RayBio[®] Label-Based (L-Series) Human Cancer screening array

Patent Pending Technology User Manual (Apr 14, 2023)

For the simultaneous detection of the relative expression of 500 Human proteins in serum, plasma, cell culture supernatants, cell/tissue lysates or other body fluids.

AAH-BLG-CAN-4 (4 Sample Kit) AAH-BLG-CAN-8 (8 Sample Kit)

Please read manual carefully before starting experiment



Your Provider of Excellent Protein Array Systems and Services

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

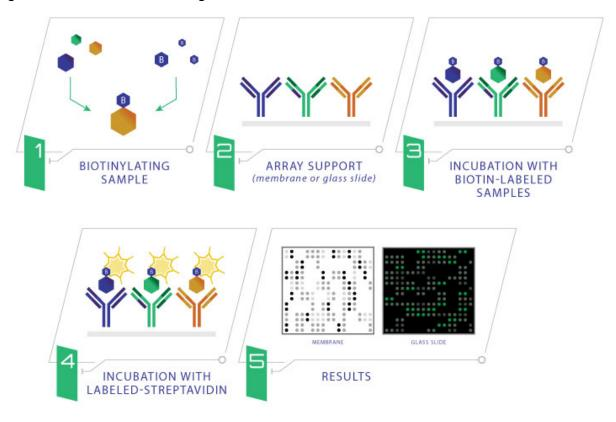
Table of Contents

I.	Introduction	3
II.	Materials Provided	4
	A. Storage Recommendations	4
	B. Additional Materials Required	4
III.	Overview and General Considerations	5
	A. Preparation and Storage of Samples	5
	B. Handling the Glass Slides	6
	C. Layout of Array Slide	7
	D. Incubations and Washes	7
IV.	Protocol	8
	A. Sample Purification	8
	B. Biotin Labeling the Sample	9
	C. Drying the Glass Slide	9
	D. Blocking and Incubations	10
	E. Fluorescence Detection	12
V.	Antibody Array Map	13
VI.	Antibody Array Target List	14
VII.	Interpretation of Results	15
	A. Explanation of Controls Spots	15
	B. Typical Results	15
	C. Background Subtraction	15
	D. Normalization of Array Data	16
	E. Threshold of Significant Difference	20
VIII.	Troubleshooting Guide	21
IX.	Selected References	22

I. Introduction

Combining direct antigen-labeling technology with our vast library of array-validated 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 glass slide arrays are then blocked, just like a western blot, and the biotin-labeled sample is added onto the glass slide, which is pre-printed with capture antibodies. The slide is incubated to allow binding of target proteins. Streptavidin-conjugated fluorescent dye (Cy3 equivalent) is then applied to the array. Finally, the glass slide is dried, and laser fluorescence scanning is used to visualize the signals.



II. Materials Provided

A. Storage Recommendations

Upon receipt, the kit should be stored at -20°C until needed. It is recommended to use the kit within 6 months of the date of shipment. After initial use, remaining reagents should be stored at 4°C and may be stored for up to 3 months. Labeling Reagent (Item B) should be prepared fresh each time before use. Unused glass slides should be kept at -20°C and repeated freeze-thaw cycles should be avoided (slides may be stored for 6 months).

ITEM	DESCRIPTION	4 SAMPLE KIT	8 SAMPLE KIT
А	Spin Columns (0.5ml)	8 columns	16 columns
В	Labeling Reagent	1 vial	2 vials
D	Stop Solution	1 vial (50 µl)	1 vial (50 µl)
E	RayBio [®] L-Series Glass Slide*	1 slide	2 slides
F	Blocking Buffer	1 bottle (8 ml)	2 bottles (8 ml)
G	20X Wash Buffer I	1 bottle (30 ml)	1 bottle (30 ml)
Н	20X Wash Buffer II	1 bottle (30 ml)	1 bottle (30 ml)
I	Cy3 equivalent-Conjugated Streptavidin	1 vial	2 vials
J	Adhesive Plastic Strips		
К	Labeling Buffer	1 bottle (30 ml)	1 bottle (30 ml)
n/a	2X Cell Lysis Buffer**	1 bottle (10 ml)	1 bottle (10 ml)
М	30 ml Centrifuge Tube	1 tube	1 tube

*Each slide contains 4 identical subarrays

**Only needed if testing cell or tissue lysates

B. Additional Materials Required

- 1 ml tube, small plastic or glass containers
- Orbital shaker or oscillating rocker
- Pipettors, pipette tips and other common lab consumables
- Laser scanner for fluorescence detection
- Aluminum foil

III. Overview and General Considerations

A. Preparation and Storage of Samples

- 1. Preparation of Cell Culture Supernatants
 - 1. Seed cells at a density of 1×10^6 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.**,+ The membrane-based array is recommended if high serum medium such as 10% FCS/FBS is used, as high background can occur on glass slide arrays with high serum containing media samples.
 - To collect supernatants, centrifuge at 1,000 x g for 10 minutes and store as less than or equal 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 fluorescent 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 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 minutes.
 - 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 2x10⁷ cells/ml in 1X Cell Lysis Buffer.

