G-Series Human Cytokine Antibody Array 640

A combination of 16 non-overlapping arrays to measure the relative expression levels of 640 human cytokines

Catalog #: GSH-CAA-640

User Manual Last revised October 1, 2021

Caution: Extraordinarily useful information enclosed



ISO 13485 Certified

3607 Parkway Lane, Suite 100 Norcross, GA 30092 Tel: 1-888-494-8555 (Toll Free) or 770-729-2992, Fax:770-206-2393 Web: www.RayBiotech.com, Email: info@raybiotech.com

Table of Contents

Section		Page #
l.	Overview	3
П.	Introduction	3
III.	How It Works	4
IV.	Materials Provided	5
V.	Storage	5
VI.	Additional Materials Required	5
VII.	General Considerations A. Sample Preparation B. Handling Glass Slides C. Incubation	6 6 6
VIII.	Protocol A. Completely Air Dry The Glass Slide B. Blocking & Incubation C. Incubation with Biotinylated Antibody Cocktail & Wash D. Incubation with Cy3 Equivalent Dye-Streptavidin & Wash E. Fluorescence Detection F. Data Analysis	7 7 7 8 8 9 10
IX.	Array Map	11
Χ.	Analysis Tool: Data Analysis Software	14
XI.	Troubleshooting Guide	15
XII.	Select Publications	16
XIII.	Experiment Record Form	17
XIV.	How To Choose An Array	18

Please read the entire manual carefully before starting your experiment

I. Overview

Assay Duration
Reproducibility
Sample Volume
Detection Method
Format
Cytokines Detected (640)

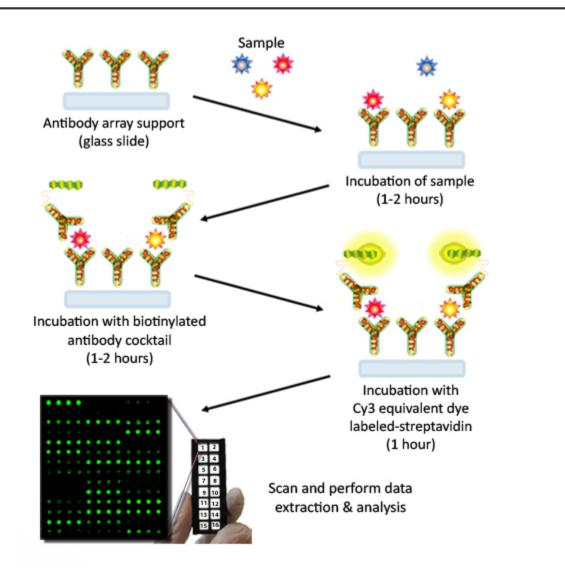
II. Introduction

Cytokines play an important role in innate immunity, apoptosis, angiogenesis, cell growth and differentiation. They are involved in interactions between different cell types, cellular responses to environmental conditions, and maintenance of homeostasis. In addition, cytokines are also involved in most disease processes, including cancer and cardiac diseases. RayBio[®] G-Series Arrays are glass slide-based antibody arrays which allow researchers to conduct rapid, accurate expression profiling of hundreds of cytokines, chemokines, growth factors, proteases, soluble receptors and other proteins from any biological fluid. Like a traditional sandwich-based ELISA, this array uses a matched pair of cytokine-specific antibodies for detection. After incubation with the sample, the target cytokines are captured by the antibodies printed on the solid surface. A second biotin-labeled detection antibody is then added, which recognizes a different epitope of the target cytokine. The cytokine-antibody-biotin complex can then be visualized through the addition of the streptavidin-conjugated Cy3 equivalent dye. Like the Quantibody[®] arrays, G-Series utilizes a highly sensitive and stable fluorescent readout which can be detected by most laser fluorescent scanner systems. After capturing the spot densities with a laser scanner,

normalization of the raw data can be easily calculated by the researcher, or by a quick copypaste into our excel-based Analysis Tool software.

This array as well as all catalog numbers beginning with 'GS' differ from the classic G-Series Arrays in a few important ways. First, each capture antibody is printed in quadruplicate instead of duplicate, delivering higher precision. Secondly, this array features the same antibody panels used in our Quantibody Arrays, allowing a seamless transition to our quantitative multiplex assay platform. Lastly, all 16 wells are spotted as sub-arrays, delivering easy handling of 16 samples simultaneously while consuming low sample volumes (10 - 100 μ l per array).

III. How It Works



IV. Materials Provided

	Catalog #	Component Name	1 Slide Box	2 Slide Box*
1	[Array-Cat-#]S	Array-specific Glass Slide	1	2
2	QA-SDB	Sample Diluent	15	ml
3	AA-WB1-30ML	20X Wash Buffer I	2 x 30 ml	3 x 30 ml
4	AA-WB2-30ML	20X Wash Buffer II	30	ml
5	[Array-Cat-#]B	<i>Array-specific</i> Biotinylated Antibody Cocktail	1-25 µl	2 x 1-25 µl
6	QA-CY3E	Cy3 equivalent dye-conjugated Streptavidin	5 μΙ	2 x 5 µl
7	QA-SWD	Slide Washer/Dryer	1 x 30 r	nl Tube
8	QA-ADH	Adhesive Film	1	2

This product is a combination of multiple arrays. Items 1, 5, & 6 are array-specific.

