

RayBio[®] Mouse Cytokine Antibody Array C Series 1000

Patent Pending Technology

User Manual (Revised June 14, 2009)

RayBio[®] Mouse Cytokine Antibody Array C series 1000
(Combination of Arrays 3 & 4 Cat# AAM-CYT-1000)

RayBio[®] Mouse Cytokine Antibody Array 3 (Cat# AAM-CYT-3)

RayBio[®] Mouse Cytokine Antibody Array 4 (Cat# AAM-CYT-4)

RayBio[®] Mouse Cytokine Antibody Custom Array (Cat# AAM-CUST)

RayBio[®] Mouse Cytokine Antibody Array Service (Cat# AAM-SERV)

Please read manual carefully before starting experiment



RayBiotech, Inc.

**A Premium Provider of
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[®] Mouse Cytokine Antibody Array 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 Array Membrane.....	6
	C. Incubation.....	6
IV.	Protocol.....	7
	A. Blocking and Incubation.....	7
	B. Detection.....	9
V.	Interpretation of Results.....	10
VI.	Troubleshooting Guide.....	14
VII.	Reference List.....	15

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 a multitude of genes and signaling pathways. New techniques such as cDNA microarrays have enabled us to analyze global gene expression¹⁻³. However, almost all cell functions are executed by proteins, which cannot be studied simply through DNA and RNA techniques. Experimental analysis clearly shows a disparity between the relative expression levels of mRNA and their corresponding proteins⁴. Therefore, analysis of the protein profile is critical. 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 their broad application. Thus, effective study of multiple protein expression levels has been complicated, costly and time-consuming until now.

Our RayBio[®] Mouse Cytokine Antibody Array is the first commercially available cytokine 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 provides a simple format and highly sensitive approach to simultaneously detect multiple cytokine expression levels from conditioned media, patient's sera, cell lysate, tissue lysates and other sources.

The RayBio[®] Mouse Cytokine Antibody Array C series 1000 can detect 96 mouse cytokines in single experiment. RayBiotech also provides RayBio[®] Human Cytokine Antibody Array C series 4000 which is the only product available in the market that can detect 274 human cytokines in single experiment.

Traditionally, cytokines are detected by using ELISA (enzyme-linked immunosorbent assays); however, RayBiotech's approach has several

advantages over ELISA. First, and most important, our approach can simultaneously detect many cytokines. Secondly, the sensitivity is higher. With this approach, most cytokines can be detected at pg/ml levels. As little as 10 pg/ml of human IL-2 can be detected in the protein array format. Furthermore, the detection range is much greater than ELISA. For example, the detection range of human IL-2 varies from 10 to 100,000 pg/ml, whereas the detection range varies only within 100-1000 fold in a typical ELISA. Therefore, the detection range with protein arrays is greater than ELISA. Additionally, variability is far lower in comparison ELISA. As determined by densitometry, the variation between two spots ranged from 0 to 10% in duplicated experiments. In contrast, variation (about 20%) in ELISA is much higher. Finally, the system is much quicker and much easier to adapt to high-throughput techniques.

Pathway-specific array systems allow investigators to focus on the specific problem and are becoming an increasingly powerful tool in cDNA microarray systems. RayBiotech's first protein array system, known as RayBio[®] Mouse Cytokine Antibody Array, is particularly useful in comparison with the mouse 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. Cytokines play an important role in innate immunity, apoptosis, angiogenesis, cell growth and differentiation¹². They 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.