3. 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 minutes 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. Determining 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 approximate 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 minutes 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

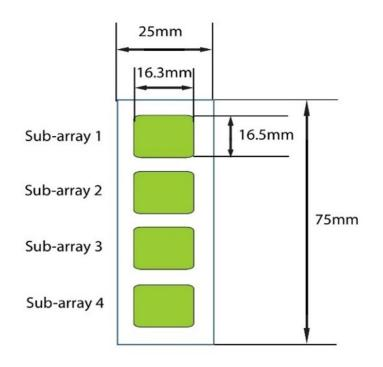
For optimal biotin labeling, it is necessary to determine the protein concentration in the cell/tissue lysate. We recommended using a BCA total protein assay (e.g., Pierce, Catalog # 23227).

B. Handling the Glass Slides

- The microarray slides are delicate. Please do not touch the array surface with pipette tips, forceps or your fingers. Hold the slides by the edges only.
- Handle the slides with powder-free gloves and in a clean environment.
- Do not remove the glass slide from the chamber assembly until step 20, and take great care not to break the glass slide when doing so.
- Remove reagents/sample by gently applying suction with a pipette to corners of each chamber. Do not touch the printed area of the array, only the sides as seen in image below.



C. Layout of Array Slide



Four identical sub-arrays on one slide

4 printed sub-arrays per glass chip

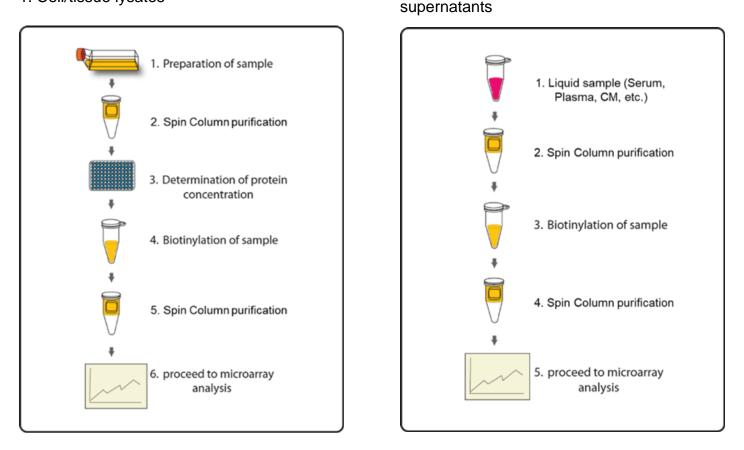
D. Incubations and Washes

- Cover incubation chamber with a Plastic Adhesive Strip (Item J) to prevent evaporation during incubation or wash steps, particularly those steps lasting 2 hours or longer.
- During incubation and wash steps avoid foaming and remove all bubbles from the sub-array surface.
- Perform all incubation and wash steps under gentle rotation or rocking motion (~0.5 to 1 cycle/sec).
- Wash steps in Wash Buffer II and all incubation steps may be performed overnight at 4°C
- Avoid cross-contamination of samples to neighboring wells. To remove Wash Buffers and other reagents from chamber wells, you may invert the Glass Slide Assembly to decant, and aspirate the remaining liquid.
- Unlike most Cy3 fluors, the streptavidin-conjugated fluor used in this kit is very stable at room temperature (RT) and resistant to photobleaching on the hybridized glass slides. However, please protect glass slides from direct, strong light and temperatures above RT.

IV. Protocol

Assay Diagram

1. Cell/tissue lysates



2. Serum, plasma, or Cell culture

A. Sample Purification

Note: This step removes low molecular weight amine derivatives or unwanted buffer from samples to ensure quality biotinylation in Steps 5-7.

