RayBio[®] Label-Based (L-Series) Rat Antibody Array L-1500 Membrane Kit

A combination of Rat L-2, L-3, and L-4 arrays

Patent Pending Technology User Manual (January 1, 2022)

For the simultaneous detection of the relative expression of 1500 rat proteins in serum, plasma, cell culture supernatants, cell/tissue lysates and other body fluids.

L-Series Rat Antibody Array L-1500 Cat# AAR-BLM-1500-2 (2 Sample Kit) Cat# AAR-BLM-1500-4 (4 Sample Kit)

Please read manual carefully before starting experiment



Your Provider of Excellent Protein Array Systems and Services

Tel: (Toll Free) 1-888-494-8555 or +1-770-729-2992; Fax: +1-770-206-2393; Website: <u>www.raybiotech.com</u> Email: <u>info@raybiotech.com</u>

TABLE OF CONTENTS

١.	Introduction and How It Works	2
II.	Materials Provided	
	A. Storage Recommendations	3
	B. Additional Materials Required	
III.	Overview and General Considerations	4
	A. Preparation and Storage of Samples	4
	B. Handling the Array Membrane	
	C. Incubation of Antibody Array	
	D. Layout of Array Membrane	7
IV.	Protocol	8
	A. Sample Purification	
	B. Biotin Labeling of Sample	
	C. Blocking and Incubations	
	D. Detection	
V.	Antibody Array Map	
VI.	Antibody Array Target Lists	
VII.	Interpretation of Results	
	A. Explanation of Controls Spots	
	B. Typical Results	
	C. Background Subtraction	
	D. Normalization of Array Data	
	E. Threshold of Significant Difference	
	F. Pathway Analysis of the Array Proteins	
VIII.	Troubleshooting Guide	
IX.	Selected References	

I. Introduction

Combining direct antigen-labeling technology with our vast library of arrayvalidated antibodies, RayBiotech has created the largest commercially available antibody array to date. With the L-Series high density array platform, researchers can now detect thousands of proteins simultaneously, obtaining a broad, panoramic view of protein expression. Our newly expanded panel includes a wide variety of metabolic enzymes, structural proteins, epigenetic markers, neuroregulatory factors, in addition to our popular list of cytokines, growth factors, receptors, adipokines, proteases, and signaling proteins. Available on both glass slide and membrane formats, this array is ideally suited for biomarker discovery studies and exploratory screens.

The first step in using the RayBio[®] L-Series Antibody Array is to biotinylate the primary amine groups of the proteins in your sample (sera or plasma, cell culture supernatants, cell lysates or tissue lysates). The membrane arrays are then blocked, similar to a Western blot, and the biotin-labeled sample is added onto the membrane array which is pre-printed with capture antibodies and incubated to allow for interaction of target proteins. After incubation with HRP-Conjugated Streptavidin, the signals can be visualized by chemiluminescence.



II. Materials Provided

A. Storage Recommendations

Upon receipt, the kit should be stored at 4 °C or below and must be used within 6 months from the date of shipment. For longer period of storage, Labeling Reagent (Item B) and Array Membrane (Item E) should be stored at -20 °C and avoid repeated freeze-thaw cycles. Labeling Reagent (Item B) should be prepared fresh before use. After initial use, Labeling Buffer, Blocking Buffer, Stop Solution, HRP-Conjugated Streptavidin, and Detection Buffers C and D should be stored at 4 °C to avoid repeated freeze-thaw cycles (may be stored for up to 3 months).

ITEM	DESCRIPTION	2 SAMPLE KIT	4 SAMPLE KIT				
В	Labeling Reagent	2 vials	4 vials				
С	Labeling Buffer	1 bottle (30 ml)	1 bottle (30 ml)				
D	Stop Solution	1 vial (50 µl)	1 vial (50 μl)				
	L-series Antibody Array	2 membranes each of Rat	4 membranes each of Rat				
E	Membranes	L-2, L-3, and L-4	L-2, L-3, and L-4				
F	4X Blocking Buffer	2 bottles (30 ml)	3 bottles (30 ml)				
	500X HRP-Conjugated						
	Streptavidin Concentrate						
К	Detection Buffer C	2 bottles (10 ml)	4 bottles (10 ml)				
L	Detection Buffer D	2 bottles (10 ml)	4 bottles (10 ml)				
G	20X Wash Buffer 1	2 bottles (30 ml)	4 bottles (30 ml)				
	Concentrate						
Н	20X Wash Buffer 2	2 bottles (30 ml)	4 bottles (30 ml)				
	Concentrate						
J-2	Spin Columns (10 ml)	4 columns	8 columns				
N/A	Plastic Incubation Trays (w/lid)	6 trays	12 trays				
N/A	2X Lysis Buffer	1 bottle (10 ml)	1 bottle (10 ml)				
Other Kit Components: Plastic Sheets							

B. Additional Materials Required

- 2-5 ml tube, small plastic or glass containers
- 50 ml conical collection tubes
- Orbital shaker or oscillating rocker
- Kodak X-Omat[™] AR film (REF 165 1454) and film processor or Chemiluminescence imaging system

- Pipettors, pipette tips and other common lab consumables
- Eppendorf tube

III. Overview and General Considerations

A. Preparation and Storage of Samples

1) Preparation of Cell Culture Supernatants

- 1. Seed cells at a density of 1x10⁶ cells in 100 mm tissue culture dishes.*
- 2. Culture cells in complete culture medium for ~24-48 hours.**
- 3. Replenish with serum-free or low-serum medium such as 0.2% FCS/FBS serum, and then incubate cells again for ~48 hours.**,[†]
- To collect supernatants, centrifuge at 1,000 x g for 10 min and store as ≤1 ml aliquots at -80°C until needed.
- 5. If you want to use cell mass for inter-sample normalization, measure the total wet weight of cultured cells in the pellet and/or culture dish. You may then normalize between arrays by dividing densitometry signals by total cell mass (i.e., express results as the relative amount of protein expressed/mg total cell mass). Or you can normalize between arrays by determining the cell lysate concentration using a total protein assay (BCA Protein Assay Kit, Pierce, Prod #: 23227).
 - *The density of cells per dish used is dependent on the cell type. More or less cells may be required.
 - **Optimal culture time may vary and will depend on the cell line, treatment conditions and other factors.
 - *Bovine serum proteins produce detectable signals on the RayBio[®] L-Series Array in media containing serum concentrations as low as 0.2%. When testing serum-containing media, we strongly recommend testing an uncultured media blank for comparison with sample results.

2) Extracting Protein from Cells

- 1. Centrifuging Cells:
 - a. Adherent Cells:
 - i. Remove supernatant from cell culture and wash cells gently twice with cold 1X PBS taking care not to disturb cell layer.
 - ii. Add enough cold 1X PBS to cover cell layer and use cell scraper to detach cells.
 - b. Cells in Suspension: Pellet the cells by centrifuging using a microcentrifuge at 1500 rpm for 10 min.
- 2. Make sure to remove any remaining PBS before adding 1X Cell Lysis Buffer (2X Cell Lysis Buffer should be diluted 2-fold with ddH₂O). Solubilize the cells at 2×10^7 cells/ml in 1X Cell Lysis Buffer.
- Pipette up and down to resuspend cells and rock the lysates gently at 2–8 °C for 30 minutes. Transfer extracts to microfuge tubes and centrifuge at 13,000 rpm for 10 min at 2-8 °C.
- Note: If the lysates appear to be cloudy, transfer the lysates to a clean tube, centrifuge again at 13,000 rpm for 20 minutes at 2-8°C. If the lysates are still not clear, store them at -20°C for 20 minutes. Remove from the freezer and immediately centrifuge at 13,000 rpm for 20 minutes at 2-8°C.
 - 4. Transfer lysates to a clean tube. Determine cell lysate concentrations using a total protein assay (BCA Protein Assay Kit, Pierce, Prod# 23227). Aliquot the lysates and store at -80°C.