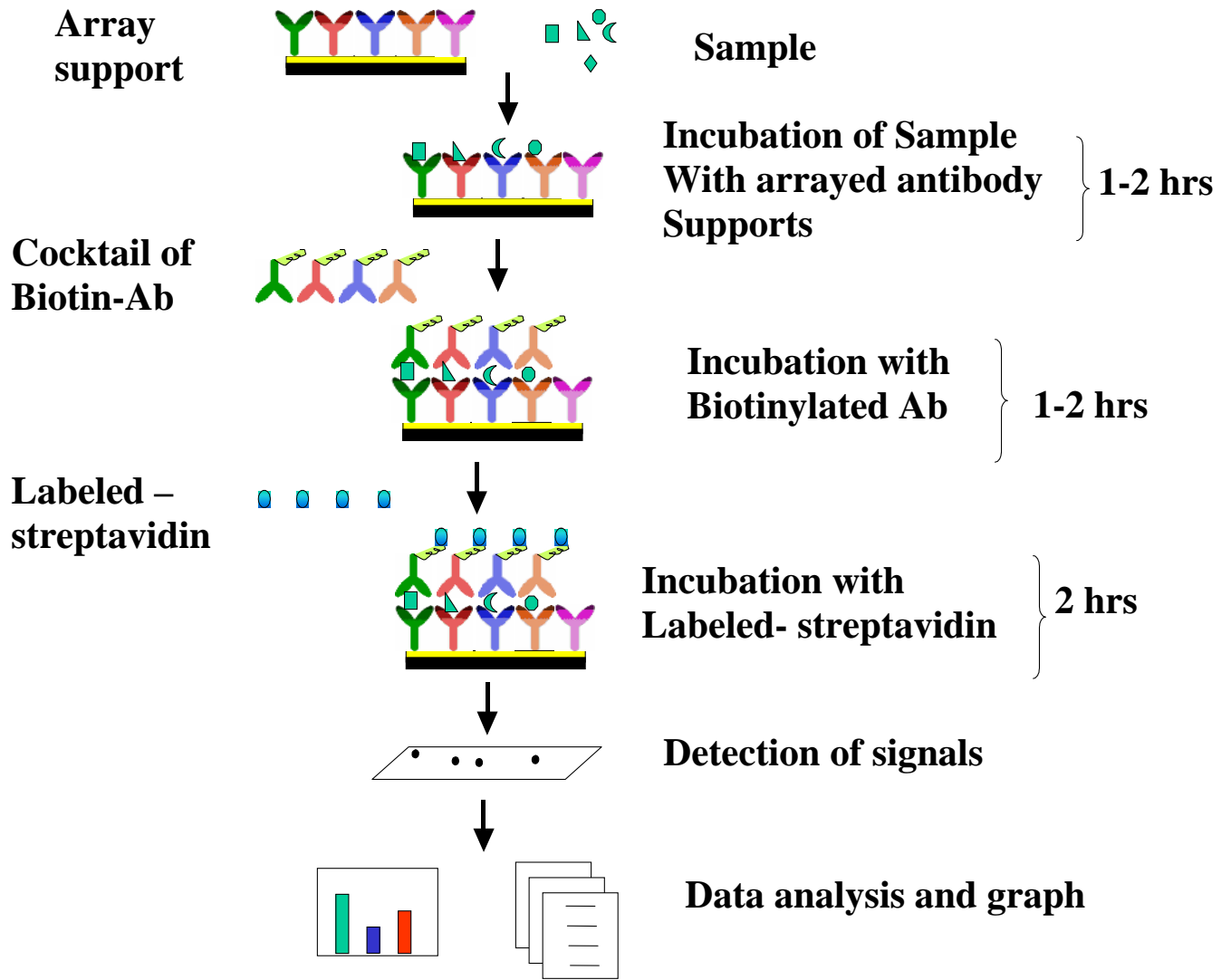
Without doubt, simultaneous detection of multiple cytokines provides a powerful tool to study cytokines.

1. LPS induces the interaction of a transcription factor, LPS-induced TNF- α factor, and STAT6(B) with effects on multiple cytokines. Tang X, Marciano DL, Leeman SE, Amar S. **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. Xu Y, Kulkoshy J, Pomerantz RJ. **PNAS**.

