

RayBio[®] Human Cytokine Antibody Array G Series

Patent Pending Technology

User Manual (Revised 032407)

RayBio[®] Human Cytokine Antibody Array G Series 3 (Cat# AAH-CYT-G3-8)
RayBio[®] Human Cytokine Antibody Array G Series 5 (Cat# AAH-CYT-G5-8)
RayBio[®] Human Cytokine Antibody Array G Series 6 (Cat# AAH-CYT-G6-8)
RayBio[®] Human Cytokine Antibody Array G Series 7 (Cat# AAH-CYT-G7-8)
RayBio[®] Human Cytokine Antibody Array G Series 8 (Cat# AAH-CYT-G8-8)
RayBio[®] Human Inflammation Antibody Array 3 (Cat# AAH-INF-G3-8)
RayBio[®] Human growth factor Antibody Array 1 (Cat# AAH-GF-G1-8)

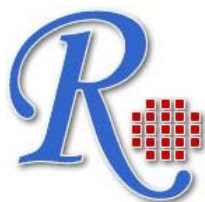
RayBio[®] Custom Human Cytokine Antibody Array (Cat# AAH-CUS-G)
RayBio[®] Human Cytokine Antibody Array Service (Cat# AAH-SER-G)



RayBiotech, Inc.

**We Provide You With Excellent
Protein Array Systems And Service**

**Tel:(Toll Free) 1-888-494-8555 or 770-729-2992; Fax: 1-888-547-0580;
Website:www.raybiotech.com Email: info@raybiotech.com**



RayBiotech, Inc.

RayBio[®] Human Cytokine Antibody Array G Series Protocol

TABLE OF CONTENTS

I.	Introduction.....	2
	How It Works.....	4
II.	Materials Provided.....	5
	Additional Materials Required.....	5
III.	Overview and General Considerations.....	6
	A. Preparation of Samples.....	6
	B. Handling Glass Chips	6
	C. Incubation.....	7
IV.	Protocol.....	7
	A. Blocking and Incubation.....	7
	B. Detection.....	10
V.	Interpretation of Results.....	11
VI.	Troubleshooting Guide.....	16
VII.	Reference List.....	17

Cytokine Antibody Arrays are RayBiotech patent-pending technology.

RayBio[®] is the trademark of RayBiotech, Inc.

I. Introduction

All cell functions, including cell proliferation, cell death and differentiation, as well as maintenance of health status and development of disease, are controlled by many genes and signaling pathways. New techniques such as cDNA microarrays have enabled us to analyze the global gene expression¹⁻³. However, almost all cell functions are executed by proteins, which cannot be studied by DNA and RNA alone. Experimental analysis clearly shows a disparity between the relative expression levels of mRNA and their corresponding proteins⁴. Therefore, it is critical to analyze the protein profile. Currently, two-dimensional polyacrylamide SDS page coupled with mass spectrometry is the mainstream approach to analyzing multiple protein expression levels^{5,6}. However, the requirement of sophisticated devices and the lack of quantitative measurements greatly limit its broad application. Thus, no simple, cost effective, and rapid method of analysis of multiple protein expression levels has been available to researchers until now.

Our RayBio[®] Human Cytokine Antibody Array is the first commercially available protein array system⁷⁻¹¹. By using the RayBiotech system, scientists can rapidly and accurately identify the expression profiles of multiple cytokines in several hours inexpensively.

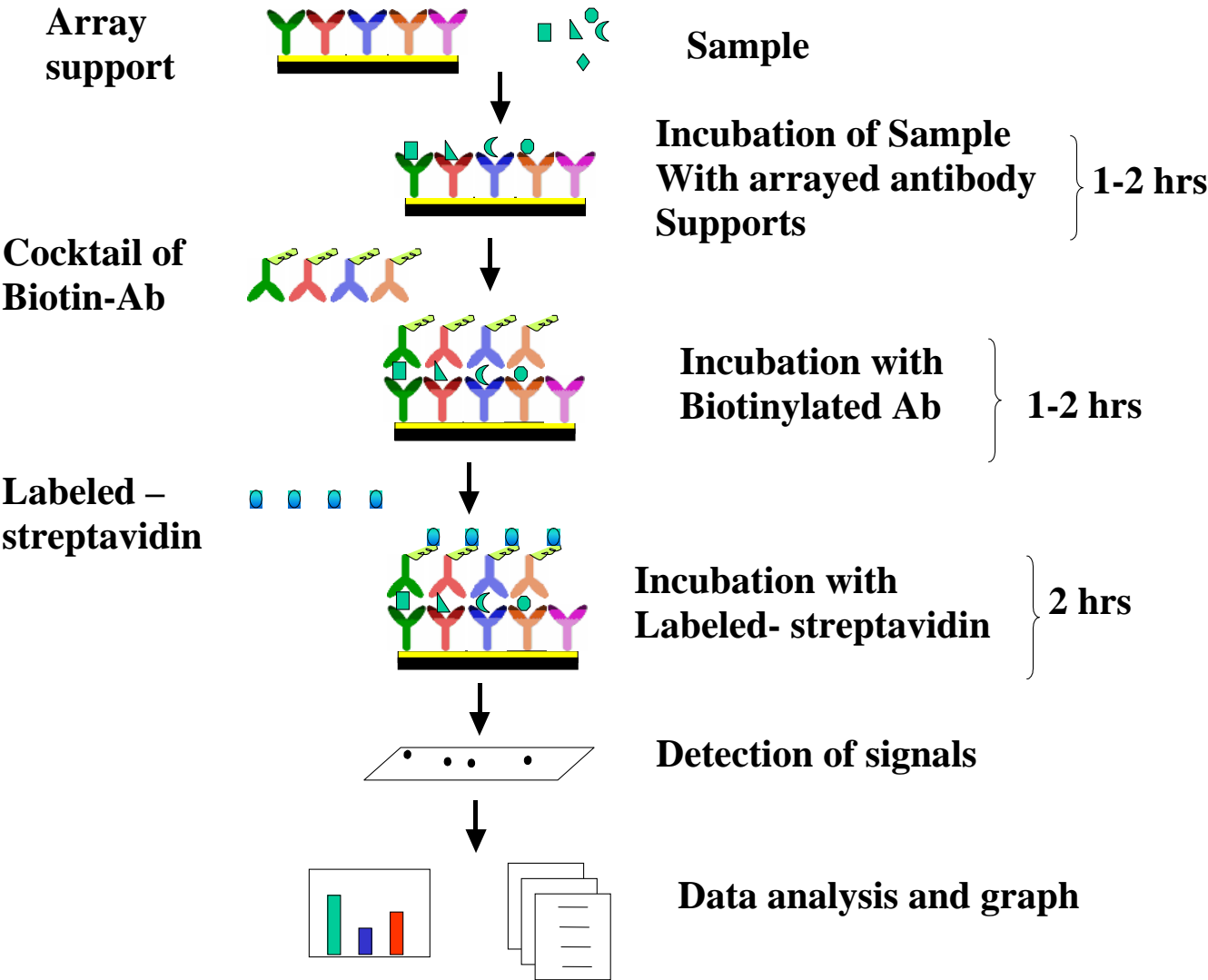
The RayBiotech kit (G series) is a glass slide format. The kit provides a highly sensitive approach to simultaneously detect multiple cytokine expression levels from cell culture supernatant, patient's serum, tissue lysate and other sources. The arrays are manufactured using non-contact arrayer. The experimental procedure is simple and can be performed in any laboratory. The signals from G series arrays are detected using a laser scanner.