* 4 slide kits are comprised of 2 separate 2 slide kits.

V. Storage

Upon receipt, all components should be stored at -20°C. The kit will retain activity for up to 6 months. Once thawed, the glass slide, antibody cocktail and dye-conjugated Streptavidin should be kept at -20°C. All other components may be stored at 4°C. The entire kit should be used within 6 months of purchase.

VI. Additional Materials Required

- Benchtop rocker or orbital rocker
- Laser scanner for fluorescence detection
- Aluminum foil
- Distilled water
- 1.5 ml Polypropylene microcentrifuge tubes

A. Preparation of Samples

- Use serum-free conditioned media if possible.
- If serum-containing conditioned media is required, it is highly recommended that complete medium be used as a control since many types of sera contains cytokines.
- Each array needs 100 µl of total sample volume. To avoid matrix effects, we recommend using a minimum of 2x dilution for serum, plasma, cell culture media, or other body fluids, or 500 µg/ml-1 mg/ml (after a 5-fold to 10-fold dilution to minimize the effects of any detergent(s)) total protein for cell and tissue lysates. Please be aware, more sample volume is required for combination arrays. For example, the minimum sample volume for a 10-array kit is 500 µl, or 500 µg cell lysate.

If you experience high background or if the fluorescent signal intensities exceed the detection range, further dilution of your sample is recommended.

B. Handling Glass Slides

- Do not touch the surface of the slides, as the microarray slides are very sensitive. Hold the slides by the edges only.
- Handle all buffers and slides with powder free gloves.
- Handle glass slide/s in clean environment.
- Permanent marker ink can significantly interfere with fluorescent signal detection. To help distinguish one slide from another, you may make a small marking (such as a number or a star) along the top or bottom edge, using a green or blue ultra-fine point Sharpie[®] brand marker. This can also serve to orient the slide. For best results during scanning, please **DO NOT**:
 - Write anywhere on the front (arrayed) side of the slide
 - $\circ\,$ Write on the slide while it is wet
 - Use red or black colored ink anywhere on the slide
 - Write over the arrayed well areas of the slide, as this interferes with scanning.

C. Incubation

- Completely cover array area with sample or buffer during incubation.
- Avoid foaming during incubation steps.

- Perform all incubation and wash steps under gentle rocking or rotation.
- Cover the incubation chamber with adhesive film during incubation, particularly when incubation is more than 2 hours or <70 µl of sample or reagent is used.
- Several incubation steps such as step 6 (blocking), step 7 (sample incubation), step 10 (detection antibody incubation), or step 13 (Cy3 equivalent dyestreptavidin incubation) may be done overnight at 4°C. Please make sure to cover the incubation chamber tightly to prevent evaporation.

VIII. Protocol

Note: This product contains sets of reagents for different arrays. Always ensure you are using the proper glass slide and biotinylated antibody cocktail for the correct corresponding array.

The following procedure is for processing any one of the arrays in the kit.

A. Completely Air Dry The Glass Slide

1. Take out the glass slide from the box, and let it equilibrate to room temperature inside the sealed plastic bag for 20-30 minutes. Remove slide from the plastic bag, peel off the cover film, and let it air dry for another 1-2 hours.

Incomplete drying of slides before use may cause the formation of "comet tails," thin directional smearing of antibody spots.

B. Blocking & Incubation

- 2. Add 100 µl Sample Diluent into each well and incubate at room temperature for 30 minutes to block slides.
- 3. Decant buffer from each well. Add 100 µl of sample to each well. Incubate arrays at room temperature for 1-2 hour.

Longer incubation time is preferable for higher signals. This step may be done overnight at 4°C.

We recommend using 50 to 100 μ l of original or diluted serum, plasma, conditioned media, or other body fluid, or 50-500 μ g/ml of protein for cell and tissue lysates. Cover the incubation chamber with adhesive film during incubation, especially if less than 70 ul of sample or reagent is used.

- 4. Wash:
 - Decant the samples from each well, and wash 5 times (5 min each) with 150 µl of 1X Wash Buffer I at room temperature with gentle shaking. Completely remove wash buffer in each wash step. Dilute 20x Wash Buffer I with H2O.
 - (Optional for Cell and Tissue Lysates) Put the glass slide with frame into a box with 1X 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.
 - Decant the 1x Wash Buffer I from each well, wash 2 times (5 min each) with 150 µl of 1X Wash Buffer II at room temperature with gentle shaking. Completely remove wash buffer in each wash step. Dilute 20X Wash Buffer II with H2O.

Incomplete removal of the wash buffer in each wash step may cause "dark spots," the background signals higher than the spots.

C. Incubation with Biotinylated Antibody Cocktail & Wash

- 5. Reconstitute the detection antibody by adding 1.4 ml of Sample Diluent to the tube. Spin briefly.
- 6. Add 80 µl of the detection antibody cocktail to each well. Incubate at room temperature for 1-2 hour.