2004 May 4, 2004 Vol. 101 No. 18.

3. Synergistic increases in intracellular Ca(2+), 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. Oh HS, Moharita A, Rameshwar P. **Cancer Res**. 64, 6327-6336.
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 **J Virol**. 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. **Infect Immunity**, Oct. 2004, p. 5868-5876.
7. Leukotriene B₄ Strongly Increases Monocyte Chemoattractant Protein-1 in Human Monocytes Li Huang, Annie Zhao, Frederick Wong, Julia M. Ayala, Jisong Cui **Arterioscler Thromb Vascul Biol**. 2004;24:1783-1788
8. Human CD1d-unrestricted NKT cells release chemokines upon Fas engagement. Giroux M 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. **Clin Cancer Res**. 2004 Oct 15; 10(20):6977-84.
10. Neuroglial activation and neuroinflammation in the brain of patients with autism (p NA). Vargas DL, Nascimbene C, Krishnan C, Zimmerman AW, Pardo CA. **Ann Neurology**. 2005 Jan 1; DOI: 10.1002/ana.20315
11. Cytokine profiling of macrophages exposed to *Porphyromonas gingivalis*, its LPS or its FimA. Zhou Q, Desta T, Graves DT, Amar S. **Infect Immunity** (IAI). 2005 Feb;73(2):935-43.

Here's how it works



II. Materials Provided

Upon receipt, all components of the RayBio[®] Mouse Cytokine Antibody Array kit should be stored at -20°C to -80°C. At -20°C to -80°C the kit will retain complete activity for up to 6 months. Once thawed, the array membranes and 1X Blocking Buffer should be kept at -20°C and all other components should be stored at 4°C. After thawing the reagents, the kit must be used within three months, and please use the kit within six months of purchase.

- RayBio[®] Mouse Cytokine Antibody Array membranes (2/4/8 array 3 membranes, 2/4/8 array 4 membranes)
- Biotin-Conjugated Anti-Cytokines (2/4/8 tubes, each tube for two membranes)
- 1,000X HRP-Conjugated Streptavidin (50 µl)
- 1X Blocking Buffer (25ml/50ml)
- 20X Wash Buffer I (10/20ml)
- 20X Wash Buffer II (10/20ml)
- 2X Cell Lysis Buffer (10/20ml)
- Detection Buffer C (1.5/2.5ml)
- Detection Buffer D (1.5/2.5ml)
- Eight-Well Tray (1 each)
- Manual

Additional Materials Required

- Small plastic boxes or containers
- Orbital shaker
- Plastic sheet protector or Saran Wrap
- Kodak X-Omat AR film (REF 165 1454) and film processor or Chemiluminescence imaging system

III. Overview and General Considerations

A. Preparation of Samples

- Use serum-free conditioned media if possible.
- If serum-containing media is required, use an uncultured media aliquot as a negative control sample, since many types of sera contain cytokines.
- For cell lysates and tissue lysates, we recommend using RayBio® Cell Lysis Buffer to extract proteins from cell or tissue (e.g. using homogenizer). Dilute 2X RayBio® Cell Lysis Buffer with H₂O (we recommend adding proteinase inhibitors to Cell Lysis Buffer before use). After extraction, spin the sample down and save the supernatant for your experiment. Determine protein concentration.
- We recommend using per membrane:
 - 1 ml of Conditioned media (undiluted), or
 - 1 ml of 2-fold to 5-fold diluted sera or plasma, or
 - 50-500 µg of total protein for cell lysates and tissue lysates (use ~200-250 µg of total protein for first experiment) ***Dilute the lysate at least 10 fold with 1 X blocking buffer.***

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

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

B. Handling Array Membranes

- Always use forceps to handle membranes, and grip the membranes by the edges only.
- Never allow the array membranes to dry during experiments.

C. Incubation

- Completely cover the membranes with sample or buffer during incubation, and cover the eight-well tray with a lid to avoid drying.

- Avoid foaming during incubation steps.
- Perform all incubation and wash steps under gentle rotation.
- Several incubation steps such as step 2 (blocking), step 3 (sample incubation), step 7 (biotin-Ab incubation) or step 10 (HRP-streptavidin incubation) may be done at 4°C for overnight.

IV. Protocol

A. Blocking and Incubation

1. Place array membrane 3 (top left corner marked with “3”) and array membrane 4 (top left corner marked with “4”) into same well of the provided eight-well tray (“3” or “4” marked side is the antibody printed side).
2. Add 2 ml 1X Blocking Buffer and incubate at room temperature for 30 min to block membranes. Add some Blocking Buffer between the two membranes. Make sure there are no bubbles between membranes.
3. Decant 1X Blocking Buffer from each container. Incubate membranes with 1.2 ml of sample at room temperature for 1 to 2 hours. Dilute sample using 1X Blocking Buffer if necessary.