Besides the products listed in this manual, RayBiotech also provides RayBio[®] Human Cytokine Antibody Array G series 1000 for detection of 120 human cytokines in single experiment and RayBio[®] Human Cytokine Antibody Array G series 2000 for detection of 174 human cytokines in single experiment.

Pathway-specific array systems allow investigators to focus on the specific problem and are becoming an increasingly powerful tool in cDNA microarray system. RayBiotech's first protein array system, known as RayBio[®] Human Cytokine Antibody Array, is particularly useful compared with the human cytokine cDNA microarray system. Besides the ability to detect protein expression, RayBiotech's system is a more accurate reflection of active cytokine levels because it only detects secreted cytokines, and no amplification step is needed. Furthermore, it is much simpler, faster, environmentally friendlier, and more sensitive.

Simultaneous detection of multiple cytokines undoubtedly provides a powerful tool to study cytokines. Cytokines play an important role in innate immunity, apoptosis, angiogenesis, cell growth and differentiation¹². Cytokines are involved in most disease processes, including cancer and cardiac diseases. The interaction between cytokines and the cellular immune system is a dynamic process. The interactions of positive and negative stimuli, and positive as well as negative regulatory loops are complex and often involve multiple cytokines.

Here's how it works



II. Materials Provided

Upon receipt, all components of the RayBio[®] Human Cytokine Antibody Array kit should be stored at -20⁰C. At -20⁰C the kit will retain complete activity for up to 6 months. Once thawed, the glass chips, Alexa-Flour 555-streptavidin, internal control and 2X Blocking Buffer should be kept at – 20⁰C and all other component should be stored at 4⁰C. Use within three months after reagents have been thawed. Please use within six months of purchase.

- RayBio[®] Human Cytokine Antibody Microarray slides (8 subarrays in each glass chip)
- Biotin-Conjugated Anti-Cytokines (2 tubes, each tube for 4 wells)
- 1,500X Alexa Flour 555-Conjugated Streptavidin (1 tube)
- 2X Blocking Buffer (8ml)
- 20X Wash Buffer I (60ml)
- 20X Wash Buffer II (40ml)
- Internal control (powder, 1 tube)
- 2X Cell Lysis Buffer (16 ml)
- Manual

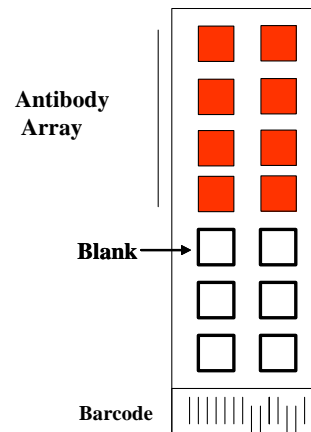
Additional Materials Required

Accessory

RayBio[®] G series antibody array accessory (including slide incubation chamber, Gasket, Protective cover, Snap-on sides and adhesive film) (Cat# G-CFF).

- Orbital shaker
- Laser scanner for fluorescence detection
- Aluminum foil
- Distilled water
- Plastic box

Layout of G series



8 arrays in one glass chip

III. Overview and General Considerations

A. Preparation of Samples

- Use serum-free conditioned media if possible.
- If serum-containing conditioned media is required, use serum as control since many types of sera contain cytokines.
- For cell lysates and tissue lysates, we recommend using 1X Cell Lysis Buffer to extract proteins from cell or tissue (e.g. using homogenizer). After extraction, spin the sample and save supernatant for experiment. Determine protein concentration. Dilute 2X Cell Lysis Buffer with H₂O (we recommend adding proteinase inhibitors to Cell Lysis Buffer before use). Prepare relative concentrated lysate since we recommend diluting lysate at least 10 folds with Blocking Buffer for array assay.
- We recommend using
 - 50 to 100 µl of cell culture supernatant
 - or 50 to 100 µl of original or 5-fold diluted serum or plasma
 - or 20-200 µg of protein for cell lysates and tissue lysates.

If you experience high background, you may further dilute your sample.

B. Handling glass chips

- The microarray slides are sensitive, do not touch the surface. Grip the slides by the edges only.
- Handle all buffers and slides with latex free gloves.
- Avoid breaking glass slide.
- Handle glass chip in clean environment.