Longer incubation time is preferable for higher signals

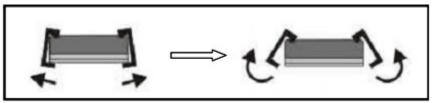
7. Decant the samples from each well, and wash 5 times (5 mins each) with 150 µl of 1X Wash Buffer I and then 2 times with 150 µl of 1x Wash Buffer II at room temperature with gentle shaking. Completely remove wash buffer in each wash step.

D. Incubation with Cy3 Equivalent Dye-Streptavidin & Wash

- 8. After briefly spinning down, add 1.4 ml of Sample Diluent to Cy3 equivalent dye-conjugated streptavidin tube. Mix gently.
- Add 80 µl of Cy3 equivalent dye-conjugated streptavidin to each well. Cover 9. the device with aluminum foil to avoid exposure to light or incubate in dark room. Incubate at room temperature for 1 hour.
- 10. Decant the samples from each well, and wash 5 times (5 mins each) with 150 µl of 1X Wash Buffer I at room temperature with gentle shaking. Completely remove wash buffer in each wash step.

E. Fluorescence Detection

11. Disassemble the device by pushing clips outward from the slide side. Carefully remove the slide from the gasket.



Be careful not to touch the surface of the array side.

- 12. Place the slide in the Slide Washer/Dryer (a 4-slide holder/centrifuge tube), add enough 1x Wash Buffer I (about 30 ml) to cover the whole slide, and then gently shake at room temperature for 15 minutes. Decant Wash Buffer I. Wash with 1x Wash Buffer II (about 30 ml) and gently shake at room temperature for 5 minutes.
- 13. Remove water droplets completely by gently applying suction with a pipette to remove water droplets. Do not touch the array, only the sides.

You may also dry the glass slide by a compressed N2 stream.

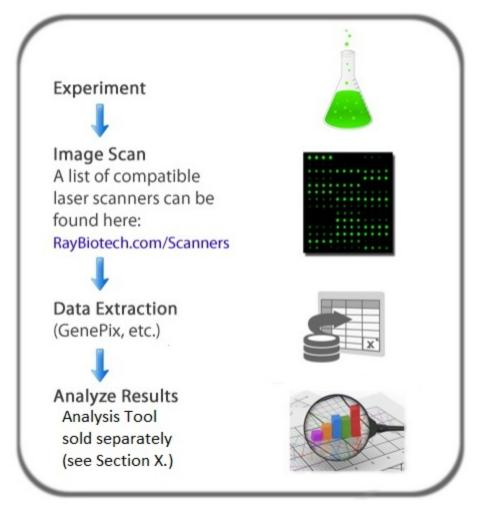
^{14.} Imaging: The signals can be visualized through use of a laser scanner equipped with a Cy3 wavelength (green channel) such as Axon GenePix or Innopsys Innoscan.

In case the signal intensity for different cytokine varies greatly in the same array, we recommend using multiple scans, with a higher PMT for low signal cytokines, and a low PMT for high signal cytokines.

F. Data Analysis

15. >Data extraction can be done using the GAL file that is specific for this array (QAH-CAA-640) along with the microarray analysis software (GenePix, ScanArray Express, ArrayVision, MicroVigene, etc.). The GAL file can be found on the product web page under the 'Files' tab.

Need help analyzing all that data? All RayBiotech array analysis tools are now free to download! Just like the GAL file, you can find this analysis tool on the product web page under the 'Files' tab. More information can be found in Section X.



QAH-INF-3

	(hINF-3 Map) Each antibody is printed in guadruplicate horizontally												
	(r	NINE-3 N	лар) Еа	ch anti	body is	printee	a in qua	arupii	cate no	rizonta	liy		
	1	2	3	4	1	2	3	4	1	2	3	4	
Α		PO	S1			PC	S2		E	BLC (C	XCL13	3)	
В	Ec	otaxin-1	(CCL'	11)	Ec	taxin-2	(MPIF	-2)	G-CSF				
C		GM-	CSF			I-3	09		ICAM-1 (CD54)				
D		IFN ga	amma			IL-1 (alpha	2		IL-1	beta	al e	
E	1	L-1ra (I	L-1 F3)		IL	-2			IL	-4		
F		IL	-5			IL	-6		IL-6sR				
G		IL	-7			IL	-8			IL-	10		
Н		IL-	11			IL-12	2p40		IL-12p70				
1		IL-	13			IL-	15			١L-	16		
J		IL-	17		1	MCP-1	(CCL2	2)		M-0	SF		
K		MIG (C	XCL9))	MIF	P-1 alpl	ha (CC	L3)	MI	P-1 be	ta (CCI	_4)	
L	MIF	P-1 delt	a (CCL	.15)		PDG	F-BB		R	ANTES	S (CCL	5)	
Μ		TIM	P-1			TIM	P-2		TNF alpha				
Ν		TNF	beta		TNF RI TNF F						RII		