- 1. Twist to remove the bottom plug of the Spin Column and loosen the cap (do not remove).
- 2. Place the Spin Column into a collection tube and centrifuge at 1,500 x g for 1 minute to remove the storage buffer. Discard the flow-through.
- 3. Wash the Spin Column three times with 300 µl Labeling Buffer each, centrifuge at 1,500 x g for 1 minute to remove the flow-through. Discard the flow-through and blot the bottom of the column to remove excess liquid. Transfer the Spin Column to a new collection tube.
- 4. Apply sample on top of the resin within the next few minutes. Centrifuge at 1,500 x g for 2 minutes. Collect the flow-through that contains the sample. The recommended sample dilutions are as follows:

- o Cell culture supernatant: 120 μl neat supernatant
- o Serum/Plasma: 2 μl serum/plasma in 100 μl Labeling Buffer
- o Cell/tissue lysate: 20 μg lysate in 100 μl Labeling Buffer

Note: Each labelled sample volume is enough for at least 3 arrays following the protocol below.

Note: The maximal sample volume is 130 μ l for each Spin Column. Do not load over 130 μ l 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 the Labeling Reagent. Briefly spin down the Labeling Reagent tube (Item B). Add 100 µI 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 8 µl of Labeling Reagent into the sample tube (for 120 µl supernatant).
 - b. For labeling serum or plasma: Add 8 µl of Labeling Reagent into the sample tube (for 2 µl serum/plasma *in 100 µl labeling buffer*).
 - c. For labeling cell or tissue lysates: Add 4 µl of 1X Labeling Reagent into the sample tube (for 20 µg lysate *in 100 µl labeling buffer*).
 - d. For all other body fluid: Add 2 μ l of Labeling Reagent Solution per 100 μ g sample to be labelled.

Note: The addition of Labeling Reagent volume is based upon the sample amount used in Step 4. If the amount of sample being labelled differs from the example in Step 6, adjust this volume proportionally.

Add 3 µl Stop Solution (Item D) to each sample 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. Drying the Glass Slide

8. Remove the package containing the Assembled Glass Slide (Item E) from the freezer. Place unopened package on the bench top for ~15 minutes, and allow the Assembled Glass Slide to equilibrate to RT.

 Open package, and take the Assembled Glass Slide out of the sleeve. Do <u>not</u> disassemble the Glass Slide from the chamber assembly. Place glass slide assembly in laminar flow hood or similar clean environment for 1-2 hours at RT.

Note: Protect the slide from dust or other contaminants.

D. Blocking and Incubations

Note: Glass slide should be completely dry before adding Blocking Buffer to wells.

- 10. Block sub-arrays by adding 400 μl of Blocking Buffer (Item F) into each well of Assembled Glass Slide and incubating at RT for 30 minutes. Ensure there are no bubbles on the array surfaces.
- 11. Dilute samples with Blocking Buffer. Recommended dilution of the biotin-labeled samples with Blocking Buffer is 10-fold for cell culture supernatants, 20-fold for serum/plasma and 100-fold for cell/tissue lysate. *Dilution for other body fluid needs to be determined by the end user. Generally, most samples can be 10-20x dilution, while tears and saliva samples may need 100x dilution.*

Note: Optimal sample dilution factor will depend on the abundance of target proteins. If the background or antigen-specific antibody signals are too strong, the sample can be diluted further in subsequent experiments. If the signal is too weak, more concentrated samples can be used.

12. Completely remove the Blocking Buffer from each well. Add 400 µl of diluted sample into appropriate wells. Remove any bubbles on array surfaces. Incubate arrays with gentle rocking or shaking for 2 hours at RT or overnight at 4°C

Note: Avoid the flow of sample into neighboring wells.

- 13. Based on number of samples and remaining protocol, calculate the amount of 1X Wash Buffer I and 1X Wash Buffer II needed to complete the experiment. Separately dilute the required amounts of 20X Wash Buffer I Concentrate (Item G) 20-fold and 20X Wash Buffer II Concentrate (Item H) with ddH₂O
- 14. Decant the samples from each well and wash 3 times with 800 μl of 1X Wash Buffer I at RT with gentle rocking or shaking for 5 minutes per wash.
- 15. Obtain a clean container (e.g., pipette tip box or slide-staining jar), place the Assembled Glass Slide into the container with enough volume of 1X Wash Buffer I to completely cover the entire assembly, and remove any bubbles in wells. Wash 2 times at RT with gentle rocking or shaking for 10 minutes per wash.
- 16. Decant the Wash Buffer I from each well, place the Assembled Glass Slide into the container with enough volume of 1X Wash Buffer II to completely cover the entire assembly, and remove any bubbles in wells. Wash 2 times at RT with gentle rocking or shaking for 5 minutes per wash.
- 17. Prepare 1X Cy3-Conjugated Streptavidin:
 - a. Briefly spin down tube containing the Cy3-Conjugated Streptavidin (Item I) immediately before use.