C. Incubation

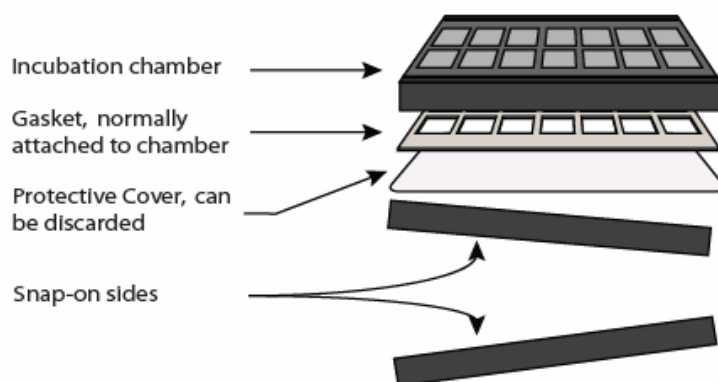
- Completely cover array area with sample or buffer during incubation, and cover the incubation chamber with adhesive film or plastic sheet protector to avoid drying.
- Avoid foaming during incubation steps.
- Perform all incubation and wash steps under gentle rotation.
- Cover the incubation chamber with adhesive film during incubation, particularly when incubation is more than 2 hours or 50 μ l of sample or reagent is used.
- Avoid cross-contamination from overflowing solution to neighboring wells.
- Several incubation steps such as step 3 (blocking), step 4 (sample incubation), step 8 (biotin-Ab incubation) or step 11 (Alexa Flour 555-streptavidin incubation) may be done at 4⁰C for overnight. Please make sure to cover the incubation chamber tightly to prevent evaporation.

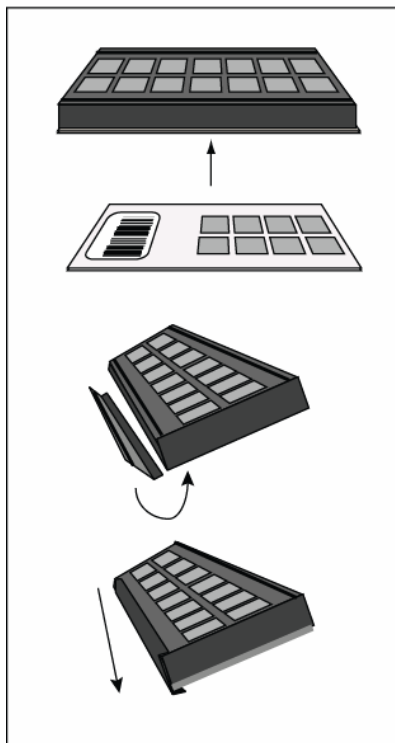
IV. Protocol

A. Blocking and Incubation

1. Take the glass chip out from the box. Let air dry for 60 minutes.
2. Assemble the glass chip into incubation chamber and incubation frame as shown below.

Instructions for incubation chamber assembly ***G Series and Q series arrays***





1 Carefully place slide at bottom of the chamber as shown. The slide will adhere somewhat to the bottom. Warning: the slide is fragile, so do not apply more than gentle force to the apparatus.

2 While gently holding chamber and slide, place side on chamber as shown, beginning with bottom flap first.

3 Then, press the top of the side into groove on chamber, and then apply even, gentle pressure from one end to the other. Repeat this procedure with the other side.

3. Add 100 μ l 1 X Blocking Buffer into each well and incubate at room temperature for 30 min to block slides. Dilute 2X Blocking Buffer with H₂O. Make sure no bubbles are in the well.

Note: only add reagents to wells printed with antibodies.

4. Decant Blocking Buffer from each well, and incubate arrays with sample at room temperature for 1 to 2 hours. Dilute sample using 1X Blocking Buffer if necessary. We strongly recommend including Internal Control in your assay, add 100 μ l of Blocking Buffer to IC tube, mix well and transfer 1 μ l of IC to each well (50 to 100 μ l of sample).

*Note: We recommend using 50 to 100 μ l of conditioned media or 50 to 100 μ l of original or 2-5 fold diluted serum or plasma or 10-200 μ g of protein for cell lysates and tissue lysates. **Dilute the lysate at least 10 folds with 1 X blocking buffer to make a total volume of 50 to 100 μ l. Make sure there is no bubble in the wells.***

Note: The amount of sample used depends on the abundance of cytokines. More of the sample can be used if signals are too weak. If signals are too strong, the sample can be diluted further.

Note: Incubation may be done at 4⁰C for overnight.

Note: when transfer IC, use 0.1 µl to 2 µl pipettor.

5. Decant the samples from each well, and wash 5 times with 150 µl of 1X Wash Buffer I at room temperature with gentle shaking. 2 min per wash. Dilute 20X Wash Buffer I with H₂O. Completely remove wash buffer I in each wash step.

Note: avoid solution flowing into neighboring wells.

6. Wash 2 times with 150 µl of 1X Wash Buffer II at room temperature with shaking. 2 min per wash. Dilute 20X Wash Buffer II with H₂O. Completely remove wash buffer II in each wash step.
7. Prepare working solution for biotin-conjugated antibodies. After brief spinning,

Add 300 µl µl of 1x blocking buffer to the Biotin-Conjugated Antibody tube. Mix gently.

Note: the diluted biotin-conjugated antibodies can be stored at 4⁰C for 2-3 days.

8. Add 70 µl of diluted biotin-conjugated antibodies to each corresponding well. Incubate at room temperature for 2 hours.

Note: incubation may be done at 4⁰C for overnight.

9. Wash as directed in steps 5 and 6.

10. Add 70 µl of 1,500 fold diluted Alexa Flour 555-conjugated streptavidin (after brief spinning, add 1.5 ml of Blocking Buffer to

Alexa Flour 555-conjugated streptavidin tube) to each subarray. Cover the incubation chamber with Adhesive film. Cover the plate with aluminum foil to avoid exposure to light or incubate in dark room.

11. Incubate at room temperature for 1 to 2 hours.

Note: incubation may be done at 4⁰C for overnight.

12. Wash with Wash Buffer I **twice** as directed in steps 5.

B. Fluorescence Detection

1. Decant excess Wash Buffer from wells.

2. Disassemble the slide out of the incubation frame and chamber.

3. Place the whole slide in 50 ml centrifuge tube, add enough Wash Buffer I (about 30 ml) to cover the whole slide and gently shake at room temperature for 10 minutes. Decant Wash Buffer I. Repeat Wash Buffer I once. Wash with Wash Buffer II (about 30 ml) with gentle shake at room temperature for 10 minutes. Or wash using slide chamber. Rinse the slide with distilled H₂O.