QAH-CHE-1

	(h	(hCHE-1 Map) Each antibody is printed in quadruplicate horizontally												
	1	2	3	4	1	2	3	4	1	2	3	4		
Α		PO	S1			PC	S2		6	Ckine	(CCL2	1)		
В		A	xl		Be	etacellu	ılin (BT	C)	CCL28 (MEC)					
C	C	TACK	(CCL2	7)		CXC	CL16		ENA-78 (CXCL5)					
D	Ec	taxin-3	(CCL	26)	G	CP-2	CXCL	6)		G	RO			
E	H	ICC-1	CCL1	4)	H	ICC-4	(CCL1	6)		IL	-9			
F		IL-1	7F			L-18 B	P alpha	a		IL-2	28A			
G		IL-	29			IL-	31		IF	P-10 (C	XCL1	0)		
Н	- I-	TAC (C	XCL1	1)		L	IF		LIGHT (TNFSF14)					
1		Lymph	otactin		1	MCP-2	(CCL8	3)	1	MCP-3	(CCL7	7)		
J	N	ICP-4 (CCL1	3)		MDC (CCL22	2)		M	IIF			
K		MIP-3	alpha			MIP-3	3 beta		N	IPIF-1	(CCL2	3)		
L		MS	SP		N	IAP-2 (CXCL	7)	09	steopor	ntin (OF	PN)		
М	F	PARC (CCL18	3)	Plate	elet Fa	ctor 4 (PF4)	SDF-1 alpha					
Ν	1	FARC (CCL17	7)	1	ECK (CCL25	5)		TS	SLP			

QAH-CYT-4

	(h	CYT-4 M	Лар) Еа	ch anti	body is	printe	d in qua	adrupli	cate ho	rizonta	lly			
	1	2	3	4	1	2	3	4	1	2	3	4		
Α		PO	S1			PC	S2		Activin A					
В		Ag	RP		Ar	ngioger	nin (AN	G)	Angiopoietin-1 (ANG-1)					
C		Angio	statin			Cathe	prin S		CD40					
D		Crip	to-1			D	AN			DK	K-1			
E		E-Ca	dherin		E	pCAM	(TROP	1)	Fas	Ligano	(TNF	SF6)		
F	F	c gamn	na RIIB	IC		Folli	statin		Galectin-7					
G	IC	CAM-2	(CD10	2)		IL-1	3 R1		IL-13 R alpha 2					
Н		IL-1	7B			IL-2 R	alpha		IL-2 R beta					
I.		IL-	23		L	AP/TG	F beta	1		NrC	CAM			
J		PA	\ -			PDG	F-AB			Res	istin			
К		SDF-	1 beta			sgp	130			Sh	h N			
L	Si	iglec-5	(CD17	0)		ST2 (II	1 R4)			TGF-	beta 2			
М		Tie	e-2		Thro	mbopo	pietin (1	(PO)	TRAIL-R4					
Ν		TRE	M-1			VEC	GF-C		VEGF-R1					

QAH-GF-1

	(hGF-1 Map) Each antibody is printed in quadruplicate horizontally													
	1	2	3	4	1	2	3	4	1	2	3	4		
Α		PO	S1			PC	S2			A	R			
В		BD	NF			bF	GF			BM	P-4			
C		BM	P-5			BM	P-7			beta	NGF			
D		EC	GF			EG	FR			EG-\	/EGF			
E		FG	F-4			FGF-7	(KGF)			GD	F-15			
F		GD	NF		Gro	wth Ho	mone	(GH)	HB-EGF					
G		HC	GF			IGFE	3P-1			IGFI	3P-2			
Н		IGFE	3P-3			IGFE	3P-4		IGFBP-6					
1		IG	F-I			Ins	ulin		MCF R					
J	N	GFR (T	NFSR1	6)		N	-3			N	- 4			
K	Oste	oprote	gerin ((OPG)		PDG	F-AA			PIGF (PLGF)	É.		
L		SC	CF		S	CFR(CD11	7)		TGF	alpha			
Μ		TGF b	oeta 1			TGF	oeta 3		VEGF-A (VEGF)					
Ν		VEG	FR2		VEGF R3 VEGI						F-D			

QAH-REC-1

	(hREC-1 Map) Each antibody is printed in quadruplicate horizontally													
	1	2	3	4	1	2	3	4	1	2	3	4		
Α		PC	S1			PC	S2		4-1BB (CD137)					
В	A	LCAM	(CD16	6)		B7-1 (CD80))	BCMA (TNFRSF17)					
C		CE)14		C	D30 (TI	VFRSF	-8)	CD40 Ligand					
D		CEAC	CAM-1		DI	R6 (TN	FRSF2	21)		D	tk			
E	Er	ndoglin	(CD10)5)		Ert	B3			E-Se	lectin			
F		Fa	as			Flt-3 L	igand		GITR (TNFRSF18)					
G	HV	EM (TN	VFRSF	14)	10	CAM-3	(CD50))	Contactin-2					
Н		IL-	1 RI			IL-2 R (gamma	a	IL-10 R beta					
1		IL-1	17R			IL-2	1 R		LIMPII					
J	Lip	ocalin	-2 (NG/	AL)	L-S	Selectir	(CD6	2L)		LYV	/E-1			
K		MI	CA			M	СВ			NRG1	-beta 1			
L		PDGF	R beta		PE	ECAM-	1 (CD:	31)		RA	GE			
М		TIM-1 ((KIM-1))		TRA	LR3		Trappin-2					
Ν		uP	AR		VCAM-1						DAR			

