

RayBio[®] Human Obesity Antibody Array 1

User Manual (Revised February 4, 2009)

RayBio[®] Human Obesity Antibody Array 1 (Cat# AAH-ADI-1)

RayBio[®] Custom Human Cytokine Antibody Array (Cat# AAH-CUST)

RayBio[®] Human Cytokine Antibody Array Service (Cat# AAH-SERV)

Please read manual carefully before starting experiment



We Provide You with Excellent
Protein Array System 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 Obesity Antibody Array Protocol

TABLE OF CONTENTS

I.	Introduction.....	2
	How It Works.....	6
II.	Materials Provided.....	7
	Additional Materials Required.....	7
III.	Overview and General Considerations.....	8
	A. Preparation of Samples.....	8
	B. Handling Array Membrane.....	8
	C. Incubation.....	9
IV.	Protocol.....	9
	A. Blocking and Incubation.....	9
	B. Detection.....	10
V.	Interpretation of Results.....	11
VI.	Troubleshooting Guide.....	17
VII.	Reference List	18

Cytokine protein 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[®] Human Cytokine Antibody Array is the first commercially available cytokine antibody array system⁷⁻¹¹. Applying similar principle, now we have developed Human Cytokine Antibody Arrays. By using the RayBiotech system, scientists can rapidly and accurately identify the expression profiles of multiple adipokines in several hours inexpensively.

The RayBiotech kit provides a simple array format, and highly sensitive approach to simultaneously detect multiple adipokine expression levels from conditioned media, patient's sera, cell lysate, tissue lysates and other sources.

Traditionally, adipokines are detected by using ELISA. However, RayBiotech's approach has several advantages over ELISA. Most importantly, our approach can detect many adipokines simultaneously with trace amount of sample.

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 antibody array system, known as RayBio[®] Human obesity antibody array, is particularly useful in comparison with the human adipokine cDNA microarray system. Besides the ability to detect protein expression, RayBiotech's system is a more accurate reflection of active adipokine levels because it only detects secreted adipokines, and no amplification step is needed. Furthermore, it is much simpler, faster, environmentally friendlier, and more sensitive.

Simultaneous detection of multiple adipokines undoubtedly provides a powerful tool to study obesity. The area of obesity research is getting hotter ever over the past years. One of the key driving factor is that adipose tissue is found no longer to be an inert energy storage organ, but is emerging as an active participant in regulating physiological and pathologic processes. Many soluble factors have been identified from the adipose tissue and are so called as adipocytokines or adipokines. But adipokines are also expressed in a number of other tissues and organs. Because all of these factors can act in an autocrine, paracrine and endocrine manner in the organisms, adipokines are thought to serve as mediators linking obesity, inflammation, immunity and other obesity related diseases¹²⁻¹⁵.