4. Remove water droplets by centrifuge at 1,000 rpm for 3 minutes and then let slide dry completely in air at least 20 minutes (protect from light). Make sure the slides are absolutely dry before the scanning procedure.

5. Image the signals using laser scanner such Axon GenePix using cy3 channel.

Note: we recommend scanning slides right after experiment. You also can store the slide at -20⁰C in dark for several days. If you do not have a laser scanner, we can provide service for you. Just simply send your slide to us and we will take care of it.

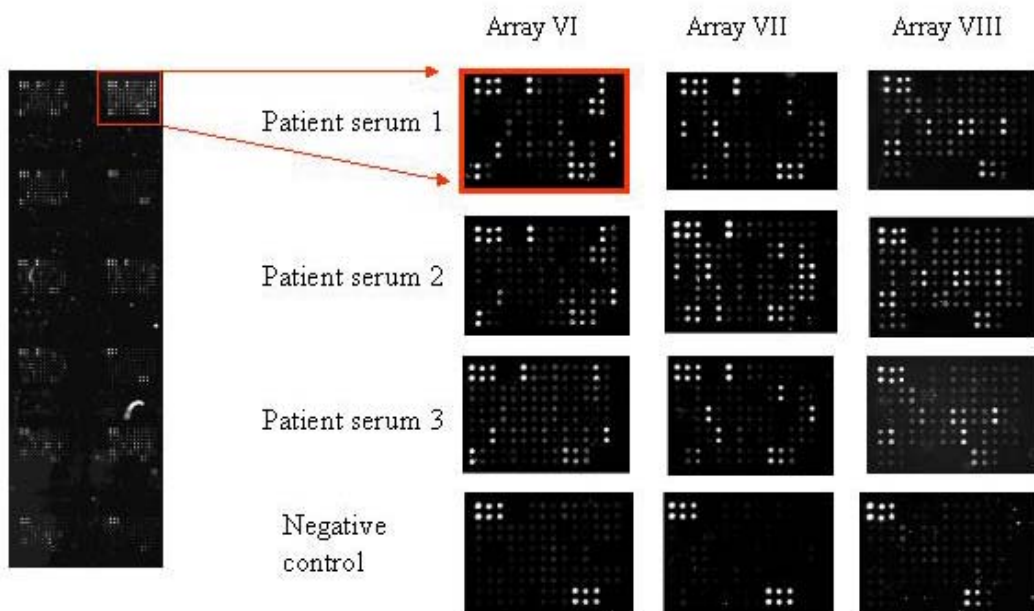
V. Interpretation of Results:

The following figure shows RayBio® Human Cytokine Antibody Array G series 2000 probed with different cell culture supernatant. The images were captured using laser scanner. The biotin-conjugated protein produces positive signals, which can be used to identify the orientation and to compare the relative expression levels among the different wells. The internal control (IC) can also be used to normalize the signal intensities among array membranes in different experiments.

The signal intensities obtained from laser scanner can simply be imported into our analysis tool. The analysis tool will help you:

- Locate your signal intensities to antibody array map
- Link the protein to website for more detailed information on the particular protein
- Protein list sorting
- Average signal intensities
- Subtract background
- Normalize the data from different samples
- Obtain protein level comparison charts among different samples

This analysis tool is very simple and affordable, which will not only assist in compiling and organizing your data, but also reduces your calculations to a “copy and paste” step.



If you do not use our **RayBio® Analysis Tool**, you can locate the cytokines by referring to corresponding RayBio® Human cytokine Antibody Array G series map.

Normalization and comparison

For biomarker discovery or for analysis of large number of arrays, great attention must be paid to the normalization. Our antibody array design includes several controls for normalization and comparison of arrays performing in different membranes and different experiments (for more information please read the reference 17).

Positive control. Positive control is biotinylated protein. It can be used to normalize the streptavidin incubation step. If the positive signals from different array membranes are similar, positive control is a simple and effective way for normalization.

Internal control. RayBio® antibody arrays also include spiking-in protein serving as internal control (IC). The spiking-in proteins do not have cross-reactivity with protein in the array. It can be used to normalize the entire process.

Negative control. Negative control is BSA. Normally, it should only give a background reading.

RayBio® Human Cytokine Antibody Array G series 3 (42)

	a	b	c	d	e	f	g	h	i	j	k	l	J
1	Pos 1	Pos 2	Pos 3	Neg	Neg	ENA-78	GCSF	GM-CSF	GRO	GRO-α	I-309	IL-1α	IL-1 β
2	Pos 1	Pos 2	Pos 3	Neg	Neg	ENA-78	GCSF	GM-CSF	GRO	GRO-α	I-309	IL-1α	IL-1 β
3	IL-2	IL-3	IL-4	IL-5	IL-6	IL-7	IL-8	IL-10	IL-12 p40p70	IL-13	IL-15	IFN-γ	MCP-1
4	IL-2	IL-3	IL-4	IL-5	IL-6	IL-7	IL-8	IL-10	IL-12 p40p70	IL-13	IL-15	IFN-γ	MCP-1
5	MCP-2	MCP-3	MCSF	MDC	MIG	MIP-1 δ	RANTES	SCF	SDF-1	TARC	TGF- β1	TNF-α	TNF-β
6	MCP-2	MCP-3	MCSF	MDC	MIG	MIP-1 δ	RANTES	SCF	SDF-1	TARC	TGF- β1	TNF-α	TNF-β
7	EGF	IGF-I	Angiogenin	Oncostatin M	Thrombopoietin	VEGF	PDGF BB	Leptin	IC 1	IC 2	IC 3	Neg	Neg
8	EGF	IGF-I	Angiogenin	Oncostatin M	Thrombopoietin	VEGF	PDGF BB	Leptin	IC 1	IC 2	IC 3	Neg	Neg

RayBio® Human Cytokine Antibody Array 5 (79)