QAH-CYT-5

	(hCYT-5 Map) Each antibody is printed in quadruplicate horizontally													
	1	2	3	4	1	2	3	4	1	2	3	4		
Α		PO	S1			PC	S2		Adiponectin (ACRP30)					
В		Adi	psin		Alpha	a-fetop	rotein (AFP)	ANGPTL4					
C	Beta-2	Microg	lobuliir	n (B2M		BC	AM			CA	125			
D		CA	15-3			CI	EA			C	RP			
E		Erb	B2			Fer	ritin			FS	SH			
F	GR	O alpha	a (CXC	:L1)	H	CG bet	a (HCC	Sb)		IGF	-IR			
G		IL-1	RI			IL	-3		IL-18 R beta					
Н		IL-	21			Le	otin		MMP-1					
L		MM	P-2			MM	P-3			MM	P-8			
J		MM	P-9			MM	P-10			MMF	P-13			
K	N	CAM-1	(CD5	6)		Nido	gen-1			NS	SE			
L	Ond	costatir	n M (OS	SM)	Pro	ocalcito	onin (P	CT)		Prola	actin			
Μ		PSA	-free			Sigl	ec-9		TACE					
Ν		Thyrog	lobulin			TIM	P-4		TSH					

QAH-CYT-6

	(hCYT-6 Map) Each antibody is printed in quadruplicate horizontally												
	1 2	3	4	1	2	3	4	1	2	3	4		
Α	PC	S1			PC	S2			2B4 (C	D244)		
В	ADA	AM-9		Angi	opoieti	n-2 (Al	IG-2)	APRIL					
С	BM	IP-2			BM	P-9		C5a					
D	Cathe	psin L			CD	200			C)97			
E	Cher	merin			Do	:R3			FA	BP2			
F	F	AP			FGI	-19		Galectin-3					
G	HG	FR		IF	N alpha	a/beta l	R2	IGF-II					
Н	IGF	-II R		IL-	-1 R6 (I	L-1 Rrp	(20	IL-24					
1	IL-33 (IL	1 F11)		Kallik	rein 14		Legumain					
J	LO	X-1			M	BL			Nepr	rilysin			
К	Not	ch-1			NOV (CCN3)		Osteoactivin					
L	PE	D-1			PG	RPs		Serpin A4					
Μ	sFR	P-3		Т	hromb	omodu	in		TL	R2			
N	TRA	IL R1			Trans	sferrin		WIF-1					

QAH-CYT-8

	(hCYT-8 Map) Each antibody is printed in quadruplicate horizontally													
	1	2	3	4	1	2	3	4	1	2	3	4		
Α		PO	S1			PC	S2		ANGPTL3					
В		beta I	G-H3			C	49		Cathepsin B					
C		CD)23			CH	3L1		CTLA4					
D		Dk	k-4			DP	PIV			ED/	A-A2			
E		Ep	o R			FG	F-6			FG	F-9			
F		Ga	is1			IGFE	3P-5		IL-1F5					
G		IL-1	1F6			IL-1	IF7			IL-1	IF8			
Н		IL-1	1F9			IL-1	F10		IL-1R5					
1		IL-1	7C			IL-	18		IL-20					
J		IL-	34			IL-5 R	alpha			IL-10 F	R alpha			
K		Lay	vilin			Lep	tin R			Mara	apsin			
L		M	er			MM	P-7		P-Cadherin					
Μ		Pros	tasin			PS	MA		SIGIRR					
Ν		TGF b	eta RIII		Tis	ssue Fa	actor (1	FF)	TWEAK					

QAH-CYT-10

	(hCYT-10 Map) Each antibody is printed in quadruplicate horizontally													
	1	2	3	4	1	2	3	4	1	2	3	4		
Α		PO	S1			PC	S2		ADAM8					
В		ADA	M12		E	37-H3 (CD276	5)	BMPR-IB					
C		Cadh	erin-4			Cadhe	erin-13		CD48 (SLAMF2)					
D		CD58 (LFA-3)	C	D84 (S	LAMF	5)		CE	99			
E		CD155	(PVR))	C	D229	(SLAM	3)		CEAC	CAM-5			
F		CF	XIV			Cysta	atin A		Cystatin B					
G		Cystat	in E/M			Desmo	glein 2	2	D	R3 (TN	FRSF2	25)		
Н	1	ErbB4	(HER4))		ES	AM		FGF-21					
1		Gale	ctin-2			Gale	ctin-9			IC	OS			
J	J	AM-A (CD321	1)	J.	AM-B (CD322	2)		Kallik	rein 5			
K		Mid	kine			Pentr	axin 3		I	Pref-1	(DLK-1)		
L		Sigle	ec-10		S	SLAM (CD150))		SF	P-D			
Μ		Synde	can-4		Tes	tican 2	(SPOC	CK2)	TIM-3 (KIM-3)					
Ν		TL	R4		TF	rail (t	NFSF1	0)	ULBP-1					