- b. Add 1000 µl of Blocking Buffer into the Cy3-Conjugated Streptavidin tube to prepare a concentrated Cy3-Conjugated Streptavidin stock solution. Pipette up and down to mix gently (do not store the stock solution for later use).
- c. To prepare 1X Cy3-Conjugated Streptavidin, add 200 µl of the concentrated Cy3-Conjugated Streptavidin stock solution into a tube with 800 µl of Blocking Buffer. Mix gently.
- Carefully remove Assembled Glass Slide from container. Remove all of Wash Buffer II from the wells. Add 400 µl of 1X Cy3-Conjugated Streptavidin to each sub-array. Cover the incubation chamber with the plastic adhesive strips.

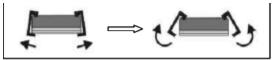
Note: Avoid exposure to light in Steps 19-25 by covering the Glass Slide Assembly with aluminum foil or incubate in a dark room.

19. Incubate with 1X Cy3-Conjugated Streptavidin at RT for 1 hour with gentle rocking or shaking.

Note: Incubation may be done overnight at 4°C

20. Decant the solution and disassemble the glass slide from the incubation frame and chamber. Disassemble the device by pushing clips outward from the side, as shown below. Carefully remove the glass slide from the gasket.

Note: Be careful not to touch the printed surface of the glass slide, which is on the same side as the barcode.



- 21. Gently place the glass slide into 30 ml Centrifuge Tube (Item M). Add enough 1X Wash Buffer I to cover the entire glass slide (about 30 ml). Wash with gentle rocking or shaking for 10 min. Remove the wash buffer. Repeat 2 times for a total of 3 washes.
- 22. Add enough 1X Wash Buffer II to cover the entire glass slide (about 30 ml). Wash with gentle rocking or shaking for 5 minutes. Remove the wash buffer. Repeat one time for a total of two washes for 5 minutes per wash.
- 23. Finally, wash the glass slide with 30 ml of ddH₂O for 5 minutes. Remove glass slide and decant water from Centrifuge Tube.
- 24. Remove buffer droplets from the slide completely by one of the following ways:
 - Put the glass slide into the Slide Washer/Dryer, and dry the glass slide by centrifuge at 1,000 rpm for 3 minutes without cap.
 - $\circ\,$ Or dry the glass slide by a compressed N2 stream.
 - Or gently apply suction with a pipette to remove buffer droplets. Do not touch the array surface, only the sides.

Note: Make sure the finished glass slide is completely dry before scanning or storage.

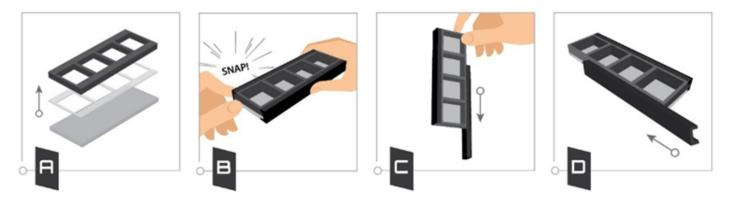
E. Fluorescence Detection

25. You may proceed immediately to scanning or you may store the slide at -20°C in the Centrifuge Tube provided or at RT to scan at a later time.

Note: <u>Please protect the finished glass slides from temperatures above RT and store them in the</u> <u>dark.</u> Do not expose glass slide to strong light, such as sunlight or a UV lamp.

Note: If you need to repeat any of the incubation steps after finishing the experiment, you must first re-assemble the glass slide into the incubation chamber by following the steps as described below. To avoid breaking the printed glass slide, you may first want to practice assembling the device with a blank glass slide.

- 1. Apply slide to incubation chamber barcode facing upward (image A).
- 2. Gently snap one edge of a snap-on side (image B).
- 3. Gently press other of side against lab bench and push in lengthwise direction (image C).
- 4. Repeat with the other side (image D)