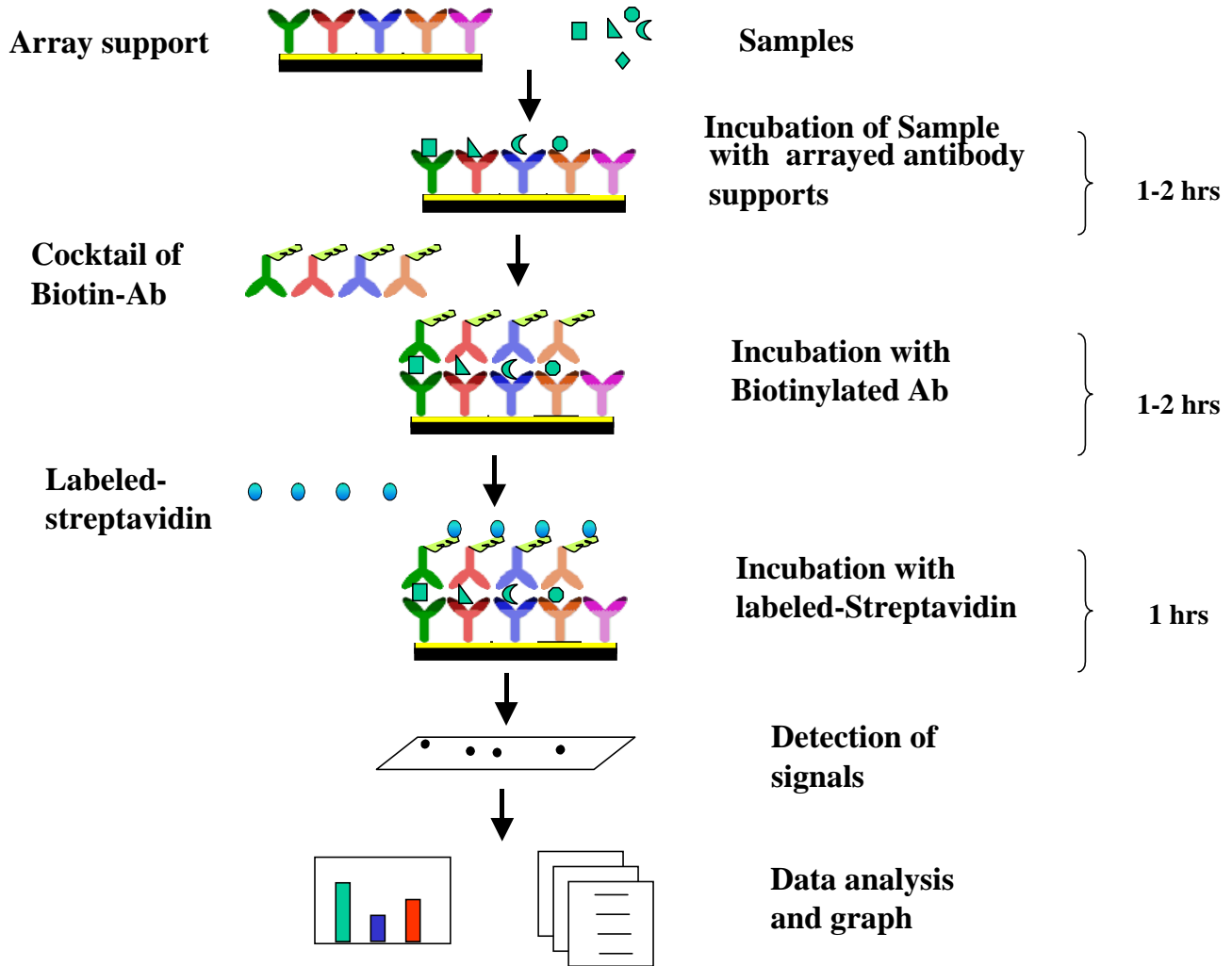
Without doubt, simultaneous detection of multiple adipokines provides a powerful tool to study adipokines.

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 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. 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 responses elicited in endothelial cells after treatment with a specific toxin. Jaya Pandey. **BioCompare Product Review**. May 13, 2004.
12. Mlinar B, et al: Molecular mechanisms of insulin resistance and associated diseases. **Clinica Chimica Acta**. 2007;375:20-35.

13. Tilg H, et al: Adipocytokines: mediators linking adipose tissue, inflammation and immunity. **Nat. Rev. Immunol.** 2006;6:772-783.
14. Trujillo ME, et al. Adipose tissue-derived factors: impact on health and disease. **Endocr Rev.** 2006;27:762-78 .
15. Van Gaal LF, et al. Mechanisms linking obesity with cardiovascular disease. **Nature.** 2006;444:875-880.
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.

Here's how it works



II. Materials Provided

Upon receipt, all components of the RayBio[®] Human Obesity 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 component 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[®] Human Obesity Antibody Array membranes (2/4/8 membranes)
- Biotin-Conjugated Anti-Adipokines (1/2/4 tubes, each tube for two membranes)
- 1,000X HRP-Conjugated Streptavidin (50 µl)
- 1X Blocking Buffer (25/50 ml)
- 20X Wash Buffer I (10/20 ml)
- 20X Wash Buffer II (10/20 ml)
- 2X Cell Lysis Buffer (10/20 ml)
- Detection Buffer C (1.5/2.5 ml)
- Detection Buffer D (1.5/2.5 ml)
- Eight-Well Tray (1 each)
- Manual

Additional Materials Required

- Small plastic boxes or containers
- Orbital shaker
- Plastic sheet protector or SaranWrap
- 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 adipokines. 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 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 8 (biotin-Ab incubation) and step 11 (HRP-streptavidin incubation) may be done at 4°C for overnight, but make sure to cover the 8 well plate tightly to prevent evaporation.

IV. Protocol

A. Blocking and Incubation

1. Place each membrane into the provided eight-well tray (“-“ means the antibody printed side).
2. Add 2 ml 1X Blocking Buffer and incubate at room temperature for 30 min to block membranes. Make sure there are no bubbles between the membranes.
3. Decant Blocking Buffer from each container, and incubate membranes with 1ml* of sample at room temperature for 1 to 2 hours. Dilute sample using 1X Blocking Buffer if necessary.

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. Prepare working solution for primary antibody.

Add 100 μ l of 1x blocking buffer to the Biotin-Conjugated Anti-Adipokines 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.