QAH-CYT-12

	(h0	CYT-12	Map) Ea	ach ant	ibody i	s printe	d in qu	adrupli	cate ho	orizonta	ally	
	1	2	3	4	1	2	3	4	1	2	3	4
Α	POS1					PC	S2			B	7-2	
В	BAFF R					Calc	itonin			Calsyn	tenin-1	
С	Cathepsin E				clA	P-2		Coa	gulatio	n Fact	or VII	
D	Complement MASP3			SP3		End	ocan			Ep	hA2	
E	EphB4					Ephr	in-A4			FG	-23	
F	FGF-5				Fl	t-3		GLP-1				
G		Glypi	can 2		G	M-CSF	Ralpl	na		GF	P73	
Н		HTF	RA2			IL-20 F	R alpha		IL-4 R alpha			
L I		JAN	N-C		Luteinizing hormone (LH)				Matrilin-3			
J		Meprir	n alpha			MS	PR		N-Cadherin			
К		Nepril	ysin-2			NK	p44			PAF	P-A	
L	Pepsinogen II					Prese	enilin 1			P	TH	
Μ	PYY					SC	X2		TFF3			
Ν		TFF	PI-2			TR/	ACP		Ubiquitin+1			

QAH-CYT-7

(hCYT-7 Map) Each antibody is printed in quadruplicate horizontally													
	1	2	3	4	1	2	3	4	1	2	3	4	
A		POS1				PC	S2		ACE-2				
В		Albu	umin			AM	ICA		Angi	opoieti	n-4 (Al	VG-4)	
C		BA	F F			CA	19-9			CD	163		
D	Clusterin				CR	TAM		C	XCL14	(BRA	K)		
E	Cystatin C				Dec	orin		Dkk-3					
F	DLL1				Fetu	in A		aFGF (FGF-1)					
G		FO	LR1			Fu	Irin			GAS	SP-1		
Н		GAS	SP-2		G-	CSF R	(CD1	14)	HAI-2				
1	IL-	17B R	(IL-17 F	RB)		IL-	27		LAG-3				
J		LD	LR		Pe	psinog	en I (P	G1)	RANK				
К	RBP4					SC	ST		Syndecan-1				
L	TACI				TF	PI		Thrombospondin 1					
Μ	TRAIL R2				TRA	NCE		Troponin I					
N	uPA			VE-	Cadhe	rin (CE)H5)	WISP-1 (CCN4)					

QAH-CYT-9

	(h	CYT-9 M	Лар) Еа	ch anti	body is	printe	d in qua	adrupli	cate ho	rizonta	lly	
	1	2	3	4	1	2	3	4	1 2 3 4			4
Α	POS1					PC	S2		ADAMTS13			
В	Aggrecan			Angi	otensin	ogen (AGT)	E	37-H1 (CD274	4)	
С	BMPR-IA (ALK-3)				BM	PR-II			Cadhe	erin-11		
D	CD27 (TNFRSF7)				CI	D6		Ck	beta 8-	1 (CCI	_23)	
E	CNTF			DI	VAM-1	(CD22	26)	EMMPRIN (CD147)				
F	FLRG			Follis	statin-li	ke 1 (F	SL1)	Fractalkine (CX3CL1)				
G		Gale	ctin-1			GITR I	igand		Gr	anulysi	n (LAG	i-2)
Н	IL-	1 R3 (IL	-1 R A	cp)	IL-15 R alpha				IL-17E (IL-25)			
1		IL-32	alpha		L1	CAM-2	2 (CHL	-1)	LRIG3			
J		LR	P-6		1	MEPE	(OF45)	Nectin-4			
K	Periostin					Pers	ephin			Re	enin	
L	RGM-B				RO	303			S10	8A00		
Μ	Siglec-7 (CD328)			Syndecan-3				Thrombospondin 2				
Ν	Th	rombos	spondi	n <mark>5</mark>		Tie	e-1		ULBP-2			

QAH-CYT-11

(hCYT-11 Map) Each antibody is printed in quadruplicate horizontally													
	1	2	3	4	1	2	3	4	1	2	3	4	
Α	POS1				POS2				ALK-1				
В	B7-H2					BLA	ME			BM	P-8		
С	CD28				Cor	mmon l	beta Cl	hain		Conta	ctin-1		
D	Desmoglein-1					Desmo	glein-3	3		ED	AR		
E	EphA1					Epl	hB6			Ephr	in-B3		
F	Epiregulin					FGF	-12		FGF-17				
G		FO	LR2			Gale	ctin-8		GHR				
Н		Glypi	can 1			Glypi	can 5		IFN-gamma R1				
L	I	L-22 R	alpha	1		IL-2	2BP		IL-23 R				
J		IL-3	1 RA			IL-7 R	alpha		Integrin alpha 5				
К	MDM2					Nec	tin-1			NK	p30		
L	Nogo Receptor				Not	ch-3		OSM R beta					
М	Prolactin R				RE	LT		Ryk					
Ν	5	Semapl	horin 6	D	9	Semapl	horin 7/	A	Siglec-11				