	A	B	C	D	E	F	G	H	I	J	K	L	M
1	Pos 1	Pos 2	Pos 3	Neg	Neg	Neg	ENA-78	GCSF	GM-CSF	GRO	GRO-α	I-309	IL-1α
2	Pos 1	Pos 2	Pos 3	Neg	Neg	Neg	ENA-78	GCSF	GM-CSF	GRO	GRO-α	I-309	IL-1α
3	IL-1 β	IL-2	IL-3	IL-4	IL-5	IL-6	IL-7	IL-8	IL-10	IL-12 p40p70	IL-13	IL-15	IFN- γ
4	IL-1 β	IL-2	IL-3	IL-4	IL-5	IL-6	IL-7	IL-8	IL-10	IL-12 p40p70	IL-13	IL-15	IFN- γ
5	MCP-1	MCP-2	MCP-3	MCSF	MDC	MIG	MIP-1β	MIP-1δ	RANTES	SCF	SDF-1	TARC	TGF-β1
6	MCP-1	MCP-2	MCP-3	MCSF	MDC	MIG	MIP-1β	MIP-1δ	RANTES	SCF	SDF-1	TARC	TGF-β1
7	TNF-α	TNF-β	EGF	IGF-I	Angiogenin	Oncostatin M	Thrombopoietin	VEGF	PDGF-BB	Leptin	BDNF	BLC	Ck β 8-1
8	TNF-α	TNF-β	EGF	IGF-I	Angiogenin	Oncostatin M	Thrombopoietin	VEGF	PDGF-BB	Leptin	BDNF	BLC	Ck β 8-1
9	Eotaxin	Eotaxin-2	Eotaxin-3	FGF-4	FGF-6	FGF-7	FGF-9	Flt-3 Ligand	Fractalkine	GCP-2	GDNF	HGF	IGFBP-1
10	Eotaxin	Eotaxin-2	Eotaxin-3	FGF-4	FGF-6	FGF-7	FGF-9	Flt-3 Ligand	Fractalkine	GCP-2	GDNF	HGF	IGFBP-1
11	IGFBP-2	IGFBP-3	IGFBP-4	IL-16	IP-10	LIF	LIGHT	MCP-4	MIF	MIP-3 α	NAP-2	NT-3	NT-4
12	IGFBP-2	IGFBP-3	IGFBP-4	IL-16	IP-10	LIF	LIGHT	MCP-4	MIF	MIP-3 α	NAP-2	NT-3	NT-4
13	Osteoprotegerin	PARC	PIGF	TGF- β 2	TGF- β 3	TIMP-1	TIMP-2	Neg	IC1	IC2	IC3	Neg	Neg
14	Osteoprotegerin	PARC	PIGF	TGF- β 2	TGF- β 3	TIMP-1	TIMP-2	Neg	IC1	IC2	IC3	Neg	Neg

RayBio® Human Cytokine Antibody Array G series 6 (60)

	a	b	c	d	e	f	g	h	i	j	k	l	m	n
1	POS 1	POS 2	POS 3	NEG	NEG	Angiogenin	BDNF	BLC	BMP-4	BMP-6	CK b 8-1	CNTF	EGF	Eotaxin
2	POS 1	POS 2	POS 3	NEG	NEG	Angiogenin	BDNF	BLC	BMP-4	BMP-6	CK b 8-1	CNTF	EGF	Eotaxin
3	Eotaxin-2	Eotaxin-3	FGF-6	FGF-7	Fit-3 Ligand	Fractalkine	GCP-2	GDNF	GM-CSF	I-309	IFN-gamma	IGFBP-1	IGFBP-2	IGFBP-4
4	Eotaxin-2	Eotaxin-3	FGF-6	FGF-7	Fit-3 Ligand	Fractalkine	GCP-2	GDNF	GM-CSF	I-309	IFN-gamma	IGFBP-1	IGFBP-2	IGFBP-4
5	IGF-I	IL-10	IL-13	IL-15	IL-16	IL-1α	IL-1β	IL-1ra	IL-2	IL-3	IL-4	IL-5	IL-6	IL-7
6	IGF-I	IL-10	IL-13	IL-15	IL-16	IL-1α	IL-1β	IL-1ra	IL-2	IL-3	IL-4	IL-5	IL-6	IL-7
7	Leptin	LIGHT	MCP-1	MCP-2	MCP-3	MCP-4	M-CSF	MDC	MIG	MIP-1delta	MIP-3α	NAP-2	NT-3	PARC
8	Leptin	LIGHT	MCP-1	MCP-2	MCP-3	MCP-4	M-CSF	MDC	MIG	MIP-1delta	MIP-3α	NAP-2	NT-3	PARC
9	PDGF-BB	RANTES	SCF	SDF-1	TARC	TGF-beta1	TGF-beta 3	TNF-alpha	TNF-beta	IC1	IC2	IC3	NEG	NEG
10	PDGF-BB	RANTES	SCF	SDF-1	TARC	TGF-beta1	TGF-beta 3	TNF-alpha	TNF-beta	IC1	IC2	IC3	NEG	NEG

RayBio® Human Cytokine Antibody Array G series 7 (60)

	a	b	c	d	e	f	g	h	i	j	k	l	m	n
1	POS1	POS2	POS3	NEG	NEG	Aarp30	AgRP	Angiopoietin-2	Amphiregulin	Axl	bFGF	b-NGF	BTC	CCL-28
2	POS1	POS2	POS3	NEG	NEG	Aarp30	AgRP	Angiopoietin-2	Amphiregulin	Axl	bFGF	b-NGF	BTC	CCL-28
3	CTACK	Dtk	EGF-R	ENA-78	Fas/TNFRSF6	FGF-4	FGF-9	GCSF	GTR-Ligand	GTR	GRO	GRO-alpha	HCC-4	HGF
4	CTACK	Dtk	EGF-R	ENA-78	Fas/TNFRSF6	FGF-4	FGF-9	GCSF	GTR-Ligand	GTR	GRO	GRO-alpha	HCC-4	HGF
5	ICAM1	ICAM3	IGFBP-3	IGFBP-6	IGF-1 SR	IL-1 R4/ST2	IL-1 RI	IL-11	IL-12 p40	IL-12 p70	IL-17	IL-2 Ralpha	IL-6 R	IL-8
6	ICAM1	ICAM3	IGFBP-3	IGFBP-6	IGF-1 SR	IL-1 R4/ST2	IL-1 RI	IL-11	IL-12 p40	IL-12 p70	IL-17	IL-2 Ralpha	IL-6 R	IL-8
7	I-TAC	Lymphotoxin	MF	MP-1alpha	MP-1beta	MP-3beta	MSP-alpha	NT-4	Osteoprotegerin	Oncostatin M	PIGF	sgp130	sTNF RI	sTNF-RI
8	I-TAC	Lymphotoxin	MF	MP-1alpha	MP-1beta	MP-3beta	MSP-alpha	NT-4	Osteoprotegerin	Oncostatin M	PIGF	sgp130	sTNF RI	sTNF-RI
9	TECK	TIMP-1	TIMP-2	Thrombopoietin	TRAIL R3	TRAIL R4	uPAR	VEGF	VEGF-D	IC1	IC2	IC3	NEG	NEG
10	TECK	TIMP-1	TIMP-2	Thrombopoietin	TRAIL R3	TRAIL R4	uPAR	VEGF	VEGF-D	IC1	IC2	IC3	NEG	NEG