3) Extracting Protein from Crude Tissue

1. Transfer approximately 100 mg crude tissue into a tube with 1 ml 1X Cell Lysis Buffer (2X Cell Lysis Buffer should be diluted 2-fold with ddH₂O).

- 2. Homogenize the tissue according to homogenizer manufacturer instructions.
- 3. Transfer extracts to microcentrifuge tubes and centrifuge for 20 min at 13,000 rpm (4°C).
- Note: If the supernatant appears to be cloudy, transfer the supernatants to a clean tube, centrifuge again at 13,000 rpm for 20 minutes at 2-8°C. If the supernatant is still not clear, store the lysate at -20°C for 20 minutes. Remove from the freezer, immediately centrifuge at 13,000 rpm for 20 minutes at 2-8°C.
 - 4. Transfer supernatant to a clean tube and store at -80°C.
- 4) Determine the total protein concentration

For optimal biotin labeling, it is necessary to determine the protein concentration in the cell/tissue lysate. We recommended using a BCA total protein assay (e.g., Pierce, Catalog # 23227).

B. Handling the Array Membranes

- Always use forceps to handle membranes and grip the membranes by the edges only.
- Never allow membranes to dry during the experiment.
- Avoid touching membranes with hands or any sharp tools.

C. Incubations of Antibody Array

- Completely cover membranes with sample or buffer during incubation and cover the Plastic Incubation Tray with the lid to avoid drying.
- Avoid foaming during incubation steps.
- Perform all incubation and wash steps under gentle rotation.

 Several incubation steps such as step 3 (sample incubation) or step 7 (HRP-Conjugated Streptavidin incubation) may be done at 4 °C overnight.

> 7 cm 9 cm

D. Layout of Array Membrane

30 columns x 36 rows

IV. Protocol

Assay Diagram

1. Cell/tissue lysates





A. Sample purification

- Note: This step removes the low molecular weight amine derivatives or unwanted buffer from samples to ensure the quality biotinylation in Steps 5–7.
 - 1. Twist to remove the bottom closure of the Spin column and loosen the cap (Do not remove).
 - Place the Spin column into a 50 ml conical collection tube, centrifuge at 1,000 x g for 3 minutes to remove the storage buffer. Discard the flowthrough.
 - 3. Wash the column three times with 5 ml ultrapure water (ddH2O) or 1xPBS (pH8.0) each, centrifuge 1,000 x g for 3 minutes to remove the

flow-through. Blot the bottom of the column to remove excess liquid, and transfer device to a new collection tube.

- 4. Apply sample on top of the resin within the next few minutes. Centrifuge at 1,000 x g for 3 minutes to collect the flow-through that contains sample. The recommended sample dilution as following:
 - Culture Media (CM): 3.5 ml neat supernatant
 - Serum/Plasma: 50 µl serum/plasma in 3.5 ml labeling buffer
 - Cell/tissue lysate: 100 µg lysate in 1 ml labeling buffer
- *Note: Each labelled sample volume is enough for at least 6 membranes following the protocol below.*
- Note: The maximal sample volume is 4 ml for each Spin Column. Do not load over 4 ml of sample into a Spin Column.

B. Biotin-Labeling the Sample

- Note: Amines (e.g., Tris, glycine) and azides quench the biotinylation reaction. Avoid contaminating samples with these chemicals prior to biotinylation.
 - 5. Immediately before use, prepare Labeling Reagent. Briefly spin down the Labeling Reagent tube (Item B). Add 100 μ I Labeling Buffer into the tube, then pipette up and down or vortex slightly to dissolve the lyophilized reagent.
 - 6. Add Labeling Reagent to the sample tube. Incubate the reaction solution at RT with gentle rocking or shaking for 30 min. Mix the reaction solution by gently tapping the tube every 5 minutes.
 - a. For labeling cell culture supernatants: Add 80 μ l of Labeling Reagent into the sample tube (for 3.5 ml supernatant).
 - b. For labeling serum or plasma: Add 80 μ l of Labeling Reagent Solution into the sample tube (for 50 μ l serum/plasma *in 3.5 ml labeling buffer*).

- c. For labeling cell or tissue lysates: Add 8 μ l of Labeling Reagent Solution into the sample tube (for 100 μ g lysate *in 1 ml labeling buffer*).
- d. For all other body fluid: Add 2 μl of Labeling Reagent Solution per 100 μg sample to be labelled.
- Note: The addition of Labeling Reagent volume is based upon the sample amount used in Step 4. If more or less amount sample is labelled, adjust this volume proportionally.
 - 7. Add 5 μ l Stop Solution (Item D) into each reaction tube. Using a new spin column, repeat Steps 1-4 of section A. Sample Purification to remove the excess non-reacted biotin reagent from each sample.
- Note: Biotinylated samples can be stored at -20°C or -80°C until you are ready to proceed with the assay.

C. Blocking and Incubations

8. Place each membrane printed side up into a Plastic Incubation Tray (provided). 1 membrane per tray.

Note: The printed membrane will have a "-" mark in the upper left corner of the membrane.

Note: Up to 4 membranes can be incubated together within one tray with proportional amount of reaction buffer. Rotate the membrane sequence at least once during sample incubation if more than one membrane is incubated in one tray.

9. Dilute 4X Blocking Buffer (Item F) with deionized or distilled water to prepare the 1X Blocking Buffer. Add 6 ml of 1X Blocking Buffer to each

membrane and cover with the lid. Incubate at room temperature with gentle shaking for 1 hour.

10. Aspirate the Blocking Buffer from each tray. Add 6 ml of diluted sample onto each membrane and cover with the lid. Incubate at room temperature with gentle shaking for 2 hours.

Note: It is recommended to use 10-20 folds diluted biotin-labeled culture supernatant, 10-20 folds diluted biotin-labeled serum/plasma, 100 folds diluted biotin-labeled cell/tissue lysate, or 10-20 folds for other body fluids. Dilute sample using 1X Blocking Buffer. The optimal concentration of sample used will depend on the abundance of target proteins. The samples can be concentrated if the overall signals are too weak. If the overall signals are too strong, the sample can be diluted further.

Note: Incubation may be done at room temperature with gentle shaking for 2 hours or overnight at 4°C.

- 11. Dilute 20X Wash Buffer 1 (Item G) with deionized or distilled water to prepare the 1X Wash Buffer 1. Aspirate the samples from each tray and then wash by adding 20 ml of 1X Wash Buffer I at room temperature with gentle shaking (5 min per wash). Repeat the wash 2 more times for a total of 3 washes.
- Aspirate the 1X Wash Buffer 1 from each tray. Dilute 20X Wash Buffer 2 (Item H) with deionized or distilled water to prepare the 1X Wash Buffer 2. Wash 3 times with 20 ml of 1X Wash Buffer 2 at room temperature with gentle shaking.
- 13. Aspirate the 1X Wash Buffer 2 from each tray.
- 14. Prepare the HRP-Conjugated Streptavidin. Briefly spin down the tube containing the 500X HRP-Conjugated Streptavidin (Item I)

immediately before use. Dilute the 500X HRP-Conjugated Streptavidin with 1X Blocking Buffer to prepare the 1X HRP-Conjugated Streptavidin. Pipette up and down to mix gently. Add 6 ml of 1X HRP-Conjugated Streptavidin to each membrane.