*Note: We recommend using 1.2 ml of undiluted conditioned media or 1.2 ml of 2-fold to 5-fold diluted sera or plasma or ~200-250 ug (range: 50-500 ug) of total protein for cell lysates and tissue lysates. **Dilute the lysate at least 10 fold with 1X blocking buffer. Add some samples between array membrane 3 and 4. Make sure there are no bubbles between membranes.***

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

4. Decant the samples from each container, and wash 3 times with 2 ml of 1X Wash Buffer I at room temperature with shaking. Please allow 5 min per wash. Dilute 20X Wash Buffer I with H₂O.

5. Wash 2 times with 2 ml of 1X Wash Buffer II at room temperature with shaking. Allow 5 min per wash. Dilute 20X Wash Buffer II with H₂O.
6. **From this step, place array membrane 3 into one well and array membrane 4 into another well.**
7. Prepare working solution for primary antibody.

Add 100 µl of 1X blocking buffer to the Biotin-Conjugated Anti-Cytokines tube. Mix gently and transfer all mixture to a tube containing 2 ml of 1X blocking buffer.

Note: the diluted biotin-conjugated antibodies can be stored at 4°C for 2-3 days.

8. Add 1 ml of diluted biotin-conjugated antibodies to each membrane. Incubate at room temperature for 1-2 hours.

Note: incubation may be done at 4°C for overnight.

9. Wash as directed in steps 4 and 5.
10. Add 2 ml of **1,000** fold diluted HRP-conjugated streptavidin (e.g. add 2 µl of HRP-conjugated streptavidin to **1998** µl 1X Blocking Buffer) to each membrane.

Note: mix the tube containing 1,000X HRP-Conjugated Streptavidin well before use since precipitation may form during storage.

11. Incubate at room temperature for 2 hours.

Note: incubation may be done at 4°C for overnight.

12. Wash as directed in steps 4 and 5.

B. Detection

*** Do not let the membrane dry out during detection. The detection process must be completed within 40 minutes without stopping.**

1. Proceed with the detection reaction.

Add 250 μ l of 1X Detection Buffer *C* and 250 μ l of 1X Detection Buffer *D* for one membrane; mix both solutions. Drain off excess wash buffer by holding the membrane vertically with forceps. Place membrane protein side up (“3“ or “4” mark is on the protein side top left corner) on a clean plastic sheet (provided in the kit). Pipette the mixed Detection Buffer onto the membrane and incubate at room temperature for 2 minutes. Ensure that the detection mixture completely and evenly covers the membrane without any air bubbles.

2. Drain off any excess detection reagent by holding the membrane vertically with forceps and touching the edge against a tissue. Gently place the membrane, protein side up, on a piece of plastic sheet (“3“ or “4” mark is on the protein side top left corner). Cover the array with another piece of plastic sheet. Gently smooth out any air bubbles. Avoid exerting any pressure on the membrane.

3. Expose the array to x-ray film (we recommend to use Kodak X-Omat AR film) and detect the signal using film developer, or the signal can be detected directly from the membrane using a chemiluminescence imaging system.

Expose the membranes for 40 Seconds. Then re-expose the film according to the intensity of signals. If the signals are too strong (background too high), reduce exposure time (e.g. 5-30 seconds). If the signals are too weak, increase exposure time (e.g. 5-20 min or overnight). Or re-incubate membranes overnight with 1X HRP-conjugated streptavidin, and redo detection in the second day.

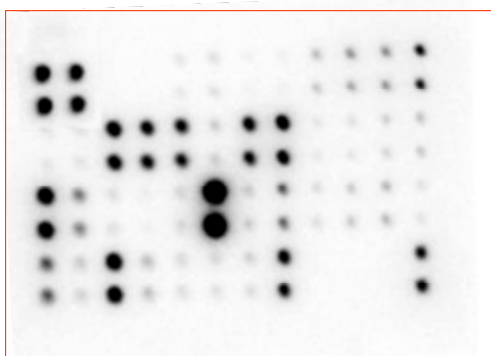
4. Save membranes in -20°C to -80°C for future reference.