V. Antibody Array Map

	1	2	3	4	5	6	7	8	9	10	11	12	13	14	15	16	17	18	19	20	21	22	23	24	25	26	27	28	29	30
1	POS1	POS1	POS2	POS2	POS3	POS3	Neg	Neg	1	1	2	2	3	3	4	4	5	5	6	6	7	7	8	8	9	9	10	10	11	11
2	12	12	13	13	14	14	15	15	16	16	17	17	18	18	19	19	20	20	21	21	22	22	23	23	24	24	25	25	26	26
3	27	27	28	28	29	29	30	30	31	31	32	32	33	33	34	34	35	35	36	36	37	37	38	38	39	39	40	40	41	41
4	42	42	43	43	44	44	45	45	46	46	47	47	48	48	49	49	50	50	51	51	52	52	53	53	54	54	55	55	56	56
5	57	57	58	58	59	59	60	60	61	61	62	62	63	63	64	64	65	65	66	66	67	67	68	68	69	69	70	70	71	71
6	72	72	73	73	74	74	75	75	76	76	77	77	78	78	79	79	80	80	81	81	82	82	83	83	84	84	85	85	86	86
7	87	87	88	88	89	89	90	90	91	91	92	92	93	93	94	94	95	95	96	96	97	97	98	98	99	99	100	100	101	101
8	102	102	103	103	104	104	105	105	106	106	107	107	108	108	109	109	110	110	111	111	112	112	113	113	114	114	115	115	116	116
9	117	117	118	118	119	119	120	120	121	121	122	122	123	123	124	124	125	125	126	126	127	127	128	128	129	129	130	130	131	131
10	132	132	133	133	134	134	135	135	136	136	137	137	138	138	139	139	140	140	141	141	142	142	143	143	144	144	145	145	146	146
11	147	147	148	148	149	149	150	150	151	151	152	152	153	153	154	154	155	155	156	156	157	157	158	158	159	159	160	160	161	161
12	162	162	163	163	164	164	165	165	166	166	167	167	168	168	169	169	170	170	171	171	172	172	173	173	174	174	175	175	176	176
13	177	177	178	178	179	179	180	180	181	181	182	182	183	183	184	184	185	185	186	186	187	187	188	188	189	189	190	190	191	191
14	192	192	193	193	194	194	195	195	196	196	197	197	198	198	199	199	200	200	201	201	202	202	203	203	204	204	205	205	206	206
15	207	207	208	208	209	209	210	210	211	211	212	212	213	213	214	214	215	215	216	216	217	217	218	218	219	219	220	220	221	221
16	222	222	223	223	224	224	225	225	226	226	227	227	228	228	229	229	230	230	231	231	232	232	233	233	234	234	235	235	236	236
17	237	237	238	238	239	239	240	240	241	241	242	242	243	243	244	244	245	245	246	246	247	247	248	248	249	249	250	250	251	251
18	252	252	253	253	254	254	255	255	256	256	257	257	258	258	259	259	260	260	261	261	262	262	263	263	264	264	265	265	266	266
19	267	267	268	268	269	269	270	270	271	271	272	272	273	273	274	274	275	275	276	276	277	277	278	278	279	279	280	280	281	281
20	POS1	POS1	POS2	POS2	POS3	POS3	Neg	Neg	282	282	283	283	284	284	285	285	286	286	287	287	288	288	289	289	290	290	291	291	292	292
21	293	293	294	294	295	295	296	296	297	297	298	298	299	299	300	300	301	301	302	302	303	303	304	304	305	305	306	306	307	307
22	308	308	309	309	310	310	311	311	312	312	313	313	314	314	315	315	316	316	317	317	318	318	319	319	320	320	321	321	322	322
23	323	323	324	324	325	325	326	326	327	327	328	328	329	329	330	330	331	331	332	332	333	333	334	334	335	335	336	336	337	337
24	338	338	339	339	340	340	341	341	342	342	343	343	344	344	345	345	346	346	347	347	348	348	349	349	350	350	351	351	352	352
25	353	353	354	354	355	355	356	356	357	357	358	358	359	359	360	360	361	361	362	362	363	363	364	364	365	365	366	366	367	367
26	368	368	369	369	370	370	371	371	372	372	373	373	374	374	375	375	376	376	377	377	378	378	379	379	380	380	381	381	382	382
27	383	383 398	384	384	385	385 400	386	386 401	387 402	387	388 403	388 403	389 404	389 404	390	390 405	391 406	391 406	392 407	392 407	393 408	393 408	394 409	394 409	395	395 410	396 411	396 411	397	397 412
28	398 413	413	399 414	399 414	400	400	401	401	402	402	403	403	404	404	405	405	406	406	407	407	408	408	409	409	410	410	411	411	412	412
29	415	415	414	414	415	415	410	410	417	417	418	418	419	419	420	420	421	421	422	422	425	425	424	424	425	440	420	420	442	427
30	443	428	429	429	450	450	451	451	452	452	455	455	434	454	450	450	450	450	457	457	458	458	459	459	440	440	441	441	442	442
31	443	443	459	459	445	445	440	440	447	447	448	440	449	449	465	450	466	466	452	452	455	455	454	454	433	433	436	430	472	437
32	400	456	439	439	400	400	401	401	402	402	405	405	404	404	465	465	400	400	482	482	408	408	484	484	470	470	471	471	472	472
33	488	473	489	474	490	475	470	491	492	492	493	478	494	494	495	495	496	496	497	497	498	498	499	494	500	500	Neg	Neg	Neg	Neg
34	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	Neg	POS3	POS3	POS2	POS2	POS1	POS1
35	HCB	INCE	Heg	HICE	neg	нев	Neg	Heg	NCE	Heg	HACE	Neg	HICE	INCE	HCB	INCE	HCB	HICE	HCE	HICE	INCE	HEE	INCE	Heg	2033	2033	2032	2032	.031	.031

