluorescent dye-Conjugated Streptavidin, the signals were visualized by laser scanning using the cy3 channel. **(B)** Immunoprecipitation was done using anti-EGFR, ErbB2 and ErbB3 monoclonal antibodies and Protein A. Immunoblots were incubated with a biotinylated anti-phosphotyrosine monoclonal antibody to detect phosphorylated target protein receptors. Bands were visualized with Streptavidin-HRP followed by a chemiluminescent detection substrate.

RayBio® G-Series Human RTK Phosphorylation Antibody Array 1

Array Map

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

VIII. Troubleshooting Guide

Problem	Cause	Recommendation
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	Reduce sample dilution or concentrate sample
	Improper storage of kit	Store kit at suggested temperature
High background	Excess of biotinylated antibodies	Make sure to use the correct amount of antibodies
	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

IX. Reference List

1. Profiling receptor tyrosine kinase activation by using Ab microarrays. Nielsen UB, Cardone MH, Sinskey AJ, MacBeath G, and Sorger PK. **PNAS**. 2003;100(16):9330-9335.
2. A Prototype Antibody Microarray Platform to Monitor Changes in Protein Tyrosine Phosphorylation. Gembitsky DS, Lawlor K, Jacovina A, Yaneva M, and Tempst P. **Mol Cell Proteomics**. 2004;3:1102–1118.
3. Analysis of receptor signaling pathways by mass spectrometry: Identification of Vav-2 as a substrate of the epidermal and platelet derived growth factor receptors. Pandey A, Podtelejnikov AV, Blagoev B, Bustelo XR, Mann M, and Lodish HF. **PNAS**. 2000; 97(1);179–184.
4. Reduced T-cell and dendritic cell function is related to cyclooxygenase-2 overexpression and prostaglandin e(2) secretion in patients with breast cancer". Pockaj BA, Basu GD. **Annal Surg Oncol**. 2004;3:327-344.
5. Cytokine Antibody Arrays: A Promising Tool to Identify Molecular Targets for Drug Discovery. Huang RP. **Comb Chem High Throughput Screen**. 2003,;6:79-99.
6. Connexin suppresses human glioblastoma cell growth by down-regulation of monocyte chemotactic protein 1, as discovered using protein array technology. Huang R, Lin Y, Wang CC, J et al. **Cancer Res**. 2002;62:2806-2812.
7. Profiling of cytokine expression by biotin-labeled-based protein arrays. Lin Y, Huang R, Chen L-P, et al. **Proteomics**. 2003, 3: 1750-1757.
8. A novel method for high- throughput protein profiling from conditioned media and patient's sera. Huang RP, Huang R, Fan Y, and Lin Y. **Ana. Biochem**. 2001;294(1):55-62.

RayBio[®] Cytokine 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 Life, 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.

RayBio[®] is a registered trademark of RayBiotech Life, Inc.

This product is for research use only.



©2012 RayBiotech Life, Inc.