V. Interpretation of Results:

The following figure shows RayBio[®] Mouse Cytokine Antibody Array membranes probed with conditioned media. Membranes were exposed to Kodak X-Omat film at room temperature for 1 minute. The biotin-conjugated IgG produces positive signals, which can be used to identify the orientation and compare relative expression levels among the different membranes.

One important parameter is background. To obtain the best results, we suggest that several exposures be attempted. We also strongly recommend using a negative control in which the sample is replaced with an appropriate mock buffer according to the array protocol, particularly during your first experiment.

Typical results using RayBio[®] Cytokine Antibody arrays



By comparing the signal intensities, relative expression levels of cytokines can be made. The intensities of signals can be quantified by densitometry. The positive control can be used to normalize the results from the different membranes being compared. The signals also can be detected and quantified by using a chemiluminescence-imaging device.

The **RayBio[®] Analysis Tool** is a program specifically designed for analysis of RayBio[®] Antibody Arrays. This tool will not only assist in compiling and organizing your data, but reduces your calculations to a “copy and paste.” Call RayBiotech, Inc. at 770-729-2992 for ordering information.

RayBio[®] Mouse Cytokine Antibody Array 3 (62)

	A	B	C	D	E	F	G	H	I	J	K	L	M	N
1	POS	POS	NEG	NEG	Blank	Axl	BLC	CD30 L	CD30 T	CD40	CRG-2	CTACK	CXCL16	Eotaxin
2	POS	POS	NEG	NEG	Blank	Axl	BLC	CD30 L	CD30 T	CD40	CRG-2	CTACK	CXCL16	Eotaxin
3	Eotaxin-2	Fas Ligand	Fractalkine	GCSF	GM-CSF	IFN γ	IGFBP-3	IGFBP-5	IGFBP-6	IL-1 α	IL-1 beta	IL-2	IL-3	IL-3 Rb
4	Eotaxin-2	Fas Ligand	Fractalkine	GCSF	GM-CSF	IFN γ	IGFBP-3	IGFBP-5	IGFBP-6	IL-1 α	IL-1 beta	IL-2	IL-3	IL-3 Rb
5	IL-4	IL-5	IL-6	IL-9	IL-10	IL-12 p40/p70	IL-12 p70	IL-13	IL-17	KC	Leptin R	Leptin	LIX	L-Selectin
6	IL-4	IL-5	IL-6	IL-9	IL-10	IL-12 p40/p70	IL-12 p70	IL-13	IL-17	KC	Leptin R	Leptin	LIX	L-Selectin
7	Lymphotactin	MCP1	MCP-5	M-CSF	MIG	MIP-1 α	MIP-1 γ	MIP-2	MIP-3 β	MIP-3 α	PF-4	P-Selectin	RANTES	SCF
8	Lymphotactin	MCP1	MCP-5	M-CSF	MIG	MIP-1 α	MIP-1 γ	MIP-2	MIP-3 β	MIP-3 α	PF-4	P-Selectin	RANTES	SCF
9	SDF-1 α	TARC	TCA-3	TECK	TIMP-1	TNF α	sTNF RI	sTNF RII	TPO	VCAM-1	VEGF	Blank	Blank	POS
#	SDF-1 α	TARC	TCA-3	TECK	TIMP-1	TNF α	sTNF RI	sTNF RII	TPO	VCAM-1	VEGF	Blank	Blank	POS

RayBio[®] Mouse Cytokine Antibody Array 4 (34)