VI. Antibody Array Target List

Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name
1	ABL1	73	CDK4	145	FGF19	217	IGF2	289	MMP9	361	Prostasin	433	SULT2A1
2	ACTN1	74	CDK6	146	FGF2	218	IGFBP2	290	MPO	362	PSA	434	SYK
3	AFP	75	CDKN1A	147	FGF20	219	IGFBP3	291	MRAS	363	PTCH1	435	TCF3
4	AKT1	76	CDKN1B	148	FGF21	220	IKBKB	292	MSH2	364	PTCH2	436	TCF7
5	AKT2	77	CDKN2A	149	FGF22	221	IKBKG	293	MSLN	365	PTEN	437	TCF7L1
6	AKT3	78	CDKN2B	150	FGF23	222	IL12B	294	MYC	366	PTGS2	438	TGFA
7	ALDH3A1	79	CEA	151	FGF3	223	IL1R2	295	NANOG	367	PTK2	439	TGFB1
8	ALK	80	CEBPA	152	FGF4	224	IL-2	296	NCOA3	368	PTPN11	440	TGFB2
9	APC	81	CEBPE	153	FGF5	225	IL2RB	297	NCOR1	369	PTPN6	441	TGFB3
10	ARAF	82	CHGA	154	FGF6	226	IL3	298	NFKB1	370	PXN	442	TGFBR1
11	ARHGEF1	83	CHUK	155	FGF7	227	IL6ST	299	NFKB2	371	RAC1	443	TGFBR2
12	ARHGEF6	84	CKMB	156	FGF8	228	IL-8	300	NFKBIA	372	RAC3	444	THBS1
13	ARNT	85	COL1A1	157	FGF9	229	ILK	301	NGFR	373	RAD51	445	TIMP1
14	ATF1 ATF2	86 87	COL4A1 CREB1	158 159	FGFR1 FGFR2	230 231	INSRR IRF3	302 303	NKX3-1 NOS2	374	RAF1 RALA	446 447	TIMP3 TLR2
15	ATF4	88	CREB1	160	FGFR2	231	IRS1	304	NOTCH1	375	RALA	447	TLR2
10	ATM	89	CREB3L2	161	FHIT	232	ITGA2	305	NOTCH2	377	RANK	449	TMPRSS2
18	AXIN1	90	CREBBP	162	FLT1	233	ITGA2	306	NOTCH3	378	RAP1A	450	TNC
19	AXIN2	91	CRK	163	FLT3	235	ITGA5	307	NOTCH4	379	RARA	451	TNFSF11
20	B2M	92	CRKL	164	FLT4	236	ITGA6	308	NR4A3	380	RARB	452	TNR
21	B7-H4	93	CSF1R	165	FN1	237	ITGAM	309	NRAS	381	RASSF1	453	TNXB
22	BAD	94	CSF2	166	FOS	238	ITGAV	310	NSE	382	RB1	454	TP53
23	BAK1	95	CTNNB1	167	FOXO1	239	ITGB1	311	NTRK1	383	RBPJ	455	TP63
24	BAX	96	CYCS	168	FOXP1	240	ITGB3	312	NTRK3	384	RECK	456	TRADD
25	BCL2	97	CYFRA21-1	169	FUT8	241	ITGB5	313	OPN	385	REL	457	TRAF1
26	BCL2L1	98	CYP1B1	170	FZD1	242	ITGB7	314	PAK1	386	RELA	458	TRAF2
27	BCL2L11	99	DDB2	171	FZD2	243	JAG1	315	PAK2	387	RET	459	TRAF3
28	BCL6	100	DDX5	172	FZD3	244	JAG2	316	PAK3	388	RHEB	460	TRAF4
29	BCR	101	DGKB	173	FZD4	245	JAK1	317	PAK4	389	RHOA	461	TRAF5
30	BIRC2	102	DGKD	174	FZD5	246	JAK3	318	PAK6	390	ROCK1	462	TRAF6
31	BIRC3	103	DGKE	175	FZD6	247	JUN	319	PAX3	391	ROCK2	463	TRIM71
32	BMI1	104	DGKG	176	FZD7	248	JUP	320	PAX5	392	RPS6	464	TSC1
33 34	BMP2 BMP4	105 106	DGKI	177 178	FZD8 GADD45A	249 250	KDR KIAA1303	321 322	PAX7 PAX8	393 394	RPS6KA5 RPS6KB1	465 466	TSC2
35	BMPR2	108	DGKQ DGKZ	178	GADD45A GADD45B	250	KLK10	323	PAX6 PDCD4	395	RRAS2	466	TWIST1 TWIST2
36	BRAF	107	DULL1	1/5	GADD456	252	KLK10	323	PDGFA	396	RUNX1	468	TYMP
37	BRCA1	100	DLL3	181	GCK	253	KRAS	325	PDGFB	397	RUNX2	469	UGT1A1