Note: Ensure that the vial containing the 500X HRP-Conjugated Streptavidin is mixed well before use, as precipitation can form during storage.

15. Incubate at room temperature with gentle shaking for 2 hours.

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

16. Wash as directed in steps 11 through 13.

D. Detection

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

- 17. For detection of 2 membranes, add 4.2 ml of Detection Buffer C and 4.2 ml of Detection buffer D into a tube and mix both solutions. Drain off excess wash buffer. Place membrane antibody side up (There is a "-" symbol on the top left corner of each membrane) on a clean plastic plate or its cover (provided in the kit). Pipette 4 ml of the mixed Detection Buffers onto each membrane and incubate at room temperature for 2 minutes with gentle shaking. Ensure that the detection mixture is evenly covering the membrane without any air bubbles.
- 18. Gently place the membrane with forceps (antibody side up) on a plastic sheet (provided) and cover the membrane with another plastic sheet. Gently smooth out any air bubbles. Avoid using pressure on the membrane. Work as quickly as possible.

- 19. The signal can be detected directly from the membrane using a chemiluminescence imaging system or by exposing the array to x-ray film (we recommend using Kodak X-Omat[™] AR film) with subsequent development. 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 the exposure time (e.g., 5–30 seconds). If the signals are too weak, increase the exposure time (e.g., 5–20 min or overnight) or re-incubate membranes overnight with 1X HRP-Conjugated Streptavidin, and repeat detection on the second day.
- 20. Save membranes at -20 °C to -80 °C for future reference.

V. Antibody Array Maps

A. RayBio[®] Rat Antibody Array L-2, L-3, and L-4 Array Map



VI. Antibody Array Target Lists

A. RayBio[®] Rat Antibody Array L-2 Target List

Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name
1	11b-HSD1	73	CD1d1	145	DYRK2	217	GLRX3	289	IPP2	361	Notch-1	433	SEMA4 D
2	14-3-3 ensilon	74	CD200	146	DVPK 3	21.8	Givovalase1	290	lc let-1	362	Notch-2	434	SEM 45.4
2	14.2.2 mm	75	00200	147	EC M1	210	Cluevelese 2	201	lagged 1	262	Notch 2	4.25	SEMAZA
5	14-3-3 6.8	75	0022	147	ECIVI1	215	Glyoxalase2	291	1988ECLT	303	NDCO	435	SLIVIA/A
4	14-5-5 Sigma	/0	CD226	148	EEAL	220	Giypican 5	292	JAIVI-A	304	INPU2	450	Serpinirz
5	14-3-3 theta	77	CD27	149	EFNA1	221	gp130	293	JAM-C	365	NRAGE	437	SerpinA1
6	4-1BB	78	CD276	150	EFNB1	222	GPT	294	JNK1	366	Nrf2	438	SerpinE1
7	A2 B5	79	CD300f	151	EGF	223	Gpt2	295	JNK2	367	NRXN1 beta	439	SerpinF1
8	ACACA	80	CD300LG	152	EGFR	224	GP X1	296	K DR	368	Olfactomedin-1	440	SH2 B1
9	ACTC1	81	CD31	153	el F5 A	225	GP X3	297	Keap1	369	OLR1	441	SHIP2
10	Actin	82	CD34	154	FMP	226	GranzymeB	29.8	Kirrel 3	370	Osteocalcin	442	SHP-1
11	Activia D2A	02	CD3.9	155	ENO1	220	CDB2	200	KITCIS VIVB1	371	03/00/01/01	442	SIGND1
12	ADAM10	0.0	00001	155	Entruin	227	CDINDA	235	KENDI KNC1	371	- 27	443	SIGNAL SIDD alaba
12	ADAMIU	84	CD39L1	156	Eotaxin	228	GRINZA	300	KINGI	372	p27	444	SIRP alpha
13	ADAMTS1	85	CD4	157	EphA5	229	GRK1	301	LAIR1	373	p38 gamma	445	SLAMF1
14	Adiponectin	86	CD47	158	EphB1	230	GRK2	302	LAR	374	p53	446	SLC4A1
15	aFGF	87	CD48	159	EphB6	231	GRK5	303	LAYN	375	p55PIK	447	Slit3
16	Agrin	88	CD5L	160	Ephr in-A2	232	GRO al pha	304	LDHA	376	PAK1	448	Smad 3
17	AIF	89	CD6	161	Ephrin-B2	233	GRP75	305	Legumain	377	PAK7	449	Smad 7
18	ΔΚ1	00	CD63	162	FP alpha	23.4	HAAO	306	Lentin	378	Pay 7	450	SMC1
10		01	0000	162	Enepo	1254		207	Loptin Decentor	270	D Codborin	450	Sortilin
19	ALCAIVI	91	00700	105	ERDD2	235	HADPI	307	Leptin Receptor	379	P-Caunerin	451	SOITHIN
20	ALK-7	92	CD79B	164	ERBB3	236	HGF	308	UF	380	PCDH-17	452	SOST
21	Alpha-Actinin 1	93	CD8 alpha	165	Erythropoietin	237	HIF-1 alpha	309	LIFR	381	PCK1	453	SOX1
22	Alpha-Synuclein	94	CD83	166	Ets-1	238	HO-2	310	LI LRA5	382	P DGF-BB	454	SOX10
23	Ameloblastin	95	CD86	167	Ezrin	239	HP RG	311	LILRC2	383	P DGFRA	455	SOX2
24	AMPK alpha 2	96	CD93	168	F2	240	HPX	312	Lipocalin-2	384	Pentraxin 2	456	SP-D
25	Androgen R	97	CDC25B	169	F3	241	HSP20	313	LMW-PTP	385	Peroxiredoxin 6	457	Src