7. 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.

8. Wash as directed in steps 4 and 5.

9. Add 2 ml of **1,000** fold diluted HRP-conjugated streptavidin (e.g. add **2** μ l of HRP-conjugated streptavidin to **1998** μ l 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.

10. Incubate at room temperature for 2 hours.

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

11. 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 (“-“ mark is on the protein side top left corner) on a clean plastic sheet (provided in the kit). Pipette the mixed Detection Buffer on to the membrane and incubated at room temperature for 2 minute. Ensure that the detection mixture is 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 (“-“ mark is on the protein side top left corner). Cover another piece of plastic sheet on the array. Gently smooth out any air bubble. Avoid using 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.

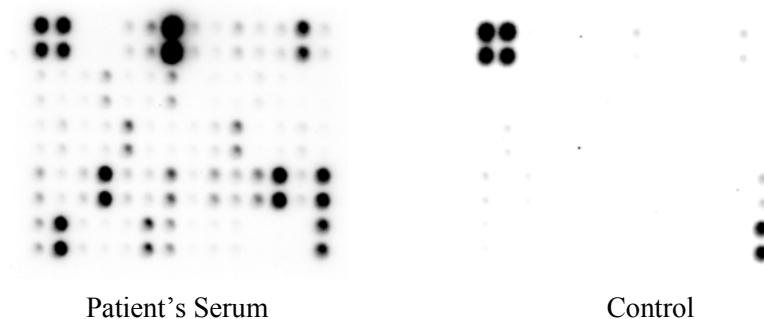
V. Interpretation of Results:

The following figure shows RayBio[®] Human Obesity Antibody Array membranes probed with conditioned media from two different cell lines. 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 to compare the 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[®] Obesity Antibody arrays

Human Adipokine Antibody Arrays 1



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

The **RayBio[®] Analysis Tool** is a program specifically designed for analysis of RayBio[®] Obesity Antibody Arrays. This tool will not only assist in compiling and organizing your data, but also reduces your calculations to

a “copy and paste.” Call RayBiotech, Inc. at 770-729-2992 for ordering information.

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 16 in introduction).

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

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

Human Adipokine Antibody Arrays 1
(for simultaneous detection of 62 human adipokine s)

	1	2	3	4	5	6	7	8	9	10	11	12	13	14
a	POS	POS	NEG	NEG	4-1BB	ACE-2	Acrp30	Adipsin	AgRP	Ang-2	Angiopoietin-1	ANGPTL4	CRP	ENA-78
b	POS	POS	NEG	NEG	4-1BB	ACE-2	Acrp30	Adipsin	AgRP	Ang-2	Angiopoietin-1	ANGPTL4	CRP	ENA-78
c	Fas	FGF-6	Growth Hormone	HCC-4	IFN-gamma	IGFBP-1	IGFBP-2	IGFBP-3	IGF-I	IGF-I SR	IL-1 R4/ST2	IL-1 sRI	IL-10	IL-11
d	Fas	FGF-6	Growth Hormone	HCC-4	IFN-gamma	IGFBP-1	IGFBP-2	IGFBP-3	IGF-I	IGF-I SR	IL-1 R4/ST2	IL-1 sRI	IL-10	IL-11
e	IL-12	IL-1alpha	IL-1beta	IL-6	IL-6 sR	IL-8	Insulin	IP-10	Leptin R	Leptin	LIF	Lymphotoctin	MCP-1	MCP-3
f	IL-12	IL-1alpha	IL-1beta	IL-6	IL-6 sR	IL-8	Insulin	IP-10	Leptin R	Leptin	LIF	Lymphotoctin	MCP-1	MCP-3
g	MCSF	MIF	MIP-1beta	MSP alpha	OPG	OSM	PAI-I	PARC	PDGF-AA	PDGF-AB	PDGF-BB	RANTES	Resistin	Serum Amyloid A
h	MCSF	MIF	MIP-1beta	MSP alpha	OPG	OSM	PAI-I	PARC	PDGF-AA	PDGF-AB	PDGF-BB	RANTES	Resistin	Serum Amyloid A
i	SDF-1	sTNF RII	sTNT RI	TECK	TGF-beta	TIMP-1	TIMP-2	TNF-alpha	VEGF	XEDAR	BLANK	BLANK	BLANK	POS
j	SDF-1	sTNF RII	sTNT RI	TECK	TGF-beta	TIMP-1	TIMP-2	TNF-alpha	VEGF	XEDAR	BLANK	BLANK	BLANK	POS

Abbreviations: OSM, oncostatin M; TPO, thrombopoietin; Pos, positive control; Neg, negative control. All others use standard abbreviations.