RayBio® Human Cytokine Antibody Array G series 8 (54)

	a	b	c	d	e	f	g	h	i	j	k	l	m	n
1	POS 1	POS 2	POS 3	NEG	NEG	Activin A	ALCAM	B7-1(CD80)	BMP-5	BMP-7	Cardiotrophin ₁	CD14	CXCL-16	DR6 (TNFRSF21)
2	POS 1	POS 2	POS 3	NEG	NEG	Activin A	ALCAM	B7-1(CD80)	BMP-5	BMP-7	Cardiotrophin ₁	CD14	CXCL-16	DR6 (TNFRSF21)
3	Endoglin	ErbB3	E-Selectin	Fas Ligand	ICAM-2	IGF-II	IL-1 R II	IL-10 Rbeta	IL-13 Ralpha2	IL-18 BPalpha	IL-18 Rbeta	MMP-3	IL-2 Rbeta	IL-2 Rgamma
4	Endoglin	ErbB3	E-Selectin	Fas Ligand	ICAM-2	IGF-II	IL-1 R II	IL-10 Rbeta	IL-13 Ralpha2	IL-18 BPalpha	IL-18 Rbeta	MMP-3	IL-2 Rbeta	IL-2 Rgamma
5	IL-21R	IL-5 Ralpha	IL-9	IP-10	LAP	Leptin R	LIF	L-Selectin	M-CSF R	MMP-1	MMP-13	MMP-9	MPIF-1	NGF R
6	IL-21R	IL-5 Ralpha	IL-9	IP-10	LAP	Leptin R	LIF	L-Selectin	M-CSF R	MMP-1	MMP-13	MMP-9	MPIF-1	NGF R
7	PDGF AA	PDGF-AB	PDGF Ralpha	PDGF Rbeta	PECAM-1	Prolactin	SCF R	SDF-1beta	Siglec-5	TGF-alpha	TGF beta2	Tie-1	Tie-2	TIMP-4
8	PDGF AA	PDGF-AB	PDGF Ralpha	PDGF Rbeta	PECAM-1	Prolactin	SCF R	SDF-1beta	Siglec-5	TGF-alpha	TGF beta2	Tie-1	Tie-2	TIMP-4
9	VE-Cadherin	VEGF R2	VEGF R3	Neg	Neg	Neg	Neg	Neg	Neg	IC1	IC2	IC3	NEG	NEG
10	VE-Cadherin	VEGF R2	VEGF R3	Neg	Neg	Neg	Neg	Neg	Neg	IC1	IC2	IC3	NEG	NEG

RayBio® Human Inflammation Antibody Array G series 3 (40)

	A	B	C	D	E	F	G	H	I	J	K	L	M
1	POS 1	POS 2	POS 3	NEG	NEG	EOTAXIN	EOTAXIN-2	GCSF	GM-CSF	ICAM-1	IFN-γ	I-309	IL-1α
2	POS 1	POS 2	POS 3	NEG	NEG	EOTAXIN	EOTAXIN-2	GCSF	GM-CSF	ICAM-1	IFN-γ	I-309	IL-1α
3	IL-1β	IL-2	IL-3	IL-4	IL-6	IL-6sR	IL-7	IL-8	IL-10	IL-11	IL-12 p40	IL-12 p70	IL-13
4	IL-1β	IL-2	IL-3	IL-4	IL-6	IL-6sR	IL-7	IL-8	IL-10	IL-11	IL-12 p40	IL-12 p70	IL-13
5	IL-15	IL-16	IL-17	IP-10	MCP-1	MCP-2	M-CSF	MG	MP-1α	MP-1β	MP-1δ	RANTES	TGF-β1
6	IL-15	IL-16	IL-17	IP-10	MCP-1	MCP-2	M-CSF	MG	MP-1α	MP-1β	MP-1δ	RANTES	TGF-β1
7	TNF-α	TNF-β	s TNF RI	s TNF RII	PDGF-BB	TIMP-2	NEG	NEG	IC 1	IC 2	IC 3	NEG	NEG
8	TNF-α	TNF-β	s TNF RI	s TNF RII	PDGF-BB	TIMP-2	NEG	NEG	IC 1	IC 2	IC 3	NEG	NEG

RayBio® Human Growth Factor Antibody Array 1

Detect 41 Growth Factors in one experiment

	a	b	c	d	e	f	g	h	i	j	k	l	m
1	Pos 1	Pos 2	Pos 3	Neg	Neg	AR	bFGF	b-NGF	EGF	EGF R	FGF-4	FGF-6	FGF-7
2	Pos 1	Pos 2	Pos 3	Neg	Neg	AR	bFGF	b-NGF	EGF	EGF R	FGF-4	FGF-6	FGF-7
3	GCSF	GDNF	GM-CSF	HB-EGF	HGF	IGFBP-1	IGFBP-2	IGFBP-3	IGFBP-4	IGFBP-6	IGF-1	IGF-1 SR	IGF-II
4	GCSF	GDNF	GM-CSF	HB-EGF	HGF	IGFBP-1	IGFBP-2	IGFBP-3	IGFBP-4	IGFBP-6	IGF-1	IGF-1 SR	IGF-II
5	M-CSF	M-CSF R	NT-3	NT-4	PDGF R a	PDGF R b	PDGF-AA	PDGF-AB	PDGF-BB	PIGF	SCF	SCF R	TGF-a
6	M-CSF	M-CSF R	NT-3	NT-4	PDGF R a	PDGF R b	PDGF-AA	PDGF-AB	PDGF-BB	PIGF	SCF	SCF R	TGF-a
7	TGF-b	TGF-b 2	TGF-b 3	VEGF	VEGF R2	VEGF R3	VEGF-D	neg	IC 1	IC 2	IC 3	Neg	Neg
8	TGF-b	TGF-b 2	TGF-b 3	VEGF	VEGF R2	VEGF R3	VEGF-D	neg	IC 1	IC 2	IC 3	Neg	Neg

We also offer Custom Human Cytokine Antibody Arrays. You can select the cytokines of interest from the following list and we will produce the customized array at an affordable price. For more information, please visit our website, www.raybiotech.com.