26	AN/CDT1	0.9	CDC37	170	EARD 1	2/12	HSP27	31.0	IDHVIS	3.86	DEKM	459	STAT2
20		30	CDC5/	174	FADP1	242	H0P40	245		200	PERMI	450	SIAIS
27	Annexin AL	99	CDHI	1/1	FABP 2	245	HSP4U	515	LKP-4	38/	PGC	459	syndecan-2
28	Annexin A4	100	CDH2	172	FABP 3	244	HSP60	316	LTBR	388	plgR	460	Syntax in 1A
29	Annexin A7	101	CDNF	173	FABP4	245	HSP70	317	LTF	389	PIM2	461	TAFA5
30	Annexin V	102	CES1	174	FABP 5	246	HSP90	318	Lyn	390	PKA C a/b	462	Talin1
31	APE	103	CF X IV	175	FAK	247	HSPA8	319	MAG	391	PKC	463	TCK-1
32	API P-1	104	CHMP2 B	176	FC AR	248	HSPH1	320	Matrilin-3	392	PKC a	464	TC-PTP
33		105	Chordin	177	ECCP1	2/0	HtrA2	321	MBL-2	3.03	PKC i/L/2	465	TDP-//3
33	APRIL Assistent 1	105	CIDI	170	FCGKI	249	IDC	321	IVIDL-2	393	P NC 1/1/2	405	104-43
54	Arginase 1	106	CIBI	1/8	FEIUB	250	IUS	522	IVICAIVI	394	PKIVIZ	400	IF
35	AR 14	107	CLEC4A2	1/9	FGF-12	251	IFNA5	323	MCP-3	395	PLAUR	467	I GE-beta RIII
36	ASAH2	108	CLEC4B2	180	FGF-21	252	IFN-alpha	324	MEK2	396	Plexin A4	468	TGM2
37	B3GNT2	109	Clusterin	181	FGFR4	253	IFN-gamma	325	MIF	397	PON3	469	THBD
38	BAFF	110	CNTF	182	Fgr	254	IFN-gamma R2	326	MIG	398	POR	470	Thioredoxin-2
39	BAK	111	CO5	183	Fibromodulin	255	IGF-1	327	MIP-1 alpha	399	PP2ACS	471	TIE-2
40	BCAM	112	COLEC12	184	EKBP12	256	IGEBP-5	328	MIP-3 beta	400	PP2C alpha	472	TIM-1
4.1	Rel-10	112	Complexip-2	105	EV 80126	250	10101	220	ME	4.01	DDA1	172	TNE allaba
41	DCI-10	113	Complexitie 1	105	FK0F12.0	257	IGGF0	325		401	PPD2D4	473	TNE D1
42	BCI-2	114	Contactin-1	180	FKBP15	258	IK D-Deta	550	IVIIVIP-2	402	PPPZR4	4/4	INF-KL
43	BCI-W	115	Contactin-2	187	FKBP25	259	IKK	331	MMP-8	403	PRDX 2	475	INFRSF11A
44	BcI-xL	116	Contactin-4	188	FKBP51	260	IL-1 beta	332	MMP-9	404	PRDX1	476	TNFSF9
45	beta 2-M	117	Cortactin	189	FKBP52	261	IL-1 RA	333	MO G	405	PRDX4	477	Tollip
46	beta IG-H3	118	CPA1	190	FLIP	262	IL-10	334	MP1	406	Pref-1	478	TPT1
47	bFGF	119	CPA2	191	FLT1	263	IL-11 R alpha	335	MPO	407	PRL-3	479	TRAF-2
48	BID	120	CPB1	192	Elt-3 Ligand	264	II-12 p70	336	MST1	408	P RI 8 44	480	TRAF-3
40	BIK	121	C DEL D1	102	Follictatio	265	IL-12 D-2	227	NCAM-1	100	PROCE	/ 01	Transgelin
45	PLVDP	121	CDELDO	10.4	EOLD4	205	IL 15 NG2	220	NCD2	405	Brolactia	401	TDEM 4
50	DLVKD	122	C RELUZ	194	FORT	200	11.1.7 Kd	338	NEDDA	410	Drepardia	402	TDUDE
51	BIVIP-2	123	UrkL	192	FRK	20/	IL-1 / RC	559	INE DD4	411	Properain	483	IKHDE
52	BMP-7	124	CRP	196	FRS2	268	IL-18	340	NEDD8	412	PSAP	484	TrkA
53	B-raf	125	CRYAB	197	GABRA4	269	IL-18 BPc	341	Nephrin	413	PSMA1	485	TrkB
54	BST1	126	CSF1R	198	GAD1	270	IL1 R1	342	Nestin	414	PSMA2	486	TrkC
55	BTLA	127	CTACK	199	Galectin-1	271	IL1 R2	343	Netrin-1	415	PTK7	487	TWEAK R
56	C4.4A	128	CTGF	200	Galectin-3	272	IL-2	344	Neurexophilin-1	416	PTP1B	488	UCH-L1
57	Cadherin-4	120	CTHPC1	201	Galectin-4	273	II-2 Pa	345	Neuritin	417	p\/p	4.80	LICH-13
50	CADM2	120	CTIAA	201	CAPDU	273	11.2 PC	3/4	Neurocan	/10	D\/DL2	400	
50		100	CILAN	202		2/4	1L-2 KG	340	Neurocari	410	P V KLZ	4.50	
59	Calcineurin A	131	Cubilin	203	Gas 1	2/5	11-21	347	Neurofascin	419	каіА	491	UNC5H2
60	Calcineurin B	132	CXCL10	204	GDF-3	276	IL-22	348	Neurogranin	420	RALT	492	VAMP-2
61	Caspr 2	133	CXCL16	205	GDF-8	277	IL-23 p19	349	Neuroligin-1	421	RANTES	493	VHR
62	Catalase	134	Cyclophilin A	206	GDNF	278	IL-31	350	Neuroligin-2	422	RBBP4	494	Vinculin
63	Cathepsin B	135	Cyclophilin B	207	GFAP	279	IL-4	351	Neuropilin-1	423	RBP4	495	VSIG1
64	Catheosin C	136	Cystatin C	208	GFRA1	280	IL-4 R	352	Neuroplastin 65	424	Reg III	496	WEDC2
65	Catheosin F	137	Outochrome C	200	GEPA2	281	1.4	35.2	NEATC3	4.25	Ren3R	407	Wet5a
60	Cathornel	120	Description	203	CERAD	201	11-0	25.4	NET	425	Degin 1	400	VILDO
00	Cathepsin L	138	Decorin	210	GERAS	262	11-7	554	INF*L	420	Keill I	498	AIAP
67	Cathepsin X	139	DEP-1	211	GGT1	283	IL-/ Ra	355	NM23-H1/H2	427	RHD	499	X PNP EP2
68	Caveolin-2	140	DGK-epsilon	212	GH	284	IL-9	356	nNOS	428	RO BO1	500	Zyxin
69	CCK-A R	141	DHFR	213	GIT1	285	IL-9 R	357	NNUP85	429	RO CK2		
70	CCL26-Like	142	Dkk-3	214	GITR	286	ILK	358	Noggin	430	SDC1		
71	CD13	143	DIL1	215	GLA	287	ILK AP	359	Nogo-A	431	Secretagogin		
72	CD14	144	D047	216	GLC1	28.8		360	None	432	SEMAAC		
12	0014	7-4-4	00117	210	0.01	200		550	nope	-12	JENNAG		