	A	B	C	D	E	F	G	H	I	J	K	L
1	POS	POS	NEG	NEG	BLANK	bFGF	DPPIV/CD26	Dtk	E-Selectin	Fc γ RIIB	It-3 Ligan	GITR
2	POS	POS	NEG	NEG	BLANK	bFGF	DPPIV/CD26	Dtk	E-Selectin	Fc γ RIIB	It-3 Ligan	GITR
3	HGF R	ICAM-1	IGFBP-2	IGF-I	IGF-II	IL-15	IL-17B R	IL-7	I-TAC	Lungkine	MDC	MMP-2
4	HGF R	ICAM-1	IGFBP-2	IGF-I	IGF-II	IL-15	IL-17B R	IL-7	I-TAC	Lungkine	MDC	MMP-2
5	MMP-3	Osteopontin	Osteoporotegerin	Pro-MMP-9	Resistin	Shh-N	Thymus CK-1	TIMP-2	TRANCE	TROY	TSLP	VEGF R1
6	MMP-3	Osteopontin	Osteoporotegerin	Pro-MMP-9	Resistin	Shh-N	Thymus CK-1	TIMP-2	TRANCE	TROY	TSLP	VEGF R1
7	VEGF R2	VEGF R3	VEGF-D	BLANK	BLANK	BLANK	BLANK	BLANK	BLANK	BLANK	BLANK	POS
8	VEGF R2	VEGF R3	VEGF-D	BLANK	BLANK	BLANK	BLANK	BLANK	BLANK	BLANK	BLANK	POS

Abbreviations: Pos: positive control; Neg: negative control. All others use standard abbreviations.

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

We also offer Custom Mouse 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.

RayBio® Mouse Custom Array Antibody List

Choose from 146 cytokines and other proteins

4-1BB	6Ckine	ACE	ALK-1	Amphiregulin	Axl
bFGF	BLC	Cardiotrophin-1	CD27	CD27 Ligand	CD30
CD30 Ligand	CD36	CD40	CD40 Ligand	Chordin	CRG-2
CTACK	CTLA-4	CXCL16	Decorin	DKK-1	DPPIV
Dtk	E-Cadherin	EGF	Endoglin	Eotaxin	Eotaxin-2
Epigen	Epiregulin	E-Selectin	FasLigand	Fc gamma RIIB	Flt-3 Ligand
Fractalkine	Galectin-1	Galectin-3	Galectin-7	GCSF	GITR
GITR Ligand	GM-CFS	Granzyme B	Growth arrest specific 1	Growth arrest specific 6	HAI-1
HGF	HGFR	ICAM-1	IFN gamma	IGFBP-2	IGFBP-3
IGFBP-5	IGFBP-6	IGF-I	IGF-II	IL-1 alpha	IL-1 beta
IL-1 R4/ST2L	IL-10	IL-11	IL-12 p40	IL-12 p70	IL-13
IL-15	IL-17	IL-17 B	IL-17 BR	IL-17 E	IL-17 F
IL-1ra	IL-2	IL-2 R alpha	IL-20	IL-21	IL-28/IFN-lambda
IL-3	IL-3 R beta	IL-4	IL-5	IL-6	IL-6 R
IL-7	IL-9	I-TAC	JAM-A	KC	Leptin R
LEPTIN(OB)	LIX	L-Selectin	Lungkine	Lymphotactin	MAdCAM-1
MCP1	MCP-5	M-CSF	MDC	MFG-E8	MIG
MIP-1alpha	MIP-1gamma	MIP-2	MIP-3 alpha	MIP-3 beta	MMP-2
MMP-3	Nepriylisin	Osteopontin	Osteoporotegerin	Pentraxin 3/TSG 14	PF-4
Prolactin	Pro-MMP-9	P-Selectin	RAGE	RANTES	Resistin
SCF	SDF-1alpha	Shh-N	sTNF RI	sTNF RII	TACI
TARC	TCA-3	TECK	Thymus CK-1	TIMP-1	TIMP-2
TNF alpha	TPO	TRANCE	TREM-1	TROY	TSLP
TWEAK R	TWEAK	VCAM-1	VEGF	VEGF-D	VEGF R1
VEGF R2	VEGF R3				

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 507 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

- Quantibody arrays for quantitative measurement of cytokine and other protein concentration

- Phosphorylation antibody arrays

- Biotin label-based antibody arrays for high density antibody arrays.

- Antibody analysis tool, software

2. ELISA

3. Cell-based phosphorylation assay

4. Custom antibody arrays

5. Antibody

6. Recombinant protein

7. 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 of interest to the scientific and research community? RayBiotech can help you commercialize your technologies, reagents and dream.