Note: IL-12 reacts both IL-12p40 and IL-12p70. IL-12p70 only recognizes IL-12p70.

Abbreviations: IP-10, Interferon-inducible protein-10; LAP, latency associated peptide (TGF- β 1); LIF, leukocyte inhibitory factor. MMP, Matrix Metalloproteinase; Pos, positive control; Neg, negative control. All other are used standard abbreviations.

4-1BB/TNFRSF9	adiponectin/Acrp30	AgRP(ART)	ALCAM	ANGIOGENIN	Angiopoietin-1
Angiopoietin-2	Angiostatin	AR (amphiregulin)	Axl	B7-1(CD80)	BDNF
bFGF	BLC	BMP-4	BMP-6	BMP-7	b-NGF
BTC	Cardiotrophin-1	CCL28/VIC	CD27	CD30	CD40
CD40 Ligand	Ck beta 8-1	CNTF	CTACK/CCL27	CTLA-4	CXCL16
Dkk-4	DR6	Dtk	EGF	EGF R	ENA-78
Endostatin	Eotaxin	Eotaxin-2	Eotaxin-3	E-Selectin	Fas/TNFRSF6
FGF-4	FGF-6	FGF-7	FGF-9	Flt-3 Ligand	Follistatin
Fractalkine	GCP-2	GCSF	GNDF	GTR Ligand/TNFSF18	GTR/TNFRF18
GM-CSF	GRO	HB-EGF	HCC-4/CCL16	HGF	HVEM
I-309	ICAM-1	ICAM-2	ICAM-3	IFN-gamma	IGFBP-1
IGFBP-2	IGFBP-3	IGFBP-4	IGFBP-6	IGF-I	IGF-I SR
IGF-II	IL-1 R4/ST2	IL-1 sRI	IL-1 sRII	IL-1 alpha	IL-1 beta
IL-1ra	IL-2	IL-2 R alpha	IL-2 R beta	IL-2 R gamma	IL-2 sR alpha
IL-3	IL-4	IL-5	IL-5 R alpha	IL-6	IL-6 sR
IL-7	IL-8	IL-9	IL-9 R	IL-10	IL-10 R alpha
IL-10 R beta	IL-11	IL-12	IL-12 p40	IL-12 p70	IL-13
IL-13 R alpha	IL-15	IL-16	IL-17	IL-18 BP alpha	IL-18 R alpha
IL-18 R beta	IL-21 R	IP-10	I-TAC/CXCL11	LAP(TGF-b1)	LEPTIN(OB)
LIF	LIGHT	L-Selectin	Lymphotactin	MCP-1	MCP-2
MCP-3	MCP-4	M-CSF	M-CSF R	MDC	MF
MIG	MIP-1 alpha	MIP-1 beta	MIP-1 gamma	MIP-3 alpha	MIP-3 beta
MMP-1	MMP-2	MMP-3	MMP-9	MMP-10	MMP-13
MPIF-1	MSP alpha	NAP-2	NGF R	NT-3	NT-4
ONCOSTATIN M	Osteoprotegerin	PARC	PDGF R alpha	PDGF R beta	PDGF-AA
PDGF-BB	PECAM-1	PF4	PIGF	Prolactin	P-selectin
RANTES	SCF	SCF R	SDF-1	sgp130	ST2
sTNF RII/TNFRS1B	sTNT R/TNFRS1A	Tarc	TECK/CCL25	TGF-alpha	TGF-beta
TGF-beta 2	TGF-beta 3	Tie-1	Tie-2	TIMP-1	TIMP-2
TIMP-3	TIMP-4	TNF-alpha	TNF-beta	TPO	TRAIL R1
TRAIL R2	TRAIL s R3/TNFRS1C	TRAIL s R4/TNFRS1D	u PAR	VCAM-1	VE-Cadherin
VEGF	VEGF-D	VEGF R2	VEGF R3	VEGI	

RayBiotech, Inc., the protein array pioneer company, strives to research and develop new products to meet demands of the biomedical community. RayBio's patent-pending technology allows detection of over 180 cytokines, chemokines and other proteins in a single experiment. Our format is simple, sensitive, reliable and cost effective. Products include: Cytokine Arrays, Chemokine Arrays, ELISA kits, Phosphotyrosine kits, Recombinant Proteins, Antibodies, and custom services.

1. Antibody arrays
 - Cytokine antibody array
 - Human cytokine antibody arrays
 - Mouse cytokine antibody arrays

Rat cytokine antibody arrays
Pathway- or disease-focused antibody arrays
Inflammation antibody array
Angiogenesis antibody array
Chemokine antibody array
Growth factor antibody array
MMP antibody array
Atherosclerosis antibody array

Antibody analysis tool, software

2. ELISA
3. Cell-based phosphorylation assay
4. Custom antibody arrays
5. Antibody
6. Recombinant protein
7. Cytokine protein arrays

RayBiotech also provides excellent custom service:

1. Antibody arrays
2. Protein arrays
3. Peptide synthesis
4. Production of recombinant protein and antibody
5. Peptide arrays
6. Phosphorylation arrays
7. ELISA

Just simply send your samples and we will do the assay for you.

Technology transfer program

Have you developed technologies or reagents interested to the scientific and research community? RayBiotech can help you commercialize your technologies, reagents and dream.