B. RayBio[®] Rat Antibody Array L-3 Target List

							<u> </u>						
Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name
1	A2M	73	CHGB	145	FGFR5	217	IL-1 R4	289	Neurturin	361	Quiescin Q6	433	TCP1
2	AADE	74	Child 1	14.6	500	21.0	11.1.00	20.0	Ni dener 1	2.62	DACE	4.2.4	TDIFO
2	AAKE	/4	CIKI	140	rgg	210	IL-I KO	290	Nidoger-1	302	KAGE	404	IDIFZ
3	ABCF1	75	Chymase	147	FH	219	IL-11	291	Nidogen-2	363	Ras	435	TECK
4	ACAT1	76	CINC-2	148	Fibronectin	220	IL-12 p40	292	NIT2	364	RELM beta	436	Tenascin X
5	Activin A	77	CINC-3	140	Ficolin-2	221	II-12 PB1	203	NNT	365	Pecistin	437	TEE1
		70	Cine J	45.0	TICO111 2	221	10 12 101	200	NOV	205	0500	400	7550
ь	ADAMISIU	/8	Cinguiin	150	FLGZ	222	IL-13	294	NOV	300	REV3L	438	IFF2
7	ADAMTS15	79	CIP29	151	FOXN3	223	IL-15	295	NP B	367	Rheb	439	TGF-beta 1
8	ADAMTSI 2	80	Claudin-3	152	Fractalkine	224	II-16	296	NPTXR	368	RNASE6	440	TGE-beta 2
0	Aggregion	01	Cloudin 4	152	Eximpled 1	225	11 17 4	207	NID2/C2	260	DO PO/	4.41	TCE both 2
9	Aggrecari	01	Claudin-4	155	FI1221eu-1	225	IE17A	297	INKOCO	209	KU DU4	441	IGF-Deta 5
10	AHCY	82	CNPY2	154	Frizzled-4	226	IL-17C	298	Nrf1	370	ROR1	442	TGF-beta R1
11	AHSG	83	CNTFR	155	Frizzled-5	227	IL-17D	299	OCT3/4	371	RP1	443	TGF-beta R2
12	Δk+2	84	COI 19A1	15.6	Frizzled-6	228	11-10	300	Orevin A	372	DDI12	4.4.4	TIMP-1
12	76.02		COLIDAI	150	111221000	220	1015	500		572	111 112		TIN 1
13	Albumin	85	COLL	157	Frizzled-/	229	IL-2 R beta	301	OSCAR	3/3	RPL23A	445	TIMP-2
14	AMPKa1	86	CPE	158	FSTL1	230	IL-24	302	OSM	374	RP LP0	446	Titin
15	ANGPTI 2	87	CRADD	159	Galanin	231	11-27	303	Ostepactivin	375	RPS13	447	TK 1
10	ANOTTE2	00	CRADD	100	CACD 1	201	10.200	20.4	Ostessadhasia	270	00014	440	TID1
10	ANGPILS	00	CKED	100	GASP-1	252	1L-26 D	504	Osteoadne in	576	KP 514	440	TUKI
17	ANK RD9	89	CRF21	161	GASP-2	233	IL-3	305	Osteoprotegerin	377	RPS15A	449	TLR3
18	ANXA6	90	CRHBP	162	G-CSF R	234	IL-3 R beta	306	p130Cas	378	RPS23	450	TLR4
10	ADBA2	01	CrkPS	163	CDE-15	235	II5	30.7	p21	370	DD S3 A	// 51	TMEEE1
19	AFUAZ	51	CINNO	105	001-10	235	IL-5	307	pzi	375	NF JUA	401	INVIETTI
20	ApoA1	92	CRTAC1	164	GDF-5	236	IMP2	308	P4HB	380	RPS5	452	TMEFF2
21	ApoA2	93	CRTAM	165	GFRA4	237	INSL3	309	Pappalysin-1	381	RPS8	453	TMEM223
22	ApoB	94	CRTH-2	166	GHR	238	Inuslin	310	PCAP	382	RPS9	454	TOMM70A
22	Ac -5	05	Countin	167	CVAH	220	LTAC	211	DODE 1	202	00501	455	TRIC
25	Apot	32	cryptic	10/	GKINI	259	I-TAL	511	FUPE-1	385	RKEDI	400	1812
24	ARHGAP1	96	CSE1L	168	GLI-2	240	Jak2	312	P D-1	384	RSF1	456	TPP1
25	ATG5	97	CSK	169	GLIPR2	241	Kallikrein 10	313	P D-ECGF	385	RUSC1	457	TRADD
26	ΔΤΡΟ	0.8	CTNND1	170	Glu#1	24.2	Kallikrein 11	31.4	PDGE-AA	3.86	\$100410	458	TRAIL P2
20	Packets	20	CHANDI	174	Cluth	242	Kelliki- 7	345	DOI-MA	207	5100A10	450	TDALL
27	B3GAI1	33	CXCR2	1/1	GIUT2	243	Kai i ikrein 5	515	PDGF-C	58/	5100A11	459	IKAM
28	B4GaIT1	100	CXCR4	172	Glut4	244	Kallikrein 6	316	PDGF-D	388	S100 A9	460	TRIM14
29	B7-1	101	CXCR7	173	Glut5	245	KIF5 B	317	PDGERB	389	S-100b	461	Tropomyosin 3
20	0.1	102	Ouclin D1	174	CM2A	245		210	DDUME	200	CPD 1	460	торал
50	D/ •F12	102	Cyclin D1	1/4	GIVIZA	240	LAWAS	210	PDLINIS	290	306-1	402	IKKAP
31	BAFF R	103	Cyclophilin F	175	GM-CSF	247	LAMP	319	PDZD2	391	SCF	463	Trypsinogen-2
32	Bax	104	Cystatin A	176	GP2	248	LASP 1	320	PENK	392	SCF R	464	TSLP
33	BDNE	105	Cystatin B	177	gn340	249	1 BP	321	Pentraxin-3	393	SDF4	465	TSP-1
24	hoto NCE	105	Oustatin D	170	C0D1	250	1 - 45 - 1	222	Desilipin 2	204	Contin 7	105	TCD 2
34	Deta-Indi	100	Cystatin D	170	GPD1	250	Lay-1	322	Perinpines	3.34	Septite/	400	TOP 4
35	BLAIVE	107	Cystatin E	1/9	GPK-39	251	Letty-A	323	Peroxiredoxin-3	395	SERBPI	46/	158-4
36	BLMH	108	Cystatin S	180	GranzymeA	252	LHPP	324	PF4	396	Serpin A3	468	TTF1
37	BMP-1	109	DAK	181	Granzyme M	253	LIX	325	PFAS	397	Serpin A5	469	TUBA6
38	BMP-15	110	DCI	182	GRHPR	254	LPS	326	PEDN6	398	Serpin B5	470	TWF2
20	BMD-0	111	DCVP	102	CPD	25.5	IDC1	227	PHCDH	200	Serpin C1	471	TVNDC15
35	DIVIP-9	111	DUAN	10.5		255	1001	327	Pindon	333	Jeipinter	4/1	TXNDCID
40	BNIP2	112	DIL4	184	GSK-3 beta	256	LKP-6	328	PICCOIO	400	SEI	472	TXNDC5
41	BOLA2	113	DMGDH	185	GSN	257	L-Selectin	329	PIK3 R2	401	sFRP-4	473	TYRO10
42	BTC	114	DSCAM	186	GSR	258	LUZP 1	330	PINCH1	402	SH3BGRL3	474	UBC9
43	BTE3	115	DSG1	187	GSTM1	25.9	Lymphotactin	331	ΡΙΡΔΚ2Δ	403	SHBG	475	Ubiquitin
45	01	445	5001	107	001011	255		001	111 4827	400	3100	475	
44	Clq	116	EDA-A2	188	GS101	260	MACAM-1	332	PLA2G1B	404	SHOX	4/6	Ubiquitin+1
45	C1s	117	EDAR	189	GULP1	261	MAN1	333	PLD4	405	Siglec-1	477	UNC45A
46	C3a	118	eEF2	190	HAI-1	262	McI-1	334	Plexin B2	406	SLC38 A10	478	UNC5H4
07	(53	110	EG-VECE	101	Hantoglobic	263	MCP-1	325	PICE-2	407	SUTDV1	470	IID A
47	004	115	LG-VEGF	191	Laplogroun	203	MOP-1	335	FIGF-2	407	SUIKKI	4/5	UPD C1
48	CAI	120	eit4E	192	HB-EGF	264	MCP-5	336	PLS3	408	SPT	480	URU C1
49	CA2	121	EMAP-II	193	HEG1	265	MDC	337	P NP	409	SLURP1	481	USP2
50	CA3	122	Endothelin	194	Hepassocin	266	MEP1A	338	POMC	410	Smad 1	482	Uteroglobin
51	Calhindin D	123	Fotavin-2	105	HEYR	267	Mesothelin	320	PON1	Δ11	Smad 4	483	VΔP-1
51		125	EUGATIF2		LICE .	207	MCSOUTCHT	333	10/11	411	0	405	VAC -1
52	cardiotrophin-1	124	EphA1	196	HGFA	268	MICB	340	PP	412	Smad 5	484	VAP-A
53	Cathepsin A	125	EphA2	197	Histone H2AY	269	MIP-3 alpha	341	PPP1CC	413	Smad 8	485	VARS
54	CC128	126	EP HX2	198	hnRNPI	270	MISRU	342	PRAT4B	414	Somatostatio	486	VDAC1
5-	00.02	107	Enirmulia	100	Howb?	270	Mitofusia 2	2/2	DDELD	A1E	COVE	107	VECE
55	00.63	12/	- cpireguin	199	10x03	2/1	wittorusin 2	343	PKEUP	412	2072	+0/	VEGP
56	CCR4	128	ERRa	200	HOXD11	272	MKK3	344	Prolactin R	416	SPARC	488	VEGF-B
57	CCT3	129	E-Selectin	201	HSP10	273	MKK4	345	ProSAAS	417	SPINK7	489	VEGF-C
58	CD133	130	EVC2	20.2	HSP47	274	MMP-10	346	Prostasio	418	SPTBN5	490	VEGER3
50	00100	104	Ener- IV	202	LITOAA	274	MAMP 40	247	Dest-1-7	440	00700	4.04	MURA
59	CD23	151	Factor IX	203	TIKAL	2/5	IVIIVIP-13	547	Protein Z	419	351KZ	491	VILIP3
60	CD24	132	Factor V	204	HVEM	276	MMP-16	348	Prouroguanylin	420	STXBP2	492	Visfatin
61	C D2AP	133	Factor VII	205	ICAM-1	277	MMP-7	349	PRR4	421	SVEP1	493	Vitronectin
62	CD30	134	Factor VII	20.6	ICAM-2	27.8	MRP 1	35.0	PRRC2 A	422	SYK	494	W/ARS
62	CDIOLiceard	105	EAM2C	200	IDE	270	Multimeria 2	254	PDTNP	122	CV0.11	405	WICD 1
05	CD40 Ligand	102	FAIVISC	207	IDE	2/9	Murumerin 2	351	PRIND	425	2101	432	WISP-1
64	CD9	136	Fas	208	IFN-beta	280	MuSK	352	P-selectin	424	IACI	496	WISP-2
65	CD90	137	Fas Ligand	209	IFNGR1	281	MyBPC3	353	PSMB1	425	TAGLN2	497	XP D
66	CDC14	138	FGF-11	210	IGF BP -2	282	NACA1	354	PSMD2	426	TALDO	498	XPG
67	CEH	120	EGE-20	211	LOSE4C	283	NADK	355	PSMDD	4.27	TALDO1	400	VV1
67	ori	135	FOF 20	211	103140	200	NACO	355	PONID9	427	THEOT		702114
68	CFI	140	FGF-23	212	IL-1 alpha	284	NAGPA	356	PSIME1	428	Talin-2	500	ZC3 H4
69	CFL1	141	FGF-9	213	IL-1 F10	285	NAP RT1	357	PTHLP	429	TARC		
70	CGA	142	FGF-BP	214	IL-1 F5	286	NeuroD1	358	PTMA	430	TARS		
71	CHCHDS	1/12	EGED1	215	11-1 55	287	Neurolusio	350	pvv	A 21	TCA-3		
/1	CHICHUO	140	TOPNI	213	10-110	201	Neur Orysin	555	F 11	431	TCA-5		
/2	Chemerin	144	I FGFR2	216	I IL-1 F9	288	Neuropilin-2	360	QARS	432	Icf20		