QAH-CYT-13

(hCYT-13 Map) Each antibody is printed in quadruplicate horizontally												
	1	2	3	4	1	2	3	4	1	2	3	4
Α		PO	S1			PC	S2		ACE			
В	Activin RIB				ADA	M23			Arte	emin		
C	Cardiotrophin-1				Cathe	psin V			FA	BP1		
D	FGF-20				GD	F-8			H/	\ I-1		
E	IL-27 R alpha				Insu	lin R			Kallik	rein 7		
F	LIF R alpha				Lipoc	alin-1		LTbR				
G		Meso	thelin			MF	RP			Neuro	pilin-2	
Н		Neu	turin			Nido	gen-2		Olfactomedin-2			2
. I.		p5	53			PD-E	CGF		PDGF-CC			
J		Progr	anulin	•		R	et		ROBO4			
K	Semaphorin 6B			3		Serp	in F1		SREC-I			
L	SREC-II				TL	R1		TLR3				
Μ	TPP1					TRE	M-2		TrkC			
Ν	TROY					Urom	odulin		XIAP			

QAH-CYT-14

QAH-CYT-1

	(h	CYT-14	Map) E	ach ant	ibody i	s printe	d in qu	adrupli	icate ho	orizonta	ally	
	1	2	3	4	1	2	3	4	1	2	3	4
Α		PO	S1		POS2				4-1BB Ligand			
В	Activin RIIB				Am	inopep	otidase	P2		BA	MBI	
C	BOC					Brev	/ican		Carb	onic A	nhydras	se XII
D	Car	boxype	ptidase	e A2		CD:	300c			CD	320	
E	CDNF					CI	00		CHST1			
F	CHST4					CIL	.P-1		CNTF R alpha			
G		CR	IM1		CRTAC1				CX/	ADR		
Н	Do	pa Dec	arboxy	ase		DF	PI		DSPG3			
1		EN	IR2			FC	AR		FCRL1			
J		FC	RL2			Ga	as6			GP	R56	
K		GPVI				He	psin		ILT2			
L	Jagged 2			Kirrel3				KLF4				
Μ	LAIR1			LAMP				LAMP1				
N		MD	GA1		MIS RI				1	Neurexin 3 beta		

	(h)	CYT-15	Map) E	ach ant	ibody i	s printe	d in qu	adrupli	icate ho	orizonta	ally	
	1	2	3	4	1	2	3	4	1	2	3	4
А		PO	S1		POS2				AMIGO			
В	Aminopeptidase LRAP				Amni	onless			Arylsulf	atase /	4	
С	Bcl-w				CD	109			CD	157		
D	CD34					CE	083			CLE	-1-C	
E	CLEC10A				CM	G-2			CR	EG		
F	Cystatin SN				Cytoke	eratin-8		Dectin-1				
G		Desmo	collin-3	3		Endo	glycan			Gale	ctin-4	
Н		HAF	LN1			Jago	ged 1		Langerin			
1		Lum	ican			Matri	ptase		MEP1B			
J		Nec	tin-3			0>	(40		OX40 Ligand			
K	p27					Pappa	lysin-2	2	Plexin B3			
L	Plexin D1				pro	GRP		PSA-total				
М	Reg1B			RGM-A				ROBO2				
Ν	Spinesin			TWEAK R				ULBP-3				

X. Array Data Analysis Tool

The RayBio Analysis Tools are array specific, Excel-based program that perform sophisticated data analysis on the raw numerical data extracted from the array scan. All RayBiotech array analysis tools are now free to download! Just like the GAL file, you can find this analysis tool on the product web page under the 'Files' tab.

Key features:

- <u>Simplicity</u>: Easy to operate and requires no professional training. With a simple copy and paste process, the cytokine expression levels are determined per sample.
- <u>Outlier Marking & Removing:</u> The software can automatically mark and remove the outlier spots for more accurate data analysis
- <u>Normalization</u>: The program allows for intra- and inter-slide normalization for large numbers of samples.
- <u>Two Positive Controls</u>: The program utilizes the two positive controls in each array for normalization.
- <u>User Intervention</u>: The program allows for user manual handling of outliers and other analytical data.
- <u>Analyze Multiple Slide:</u> The data for multiple slides can be inputted for easy slideto-slide comparison.

XI. Troubleshooting Guide

Problem	Cause	Recommendation
	Inadequate detection	Increase laser power and PMT parameters
	Inadequate reagent volumes or improper dilution	Check pipettes and ensure correct preparation
Weak Signal	Short incubation time	Increase incubation time or change sample incubation step to overnight
	Too low protein concentration in sample	Lessen dilution or do not dilute sample. Concentrate sample if necessary.
	Improper storage of kit	Store kit as suggested temperature. Don't freeze/thaw the slide.
	Bubble formed during incubation	Decrease amount of rocking/shaking during incubations. check for bubble formation and remove bubbles.
Uneven signal	Arrays are not completed covered by reagent	Completely cover arrays with solution for all required steps.
	Reagent evaporation	Cover the incubation chamber with adhesive film during incubation
	Overexposure	Lower the PMT or sigmal gain.
	Dark spots	Completely remove wash buffer in each wash step.
High background	Insufficient wash	Increase wash time and use more wash buffer
	Dust	Work in clean environment
	Slide is allowed to dry out	Don't dry out slides during experiment.