VI. Troubleshooting guide

Problem	Cause	Recommendation
Weak signal or no signal	1. Taking too much time for Detection.	1. The whole Detection process must be completed in 30 min.
	2. Film developer does not work properly.	2. Fix film developer.
	3. Did not mix HRP-streptavidin well before use.	3. Mix tube containing 1,000X HRP-Conjugated Streptavidin well before use since precipitation may form during storage.
	4. Sample is too dilute.	4. Increase sample volume, (e.g. using undilute sample) or using more cells (e.g. seed 2 million cells. After 1 or 2 days, change complete medium with low serum medium and collect conditioned medium 2 day later. Use about 1 to 2 ml of conditioned medium for experiment).
	5. Other.	1. Reduce blocking concentration by diluting in 1X Wash Buffer II. 2. Slightly increase HRP concentrations. 3. Slightly increase biotin-antibody concentrations. 4. Expose film for overnight to detect weak signal.
Uneven signal	1. Bubbles formed during incubation.	1. Remove bubble during incubation.
	2. Membranes were not completely covered by solution.	2. Completely cover membranes with solution.
High background	1. Exposure to x-ray file is too long.	1. Decrease exposure time.
	2. Membranes were allowed to dry out during experiment.	2. Completely cover membranes with solution during experiment.
	3. Sample is too concentrated.	3. Use more diluted sample.

Reference List

- 12.
13. Y. Lin, Ruochun Huang, Li-Pai Chen, Henry Lisoukov, Zhen-Hai Lu, Shiyong Li, Cheng C. Wang and R.-P. Huang. (2003) Profiling of cytokine expression by biotin-labeled-based protein arrays. **Proteomics**.3: 1750-1757.
14. C. C. Wang, R.-P. Huang, H. Lisoukov, M. Sommer, R. Huang, Y. Lin and J. Burke. (2002) Array-based multiplexed screening and quantitation of human cytokines and chemokines. **J. Proteome Res.** 1:337-343.
15. R. Huang, Y. Lin, C. C. Wang, J. Gano, B. Lin, Q. Shi, A. Boynton, J. Burke and R.-P. Huang. (2002) Connexin suppresses human glioblastoma cell growth by down-regulation of monocyte chemotactic protein 1, as discovered using protein array technology. **Cancer Research.** 62 : 2806-2812.

R.-P. Huang, R. Huang, Y. Fan and Y. Lin. (2001). A novel method for high-throughput protein profiling from conditioned media and patient's sera. Ana. Biochem. 294(1):55-62.

Selected References Using RayBiotech Arrays

1. Ju Z, Jiang H, Jaworski M, Rathinam C, et al. Telomere dysfunction induces environmental alterations limiting hematopoietic stem cell function and engraftment. **Nature Med.** 2007; 13(6):742-747.
2. Yang F-C, Ingram DA, Chen S, Hingtgen CM, et al. Neurofibromin-deficient Schwann cells secrete a potent migratory stimulus for Nf1+/- mast cells. **J Clin Invest.** 2003;112(12):1851-1861.
3. Kong X-N, Yan H-X, Chen L, Dong L-W, et al. LPS-induced down-regulation of signal regulatory protein alpha contributes to innate immune activation in macrophages. **J Exper Med.** 2007; 204(11):2719-2731.
4. Young LR, Borchers MT, Allen HL, Gibbons RS, McCormack FX. Lung-Restricted Macrophage Activation in the Pearl Mouse Model of Hermansky-Pudlak Syndrome. **J Immunol.** 2006;176:4361-4368.

5. Kumar P, Wu H, McBride JL, Jung K-E, et al. Transvascular delivery of small interfering RNA to the central nervous system. **Nature**. 2007; 448:39-43.

16. 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 from 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 manufacture, RayBiotech's liability to the buyer for any claim relating to products shall be limited to replacement or refund of the purchase price.

Kodak X-Omat[™] is the trademark of Eastman Kodak Company.

This product is for research use only.



©2007 RayBiotech, Inc.