C. RayBio[®] Rat Antibody Array L-4 Target List

Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name
1	14-3-3 heta	73	CD40	145	ECGR34	217	IMPAD1	289	NDEIP1	361	PNUTS	433	SHIP
2	14.2.2	74	0010	140	FCCRT	210	IMPDU2	200	Nantia 2	202	0024	434	2 0 1 2
2	14-5-5 gamma	/4	0044	140	FUGRI	210	TMFDH2	290	Necun-5	502	FFZA	454	387-2
3	A1BG	75	CD51	147	Fen 1	219	Inhibin beta	291	Nesfatin-1	363	PPM1B	435	SIGNR3
4	A1M	76	CD59	148	Filamin A	220	iNOS	292	Nesprin2	364	PPM 1L	436	Six3
5	a Amyl ase	77	CDC25A	149	FKBP38	221	Intelectin-1	293	Neurogenin-2	365	PPP1R9B	437	SMAGP
6	ACE2	78	CDC25C	150	FoxA2	222	IRF1	294	Neuroglycan C	366	PRCP	438	SMOC-1
7	ACIR	70	CDK1	151	Eev P2	222	IDS1	205	NGER	267	PPDVE	420	SMURE2
	Act	/5	CONI	151	10/15	225	1101	255	NGIN	507	11073		5000012
8	ACTN2	80	CDK2	152	FPRP	224	IRS2	296	Nicalin	368	PRG2	440	SNAP25
9	ADAM17	81	CEACAM1	153	FSTL4	225	ITGA8	297	Ninjurin-2	369	PRNP	441	SOD1
10	ADAM 9	82	CELF1	154	FUCA1	226	ITGB4BP	298	NIPP1	370	Prohibitin	442	SOD2
11	ADNP	83	CES3	155	Evn	227	ITGB5	299	NKX2.2	371	Prss21	443	SOD-3
12	40.002	9.4	CHORDC1	155	(29P	220	ITGRE	200	NIRP10	272	DSD. GE	444	SPOCK2
12	AD1102	07	CHORDCI	157	0.00	220	17000	201	NDC1	372	DTEN	445	SI OCK2
15	AFF	65	CKBB	157	GBPD	229	TIPRS	301	NPCI	5/5	PTEN	445	30,311/11
14	AGT	86	CLEC1B	158	GABAB R1	230	JAB1	302	NR3C1	374	PTGDS	446	SR-AI
15	Akt1	87	CLEC5A	159	GABAB R2	231	Jak1	303	NrCAM	375	PTGES3	447	ST3GAL2
16	ALDH2	88	COL1A1	160	GABRA1	232	JI P1	304	NS E	376	PTP gamma	448	STAT5b
17	ALOX5	89	CO16A1	161	GABRAS	233	Kallikrein 7	305	NT5F	377	PTP-MEG2	449	STAT6
10	alaba 2u Globulia	90	COLEC10	162	GALNE2	224	KCNP2	205	NUAK1	270	DTDDM	450	STI1
10	arpna zu-dioburin	50	COLLCIO	102	GALINIZ	234	KONDZ	300	NURKI	370	FTERIVI	450	3111
19	ALPP	91	Complexin-1	163	gamma Catenin	235	KCNC1	307	Nucleostemin	3/9	PIPKU	451	SIIM1
20	AMBP	92	Contactin-3	164	GATE-16	236	KIAA1967	308	NXPH3	380	PVRL1	452	ST K3
21	AMH	93	COPZ1	165	GBL	237	Klotho beta	309	Oligodendrocyte Marker O1	381	Q DPR	453	Substance P
22	Amphiphysin	94	CPEB3	166	GDF1	238	KMO	310	Oligodendrocyte Marker O4	382	Rab11A	454	SUM 03
23	AMPK beta 1	95	CPM	167	GDF7	239	KOR	311	Oncomodulin	383	Rab27a	455	SUSD2
24	ANG 2	95	CSNV1A	160	GD11	240	KPND1	212	OPPM1	394	RAP7A	456	Synantotamin 1
24	And-2	30	CONKIA	100	3011	240	NEND1	312	OFAMI	304	nAD/A	-30	Synapio (agrinin-1
25	Angiogenin	97	CSNK1D	169	Gephyrin	241	Kynureninase	313	Osteopontin	385	RAC1	457	Syndecan-3
26	Annexin A11	98	CSNK1E	170	GLUD1	242	LAMC1	314	0V-6	386	RACK1	458	Syntaxin 1B
27	Annexin A2	99	CSNK1G	171	Glycine R	243	Laminin S	315	P20Sb3	387	Rad17	459	Syntaxin 7
28	ApoH	100	CSNK2B	172	GOLGB1	244	LC3B	316	p38 alpha	388	Raf-1	460	Svntaxin 8
29	ARC	101	CSRP1	173	GPID1	245	IHX5	317	n70.56 Kinase	389	Ran1A/R	461	Syntaxin BP1
20	ATE2	102	CVCP2	174	GPRE4	245	LING	210	PARP	200	Pap2A/P	462	T Call Resenter alpha Chain V
50	AIF2	102	CACHS	1/4	GFR04	240	UFG .	510	FABE	330	парадув	402	r cen keceptor alpha charry
31	AIF6	103	CXCRb	1/5	GPX2	247	Lipin 2	319	PAK4	391	RECQ4	463	I CEB2
32	ATG3	104	CYTL1	176	GPX4	248	LMAN1L	320	PAK6	392	REG4	464	TCP1 eta
33	ATM	105	DARC	177	GRB7	249	LMNA	321	Pannexin-1	393	Relaxin R1	465	Tenascin R
34	Axin-1	106	DARPP-32	178	GRP78	250	LOK	322	Park7	394	RELM gamma	466	TfR
35	B7-H4	107	DDC	179	GSK-3 al nha	251	IRPAP	323	PARI	395	RGM-B	467	TGN38
26	PAG4	100	DDT	100	LCDD	252	lumiene	224	Pasualhumia	205	PCM C	460	TU
27	BACC	100	DDV1	100	HADDO	252	Lunicali	225	Pavillia	307	RUN-C	400	Theorie
57	BAGB	109	DDX1	101	HABP2	255	Lysozyme	525	Paxillin	59/	KHUG	469	Themis
38	BAM BI	110	DEFA6	182	HAO-1	254	LYVE1	326	PCBP2	398	RIBP	470	Thioredoxin-1
39	BarX1	111	DGK-gamma	183	HBB	255	MAD2L1	327	PCDH12	399	RIPK1	471	Thrombopoi eti n
40	BCHE	112	DGK-theta	184	HCLS1	256	MafB	328	PCK2	400	RKIP	472	TLR7
41	Beclin 1	113	DISC 1	185	HDAC2	257	MAP4K4	329	PCNA	401	RNASE4	473	TOP2B
42	hete Actio	114	Disk 1	100	HDAC4	250	Matrilia 4	220	BCEKB	402	DNE2	474	TOP
42	Dea-Acuit	114	DKK-1	100	HDAC4	250	Mag Mag	224	10385	402	10072 D0.0%4	475	TRUMER