38	BRCA2	110	DLL4	182	GLI1	254	LAMA1	326	PDGFC	398	RXRA	470	VAV1
39	C3	111	DUSP6	183	GLI2	255	LAMA3	327	PDGFD	399	RXRB	471	VEGFA
40	CA125	112	DVL1	184	GLI3	256	LAMA4	328	PDGFRA	400	RXRG	472	WAS
41	CA15-3	113	DVL2	185	GP73	257	LDHA	329	PDGFRB	401	SCCA1	473	WASF1
42	CA19-9	114	DVL3	186	GPC1	258	LDHB	330	PDPK1	402	SCF	474	WASF3
43	CA242	115	E2F1	187	GPC3	259	LEF1	331	PFKM	403	SDC1	475	WNT1
44	CA72-4	116	E2F2	188	GRB2	260	Leptin	332	PGA3	404	SDC2	476	WNT10A
45	CAMK2A	117	EFNA3	189	GSK3B	261	LMO2	333	PGR	405	SDC4	477	WNT10B
46	CAMK2B	118	EFNA4	190	GSN	262	LRP5	334	PIK3CA	406	SERPINB5	478	WNT11
47	CAMK2D	119	EFNA5	191	GSTP1	263	LRP6	335	PIK3CB	407	SHC1	479	WNT16
48	CASP3	120	EGF	192	GZMB	264	LTBR	336	PIK3CD	408	SHC2	480	WNT2
49	CASP8	121	EGFR	193	HBEGF	265	LUM	337	PIK3R1	409	SHH	481	WNT2B
50 51	CASP9 CAV1	122 123	EGLN1 EGLN2	194 195	HCGb HDAC1	266 267	MAD1L1 MAF	338 339	PIK3R2 PIK3R3	410	SIN3A SIRT1	482 483	WNT3A WNT4
51	CAV1 CAV2	123	EGLN2 EGLN3	195	HDAC1 HDAC2	267	MAP2K1	339	PIK3R3 PIM1	411 412	SIRT1 SIRT3	483	WNT5A
53	CAV2 CAV3	124	EGLINS EIF2AK2	196	HDAC2	268	MAP2K1 MAP2K2	340	PIM1 PIM2	412	SKP2	484	WNT5A WNT5B
54	CBL	125	EIF4B	197	HE4	209	MAP2K2 MAPK1	341	PLA2G4A	413	SLC22A1	485	WNT6
55	CCNA1	120	EIF4EBP1	199	HGF	271	MAPK11	343	PLAT	415	SLC22A1	487	WNT7A
56	CCNA2	128	ELK1	200	HHEX	272	MAPK12	344	PLAU	416	SLC2A1	488	WNT7B
57	CCND1	129	EP300	201	HIF1A	273	МАРКЗ	345	PLAUR	417	SLC2A2	489	WNT8B
58	CCND2	130	EPAS1	202	HIST3H3	274	MAPK7	346	PLCG1	418	SLC7A5	490	WNT9A
59	CCND3	131	ERBB2	203	HK2	275	MAPK8	347	PLCG2	419	SMAD2	491	WNT9B
60	CCNE1	132	ERBB3	204	HMGA2	276	MAPK9	348	PLD1	420	SMAD3	492	YWHAB
61	CCNE2	133	ERBB4	205	HMOX1	277	MARCKS	349	PLD2	421	SMAD4	493	YWHAE
62	CCR3	134	EREG	206	HOXD10	278	MAX	350	PML	422	SMO	494	YWHAH
63	CCR5	135	ESR1	207	HPGD	279	MCL1	351	PPARD	423	SOS2	495	YWHAQ
64	CCR8	136	ESR2	208	HPN	280	MCSF	352	PPARG	424	SP1	496	YWHAZ
65	CD14	137	ETV5	209	HRAS	281	MDM2	353	PPP1CA	425	SPI1	497	ZBTB16
66	CD44	138	EZR	210	HSD11B1	282	MEF2C	354	PRKACA	426	SPINT1	498	ZBTB17
67	CDC25A	139	FCGR1A	211	HSP90AB1	283	MEN1	355	PRKACB	427	SPRY2	499	ZEB1
68	CDC25B	140	Ferritin	212	HSP90B1	284	MET	356	PRKACG	428	SRC	500	ZEB2
69	CDC25C	141	FGF1	213	HSPG2	285	MITE	357	PRKCA	429	ST14		
70	CDC42 CDH1	142 143	FGF10 FGF16	214 215	ID2 IGF1	286 287	MMP16 MMP2	358 359	PRKCB PRKCE	430 431	STAT3 STAT5A		
71	CDH1 CDK2	145	FGF10 FGF17	215	IGF1	287	MMP2 MMP3	360	PRKCE	431	STAT5A STAT5B	8	
12	CUNZ	144	10/1/	210	IGI'IK	200	WIVIP3	300	FRAUG	432	JIAIDD	I	1