RayBiotech also offers customized cytokine protein 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,

Human Custom Antibody Array List (285 proteins)

4-1BB/TNFRSF9	CNTF	GDNF	IL-18 R alpha	MIP-1 alpha	SCF
ACE-2	Cripto-1	GITR	IL-18 R beta	MIP-1 beta	SCF R
Activin A	CRP	GITR Ligand	IL-1ra	MIP-1 delta	SDF-1 alpha
Adiponectin/Acrp30	CTACK/CCL27	GM-CSF	IL-2	MIP-3 alpha	SDF-1 beta
Adipsin/Factor D	CTLA-4	GRO	IL-2 R alpha	MIP-3 beta	sgp130
AFP	CXCL16	GRO-a	IL-2 R beta	MMP-1	Shh N
AgRP(ART)	DAN	Growth Hormom	IL-2 R gamma	MMP-2	Siglec-5
ALCAM	Decorin	HB-EGF	IL-21 R	MMP-3	Siglec-9
Angiogenin	DKK-1	HCC-4/CCL16	IL-22	MMP-7	sTNF RII
Angiopoietin-1	DKK-3	hCGa, intact	IL-28A/IFN-lambda	MMP-8	sTNT RI
Angiopoietin-2	DKK-4	HGF	IL29/IFN-lambda 1	MMP-9	TACE
Angiostatin	DPPIV/CD26	HVEM	IL-3	MMP-10	TARC
ANGPTL4	DR6	I-309	IL-31	MMP-13	TECK/CCL25
AR (amphiregulin)	Dtk	ICAM-1	IL-4	MPIF-1	TGF-alpha
Axl	E-Cadherin	ICAM-2	IL-5	MSP a Chain	TGF-beta 1
B7-1(CD80)	EDA-A2	ICAM-3	IL-5 R alpha	NAP-2	TGF-beta 2
Bate2 M	EGF	IFN-gamma	IL-6	NCAM-1	TGF-beta 3
BCAM	EGF R	IGFBP-1	IL-6 sR	NGF R	Thyroglobulin
BCMA/TNFRSF17	EG-VEGF/PK1	IGFBP-2	IL-7	Nidogen-1/Entactin	Tie-1
BDNF	ENA-78	IGFBP-3	IL-8	NrcAM	Tie-2
beta IG-H3	Endoglin	IGFBP-4	IL-9	NRG1-beta 1/HRG1-beta 1	TIM-1
Betacellulin (BTC)	Endostatin	IGFBP-5	IL-9 R	NT-3	TIMP-1
bFGF	Eotaxin	IGFBP-6	Insulin	NT-4	TIMP-2
BLC	Eotaxin-2	IGF-I	IP-10	Oncostatin M	TIMP-4
BMP-4	Eotaxin-3	IGF-I sR	I-TAC/CXCL11	Osteopontin	TNF-alpha
BMP-5	EpCAM/TROP1	IGF-II	LAP(TGF-b1)	Osteoprotegerin	TNF-beta
BMP-6	ErbB2	IL-1 alpha	Leptin R	PAI-1	TPO
BMP-7	ErbB3	IL-1 beta	LEPTIN(OB)	PARC	TRAIL R1
b-NGF	Erythropoietin R (EPO R)	IL-1 R4/ST2	LH	P-Cadherin	TRAIL R2
BTC	E-Selectin	IL-1 sRI	LIF	PDGF R alpha	TRAIL R3
CA125	Fas Ligand	IL-1 sRII	LIGHT	PDGF R beta	TRAIL R4
CA15-3	Fas/TNFRSF6	IL-10	LIMPII/SR-B2	PDGF-AA	Trappin-2/Etafin
CA19-9	Fcr RIIB/C	IL-10 R alpha	Lipocalin-2/NGAL	PDGF-AB	TREM-1
Carbonic Anhydrase IX(CA9)	Ferritin	IL-10 R beta	L-Selectin	PDGF-BB	TROY
Cardiotrophin-1 (CT-1)	FGF-4	IL-11	Lymphotactin	PECAM-1	TSH
Cathepsin S	FGF-6	IL-12 p40	LYVE-1	Platelet Factor 4	TSLP
CCL14a/HCC-1	FGF-7	IL-12 p70	Marapsin/Pancreasin	PIGF	u PAR
CCL21/6ckine	FGF-9	IL-13	MCP-1	Procalcitonin/Calcitonin	Ubiquitin+1
CCL28/VIC	FLRG	IL-13 Ra1	MCP-2	Prolactin	VCAM-1
CD14	Flt-3 Ligand	IL-13 Ra2	MCP-3	PSA-free	VE-Cadherin
CD23/Fc epsilon RII	Follistatin	IL-15	MCP-4	PSA-total	VEGF
CD27	Fractalkine	IL-16	MCSF	P-selectin	VEGF R2
CD30	FSH	IL-17	M-CSF R	RAGE	VEGF R3
CD40	Furin	IL-17B	MDC	RANK	VEGF-C
CD40 Ligand	Galectin-7	IL-17C	MICA	RANTES	VEGF-D
CEA	GCP-2	IL-17F	MICB	Resistin	
CEACAM-1	GCSF	IL-17R	MIF	S-100b	
CK beta 8-1	GDF-15/MIC-1	IL-18 BPa	MIG	SAA	