43	beta-i lubulin	115	DKK-Z	18/	HHEX	259	мвр	331	PDAP1	403	ROCK1	4/5	TRIM63
44	BMX	116	DOCK1	188	HIBADH	260	MCHR1	332	PDCD5	404	RPL10A	476	Troponin T
45	BNI P3L	117	DOT1L	189	HIF-2 alpha	261	M-CSF	333	PDCD6	405	RPL11	477	TRP14
46	BOK	118	DRAK2	190	Histamine H3 R	262	MDGA2	334	PDHX	406	RPL22	478	TRPV1
47	Brevican	119	Draxin	191	Histone H1 3	263	MDH1	335	PDK-1	407	RPLP2	479	TRXR1
40	CA14	120	DSC2	102	Histone H2AV	262	MDM2	225	PDV 1	109	DD011	190	Truccin 2
40	Cadacia 45	120	DVDV4A	102	LIM CD4	204	MENT	330	PD744	400	DDC10	404	Taxasi- D
49	Caonerin-15	121	DIKK1A	193	HIVI GB1	265	IVIEK1	55/	PUZK1	409	KP519	481	Trypsin Pan
50	Cadherin-8	122	Dystroglycan	194	HMGN2	266	MEKK2	338	Perilipin-1	410	RPS25	482	TSC22
51	CALD1	123	EDN	195	HMOX1	267	MESDC2	339	PGAM2	411	RPS4X	483	TSH
52	Calretinin	124	EFEMP2	196	HN1	268	Metallothionein	340	PGK1	412	RPS 6	484	TXNDC4
53	CaM Kinasell	125	EGLN1	197	hnRNP G	269	mGluR1	341	PGLS	413	RRAS2	485	UBASH3B
54	CaMKK aloba	125	FIERD	192	hnRNPII	270	mGLu82/3	342	PGM 1	414	RSK1	486	LIBE2N
57	Careford Careford	120	ELAV/11	100	HOMERI	270	moranz/s	242	DCDD S	415	DEKO	407	10,000
55	capo	127	EDAVLI	133	HUMEK1	2/1	maluks	545	rukr-o	415	naK2	46/	UUCKB
56	CART	128	Endoglin	200	HP1BP3	272	MIB1	344	PIK3R1	416	RTN1-A	488	UROD
57	Cathepsin G	129	Endophilin A1	201	HPRT	273	MIOS	345	PIWIL2	417	RYK	489	VAP-B
58	Caveolin-1	130	Endorepellin	202	HS6ST3	274	MIP-1 beta	346	PKA RI beta	418	SC35	490	VE-Cadherin
59	CBP	131	ENSA	203	HSP90B1	275	MKP-3	347	PKC beta 1	419	SCGB3A1	491	Versican
60	CCB11	122	EnCAM	204	HSPA2	276	MIKA	349	PKCgamma	420	SCGE	492	Vimentin
60	00000	132	Eponin Fatter	207	100.02	270	MANA	240	nvio	424	000	402	ALL
61	CCR2	133	EphA3	205	HSPB8	2//	MN1	349	PKLR	421	SEC13	493	WNK1
62	CCR6	134	Ephrin-B3	206	IBP160	278	MPP5	350	PKN2	422	SECISBP2	494	WNK2
63	CCR8	135	ERBB4	207	ICAM-5	279	M-Ras	351	PLA2G2A	423	SEMA3F	495	WT1
64	CCR9	136	ERK1	208	IKK al pha	280	MSH6	352	PLC-beta 4	424	SENP8	496	W WOX
65	CD106	137	ERK2	209	IKK gamma	281	Musashi-1	353	PLC-gamma 1	425	Serpin A12	497	XPB
66	CD161	138	FRK4	210	II-12 R beta 2	282	MyD88	354	Plexin A1	426	Secolo A3N	498	YAP1
67	00164	129	EAIAO	211	11.175	292	MVHC	255	Playin A2	427	Samin AF	499	V
6/	00164	159	FAINS	211	16-1/1	203	IVITHC .	222	Flexin Az	427	Serpin A6		105
68	CD19	140	FANCD2	212	IL17RA	284	Myoglobin	356	Plexin A3	428	serpin D1	500	ZB184
69	CD28	141	Fascin	213	IL-20RB	285	NAP1L1	357	Plexin B3	429	SerpinE2		
70	CD29	142	FASN	214	IL-21R	286	Nbs1	358	PLO D2	430	SerRS		
71	CD36	143	FBPase1	215	IL-6R	287	NCAM2	359	PLTP	431	SGSH		
72	CD39L4	144	ECGR2B	216	IMPA1	288	NCOR1	360	PNPIA2	432	SHC1		
						~~~							i

## VII. Interpretation of Results:

#### A. Explanation of Controls Spots

To obtain optimal results using a chemiluminescence imaging system (UVP Biolmaging Systems), it is suggested to try several different exposure times until the best one is determined. Then, by comparing the signal intensities, relative expression levels of the target proteins can be made. The intensities of signals can be quantified by densitometry. There are three Positive Controls (POS1, POS2, POS3) in each array. These are three levels of standardized anti-HRP antibodies, which will produce positive control signals after incubation with HRP-conjugated Streptavidin. With all other variables being equal, the Positive Control intensities will be the same for each subarray, which allows for inter-array normalization. Antibody affinity to its target varies significantly between antibodies. The intensity detected on the array with each antibody depends on this affinity; therefore, signal intensity comparison can be performed only within the same antibody/antigen system and not between different antibodies. Some arrays may have beta-actin and GAPDH as internal controls, much as "housekeeping" genes or proteins are used to normalize results in PCR or Western blots, respectively.