- Stechova, et al. Influence of Maternal Hyperglycaemia on Cord Blood Mononuclear Cells in Response to Diabetes-associated Autoantigens. Scandinavian Journal of Immunology. 2009. 70(2):149-158
- Willingham, SB et al. NLRP3 (NALP3, Cryopyrin) facilitates in vivo caspase-1 activation, necrosis, and HMGB1 release via inflammasome-dependent and -independent pathways. J Immunol. 2009; 183(3):2008-15
- 3. El Karim et al. Neuropeptides Regulate Expression of Angiogenic Growth Factors in Human Dental Pulp Fibroblasts. Journal of Endodontics, 2009; 35(6): 829-833
- 4. Souquière S. et al. T-Cell tropism of simian T-cell leukaemia virus type 1 and cytokine profiles in relation to proviral load and immunological changes during chronic infection of naturally infected mandrills (Mandrillus sphinx). J Med Primatol. 2009; 38(4):279-89
- Sharma, et al. Induction of multiple pro-inflammatory cytokines by respiratory viruses and reversal by standardized Echinacea, a potent antiviral herbal extract. Antiviral Research. 2009; 83(2)165-170.
- 6. Altamirano-Dimas, et al. Echinacea and anti-inflammatory cytokine responses: Results of a gene and protein array analysis. Pharmacuetical Biology. 2009; 47(6): 500-508.
- 7. Cheung, et al. Cordysinocan, a polysaccharide isolated from cultured Cordyceps, activates immune responses in cultured T-lymphocytes and macrophages: Signaling cascade and induction of cytokines. Journal of Ethonopharmacology. 2009; 124(1): 61-68.
- Du, et al. P2-380: Identification and characterization of human autoantibodies that may be used for the treatment of prion diseases. Alzheimers and Dementia. 2009; 4(4): T484-T484.
- 9. Van Rossum et al. Granulocytosis and thrombocytosis in renal cell carcinoma: a proinflammatory cytokine response originating in the tumour. Neth J Med. 2009; 67(5):191-4.
- Zhai, et al. Coordinated Changes in mRNA Turnover, Translation, and RNA Processing Bodies in Bronchial Epithelial Cells following Inflammatory Stimulation. Molecular and Cellular Biology. 2008; 28(24): 7414-7426.
- 11. Gao, et al. A Chinese herbal decoction, Danggui Buxue Tang, activates extracellular signal-regulated kinase in cultured T-lymphocytes. FEBS Letters, 2007; 581(26): 5087-5093. This reference validates mulitplex ELISA results for several analytes with standard ELISA test results.
- 12. Piganelli, et al: Autoreactive T-cell responses: new technology in pursuit of an old nemesis. (*Editorial Review*) Pediatric Diabetes 2007: 8: 249–251

Note: The citations listed above are for the Quantibody® product line, which is the same as the GS-Series, but include protein standards for quantitation.

XIII. Experiment Record Form

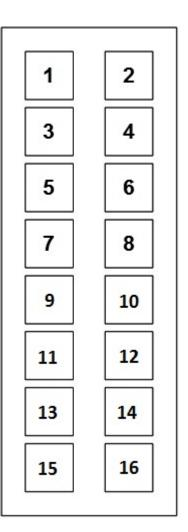
Date:_____

File Name:_____

Laser Power:_____

PMT:_____

Well No.	Sample Name	Dilution factor
1		
2		
3		
4		
5		
6		
7		
8		
9		
10		
11		
12		
13		
14		
15		
16		



XIV. How to Choose a GS-Series Array?

Species-based selection: Human (GSH-) Mouse (GSM-) Rat (GSR-) Bovine (GSB-) Canine (GSC-) Equine (GSE-) Feline (GSF-) Ovine (GSO-) Primates (GSN-) Porcine (GSP-) Rabbit (GSL-)

Function-based selection:

Adhesion Molecule Arrays	Angiogenesis Arrays	Bone Metabolism Arrays	Chemokine Arrays
Cancer Biomarker Arrays	Custom Arrays	Cytokine Arrays	Growth Factor Arrays
IGF Signaling Arrays	IL-1 Family Arrays	Immune Response Arrays	Inflammation Arrays
Interleukin Arrays	Isotyping Arrays	MMP Arrays	Obesity Arrays
Ophthalmic Arrays	Periodontal Disease Arrays	Receptor Arrays	Th1/Th2/Th17 Arrays

Cytokine Number-based selection:

Arrays are available in the GS-Series & Quantibody[®] platform to detect 660 human, 200 mouse, or 67 rat proteins. GLP-Compliant testing services are also available.

This product is for research use only.



©2015 RayBiotech, Inc