VI. Troubleshooting guide

Problem	Cause	Recommendation
Weak signal	Inadequate detection	Check laser power and PMT parameters
	Inadequate reagent volumes or improper dilution	Check pipetters 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	Excess of biotinylated antibodies	Make sure correct amount of antibodies
	Excess of streptavidin	Make sure correct amount of streptavidin
	Inadequate detection	Check laser power And PMT parameters
	dust	Work in clean environment
	Insufficient wash	Increase wash time and use more wash buffer
Uneven signal	Bubbles formed during incubation	Avoid bubble formation during incubation
	Arrays are not completed Covered by reagent	Completely cover arrays with solution

Reference List

1. LPS induces the interaction of a transcription factor, LPS-induced TNF- α factor, and STAT6(B) with effects on multiple cytokines. Xiaoren Tang, Deborah Levy Marciano, Susan E. Leeman and Salomon Amar. **PNAS**. April 5, 2005 vol. 102 no. 14 5132-5137
2. HIV-1-mediated apoptosis of neuronal cells: Proximal molecular mechanisms of HIV-1-induced encephalopathy. Yan Xu, Joseph Kulkoshy, Roger j. Pomerantz. **PNAS**. 2004 May 4, 2004 Vol. 101 No. 18.
3. Synergistic increases in intracellular Ca²⁺, and the release of MCP-1, RANTES, and IL-6 by astrocytes treated with opiates and HIV-1 Tat. **GLIA**. 2005 Apr 15;50(2):91-106.
4. Bone Marrow Stroma Influences Transforming Growth Factor- β Production in Breast Cancer Cells to Regulate c-myc Activation of the Preprotachykinin-I Gene in Breast Cancer Cells. Hyun S. Oh, Anabella Moharita... Pranela Rameshwar. **CANCER RESEARCH**. 64, 6327-6336, September 1, 2004
5. Recombinant Herpes Simplex Virus Type 1 (HSV-1) Codelivering Interleukin-12p35 as a Molecular Adjuvant Enhances the Protective Immune Response against Ocular HSV-1 Challenge **JOURNAL OF VIROLOGY**. Mar. 2005 Vol. 79, No. 6.
6. Dysregulated Inflammatory Response to *Candida albicans* in a C5-Deficient Mouse Strain. Alaka Mullick, Miria Elias, Serge Picard, Philippe Gros. **Infection and Immunity**, Oct. 2004, p. 5868-5876, DOI: 10.1128/IAI.72.10.5868-5876.2004
7. Leukotriene B₄ Strongly Increases Monocyte Chemoattractant Protein-1 in Human Monocytes Li Huang, Annie Zhao, Frederick Wong, Julia M. Ayala, Jisong Cui **Arteriosclerosis, Thrombosis, and Vascular Biology**. 2004;24:1783-1788
8. Human CD1d-unrestricted NKT cells release chemokines upon Fas engagement. Martin Giroux and François Denis. Yan Xu, Joseph Kulkoshy, Roger j. Pomerantz. **Blood**. prepublished online September 2, 2004; DOI 10.1182/blood-2004-04-1537

9. Monitoring the response of orthotopic bladder tumors to granulocyte macrophage colony-stimulating factor therapy using the prostate-specific antigen gene as a reporter. Wu Q, Esuvaranathan K, Mahendran R. **Clinical Cancer Research**. 2004 Oct 15; 10(20):6977-84.
10. Neuroglial activation and neuroinflammation in the brain of patients with autism (p NA). Diana L. Vargas, Caterina Nascimbene, Chitra Krishnan, Andrew W. Zimmerman, Carlos A. Pardo. **Annals of Neurology**. 2005 Jan 1; DOI: 10.1002/ana.20315
11. Cytokine profiling of macrophages exposed to *Porphyromonas gingivalis*, its LPS or its FimA. Qingde Zhou, Tesfahun Desta, Dana T. Graves and Salomon Amar. **Infection and Immunity (IAI)**. 2005 Feb;73(2):935-43.
12. Veto-like activity of mesenchymal stem cells: functional discrimination between cellular responses to alloantigens and recall antigens. Rameshwar P. **Journal of Immunology**. 2003 Oct 1;171(7):3426-34.
13. Cytokine responses elicited in endothelial cells after treatment with a specific toxin. Jaya Pandey. **BioCompare Product Review**. May 13, 2004
14. Proteomic Characterization of the Interstitial Fluid Perfusing the Breast Tumor Microenvironment. A Novel Resource for Biomarker and Therapeutic Target Discovery. Julio E. Celis, Pavel Gromov, Teresa Cabezón, José M. A. Moreira, Noona Ambartsumian, Kerstin Sandelin, Fritz Rank, and Irina Gromova. **Molecular Cellular Proteomics**. April 2004; 11(3):328-39.
15. Increased Expression and Secretion of Interleukin-6 in Patients with Barrett's Esophagus.. Katerina Dvorakova, Harinder Garewal **Clinical Cancer Research**. 2004 Mar 15;10(6):2020-8.
16. Antibody array-generated profiles of cytokine release from THP-1 leukemic monocytes exposed to different amphotericin B formulations. Turtinen LW, Prall DN, Bremer LA, Nauss RE, Hartsel SC. **Antimicrobial Agents Chemotherapy**. 2004 Feb;48(2):396-403.
17. The promise of cytokine antibody arrays in drug discovery process. R.-P. Huang, W. Yang, D. Yang, L. Flowers, I. R. Horowitz, X. Cao and R. Huang. Expert opinion on drug discovery. (2005) 9:601-615.

Note:

RayBio[®] is the trademark of RayBiotech, Inc.

Cytokine protein arrays are RayBiotech patent-pending technology.

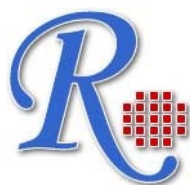
This product is intended for research only and is not to be used for clinical diagnosis. Our products may not be resold, modified for resale, or used to manufacture commercial products without written approval by RayBiotech, Inc.

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

Products are guaranteed for three 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.

ECL[™] is the trademark of Amersham Pharmacia Biotech.

This product is for research use only.



©2004 RayBiotech, Inc.