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 274 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, EIA kits, Recombinant Proteins, Antibodies, and custom services.

1. Antibody arrays
2. 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
 - Obesity antibody arrays
 - Antibody analysis tool, software
3. ELISA
4. Cell-based phosphorylation assay
5. Custom antibody arrays
6. Antibody
7. Recombinant protein
8. Cytokine protein arrays
9. Quantibody arrays for quantitative measurement of cytokine and other protein concentration.
10. Phosphorylation antibody arrays
11. Biotin label-based antibody arrays for high density antibody arrays
12. EIA
13. Peptide

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
8. EIA
9. Assay development

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 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.

VII. Selected References Using RayBiotech Antibody Arrays

1. Protein Array Reveals Differentially Expressed Proteins in Subcutaneous Adipose Tissue in Obesity. Škopková M, Penesová A, Sell H, Rádíková Z, et al. **Obesity**. 2007;15:2396-2406.
2. Adipokines oversecreted by omental adipose tissue in human obesity. Maury E, Ehala-Aleksejev K, Guiot Y, Detry R, Vandenhooft A, Brichard SM. **Am J Physiol Endocrinol Metab**. 2007; 293:E656-E665.
3. Atrial natriuretic peptide inhibits the production of adipokines and cytokines linked to inflammation and insulin resistance in human subcutaneous adipose tissue. Moro C, Klimeakova E, Lolmede K, Berlan M, et al. **Diabetologia**. 2007; 50:1038-1047.
4. Adipokine protein expression pattern in growth hormone deficiency predisposes to the increased fat cell size and the whole body metabolic derangements. Ukropec J, Penesova A, Skopkova M, Pura M, et al. **J Clin Endocrinol Metab**. 2008;93:2255-2262.
5. Cytokine secretion by human adipocytes is differentially regulated by adiponectin, AICAR, and troglitazone. Sell H, Dietze-Schroeder D, Eckardt K, Ecker J. **Biochem Biophys Res Comm**. 2006; 343(3):700-706.
6. Plasma, urine, and renal expression of adiponectin in human systemic lupus erythematosus. Rovin BH, Song H, Hebert LA, Nadasdy T, et al. **Kidney International**. 2005;68:1825-1833.
7. Adiponectin and functional adiponectin receptor 1 are expressed in airway epithelial cells in Chronic Obstructive Pulmonary Disease. Miller M, Cho JY, Pham A, Ramsdell J, Broide DH. **J Immunol**. 2009;182:684-691.
8. Sell H, Dietze-Schroeder D, Kaiser U, Eckel J. MCP-1 is a potential player in the negative crosstalk between adipose tissue and skeletal muscle. **Endocrinol**. 2005;147:(5):2458-2467.
9. Autocrine Action of Adiponectin on Human Fat Cells Prevents the Release of Insulin Resistance-Inducing Factors. Dietze-Schroeder D, Sell H, Uhlig M, Koenen M, Eckel J. **Diabetes**. 2005;54:2003-2011.

10. Leptin-mediated cytokine release and migration of eosinophils: Implications for immunopathophysiology of allergic inflammation. Wong CK, Cheung PFY, Lam CWK. **Eur J Immunol**. 2007;37:2337-2348.
11. Secreted proteins and genes in fetal and neonatal pig adipose tissue and stromal-vascular cells. Hausman GJ, Poulos SP, Richardson RL, Barb CR, et al. **J Animal Science**. 2006; 84:1666-1681.
12. Diet-induced adiposity alters the serum profile of inflammation in C57BL/6N mice as measured by antibody array. Fenton JI, Nuñez NO, Yakar S, Perkins SN, et al. **Diabetes Obesity Metab**. <http://dx.doi.org/10.1111/j.1463-1326.2008.00974.x>

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.

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

This product is for research use only.



©2009 *RayBiotech, Inc.*