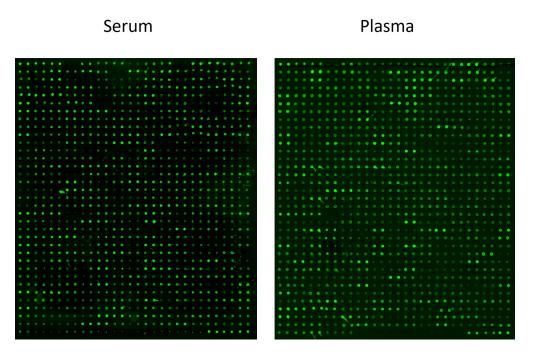
VII. Interpretation of Results

A. Explanation of Controls Spots

There are three Positive Controls (POS1, POS2, POS3) in each array. These are three levels of standardized biotinylated IgG. All other variables being equal, the Positive Control intensities will be the same for each sub-array. This allows for normalization based upon the relative fluorescence signal responses to a known control, 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 this array probed with sample(s). The images were captured using an Axon GenePix laser scanner. The Positive control signals in the upper left and lower right corners of each array can be used to identify the orientation and help normalize the results between arrays.



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 (i.e., 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 fluorescence intensity data, you should subtract the background and normalize to the Positive Control signals before proceeding to analysis.

Most laser fluorescence scanners' software has an option to automatically measure the local background around each spot. For best results, we recommend comparing signal intensities representing the MEAN signals minus local background. If your resulting fluorescence signal intensity reports do not include these values (e.g., a column labeled as "F532 Mean - B532"), you may need to subtract the background manually or change the default settings on your scanner's data report menu.

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 freely 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. Analysis Tool software can be downloaded from the product page on the RayBiotech website.

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 greater than or equal to 1.5-fold increase or less than or equal to 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 is around 95%).

VIII. 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	Ensure sufficient incubation time and change sample incubation step to overnight							
	Too low protein concentration in sample	Dilute starting sample less or concentrate sample							
	Improper storage of kit	Store kit as suggested temperature. Don't freeze/thaw the slide.							
	Bubble formed during incubation	Handle and pipette solutions more gently; De-gas solutions prior to use							
Uneven Signal	Arrays are not completed covered by reagent	Prepare more reagent and completely cover arrays with solution							
	Reagent evaporation	Cover the incubation chamber with adhesive film during incubation							
	Cross-contamination from neighboring wells	Avoid overflowing wash buffer between wells							
General	Comet tail formation	Air dry the slide for at least 1 hour before usage							
	Inadequate detection	Increase laser power so the highest standard concentration for each cytokine receives the highest possible reading yet remains unsaturated							
	Overexposure	Lower the laser power							
	Dark spots	Completely remove wash buffer in each wash step							
High	Insufficient wash	Increase wash time and use more wash buffer							
Background	Dust	Minimize dust in work environment before starting experiment							
	Slide is allowed to dry out	Take additional precautions to prevent slides from dying out during experiment							

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 biotin-labeled-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: relationship 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. GenePix[®] is a registered trademark of Molecular Devices, Inc.

This product is for research use only.



©2022 RayBiotech, Inc