#### **B.** Typical Results

The following figure shows the typical result of arrays probed with rat sample.



RayBio[®] L-Series Membrane Protocol – Rat Antibody Array L-1500

Note: In the absence of an external standard curve for each protein detected, there is no means of assessing absolute or relative concentrations of different proteins in the same sample using immunoassays. If you wish to obtain quantitative data (ie, concentrations of the various analytes in your samples), try using our Quantibody® Arrays as a targeted follow up experiment.

### C. Background Subtraction

Once you have obtained densitometry data, it is recommended to subtract the local background and normalize to the Positive Control signals before proceeding to analysis.

#### D. Normalization of Array Data

To normalize signal intensity data, one sub-array is defined as "reference" to which the other arrays are normalized. This choice is arbitrary. For example, in our Analysis Tool Software (described below), the array represented by data entered in the left-most column each worksheet is the default "reference array."

You can calculate the normalized values as follows:

X(Ny) = X(y) * P1/P(y)

Where:

P1 = mean signal intensity of POS spots on reference array

P(y) = mean signal intensity of POS spots on Array "y"

X(y) = mean signal intensity for spot "X" on Array "y"

X(Ny) = normalized signal intensity for spot "X" on Array "y"

The RayBio[®] Analysis Tool software is available for use with data obtained using RayBio[®] Biotin Label-based Antibody Arrays. You can copy and paste your signal intensity data (with and without background) into the Analysis Tool, and it will automatically normalize signal intensities to the Positive Controls.

## E. Threshold of Significant Difference

After subtracting background signals and normalization to Positive Controls, comparison of signal intensities between and among array images can be used to determine relative differences in expression levels of each protein between samples or groups.

Any  $\geq$ 1.5-fold increase or  $\leq$ 0.65-fold decrease in signal intensity for a single analyte between samples or groups may be considered a measurable and significant difference in expression, provided that both sets of signals are well above background (Mean background + 2 standard deviations, accuracy  $\approx$ 95%).

#### F. Pathway Analysis of the Array Proteins

Rat antibody array L-1500 detects 1500 unique rat proteins, including most analyzed cytokines, chemokines, adipokines, extracellular matrix proteins, growth factors, angiogenic factors, proteases, enzymes, soluble and transmembrane receptors and transport proteins, adhesion molecules and other proteins. All the array proteins are provided with their Uniprot number and GeneID, which are essential for further data mining. Raybiotech offers affordable biostatistics and bioinformatics service, including data clean-up, differential expression analysis, cluster analysis, biomarker selection, pathway analysis and experimental design. See more details on the website: https://www.raybiotech.com/biostatistics-and-bioinformatics-services

## VIII. Troubleshooting Guide

Problem	Cause	Recommendation					
	Taking too much time for detection	The whole detection process must be completed within 30 min.					
	Film developer does not work properly.	Fix film developer.					
	Did not mix HRP-Streptavidin well before use.	Mix tube containing HRP-Conjugated Streptavidin well before use since precipitates may form during storage.					
	Sample is too diluted	Increase sample concentration					
Weak Signal	Labeling reagent does not function well	Labeling reagent needs to be saved in - 20C and avoid free thaw cycle. Always use fresh labeling reagent for sample labelling.					
	Other	Check if there were any contamination with any solution containing amines in biotin-labeling step.					
		Slightly increase HRP concentrations.					
		Work as quickly as possible after mix Detection Buffer C and D.					
	Bubble formed during incubation	Remove bubbles during incubation.					
Uneven signal	Membranes were not completely covered with solution	Completely cover membranes with solution.					
	Insufficient wash	Use more stringent wash.					
	Exposure time is too long	Decrease exposure time.					
High background	Membranes dry out during experiment.	Completely cover membranes with solution during experiment. Cover tray with lid.					
	Sample is too concentrated.	Dilute sample.					

## **IX. Selected References**

Christina Scheel et all., *Paracrine and Autocrine Signals Induce and Maintain Mesenchymal and Stem Cell States in the Breast*. Cell. 2011;145, 926–940

Lin Y, Huang R, Chen L, et al., *Profiling of cytokine expression by biotinlabeled-based protein arrays*. Proteomics. 2003, 3: 1750–1757.

Huang R, Jiang W, Yang J, et al., *A Biotin Label-based Antibody Array for High-content Profiling of Protein Expression*. Cancer Genomics Proteomics. 2010; 7(3):129–141.

Liu T, Xue R, Dong L, et al., *Rapid determination of serological cytokine biomarkers for hepatitis B–virus-related hepatocellulare carcinoma using antibody arrays*. Acta Biochim Biophys Sin. 2011; 43(1):45–51.

Cui J, Chen Y, Chou W-C, et al., *An integrated transcriptomic and computational analysis for biomarker identification in gastric cancer*. Nucl Acids Res. 2011; 39(4):1197–1207.

Jun Zhong et all., *Temporal Profiling of the Secretome during Adipogenesis in Humans*. Journal of Proteome Research. 2010, 9, 5228–5238

Chowdury UR, Madden BJ, Charlesworth MC, Fautsch MP., *Proteomic Analysis of Human Aqueous Humor*. Invest Ophthalmol Visual Sci. 2010; 51(10):4921–4931.

Wei Y, Cui C, Lainscak M, et al., *Type-specific dysregulation of matrix metalloproteinases and their tissue inhibitors in end-stage heart failure patients: relationshp between MMP-10 andLV remodeling*. J Cell Mol Med. 2011; 15(4):773–782.

Kuranda K, Berthon C, Lepêtre F, et al., *Expression of CD34 in hematopoietic cancer cell lines reflects tightly regulated stem/progenitor-like state*. J Cell Biochem. 2011; 112(5):1277–1285.

Toh HC, Wang W-W, Chia WK, et al., *Clinical Benefit of Allogenic Melanoma Cell Lysate-Pulsed Autologous Dendritic Cell Vaccine in MAGE-Positive Colorectal Cancer Patients*. Clin Chem Res. 2009; 15:7726–7736. Zhen Hou, Cytokine array analysis of peritoneal fluid between women with endometriosis of different stages and those without endometriosi. Biomarkers. 2009;14(8): 604-618.

Yao Liang Tang, et al., *Hypoxic Preconditioning Enhances the Benefit of Cardiac Progenitor Cell Therapy for Treatment of Myocardial Infarction by Inducing CXCR4*. Circ Res. 2009;109:197723

RayBio[®] L-series Antibody Arrays are patent-pending technology developed by RayBiotech.

This product is intended for research only and is not to be used for clinical diagnosis. Our produces 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 six months from the date of shipment when handled and stored properly. In the event of any defect in quality or merchantability, RayBiotech's liability to buyer for any claim relating to products shall be limited to replacement or refund of the purchase price.

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

This product is for research use only.



©2021 RayBiotech, Inc.