RayBio[®] Label-Based (L-Series) Human Neuro disease 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-NEU-4 (4 Sample Kit) AAH-BLG-NEU-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

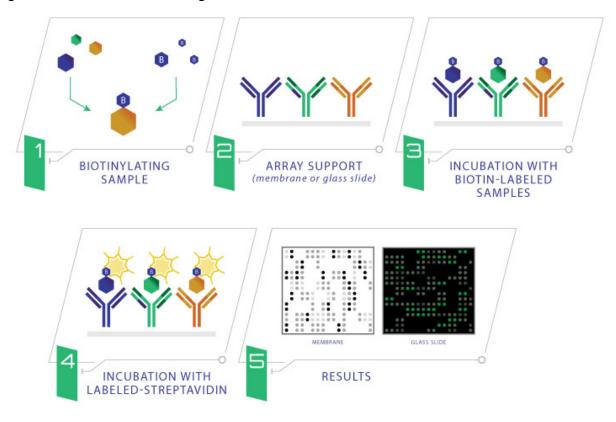
Table of Contents

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

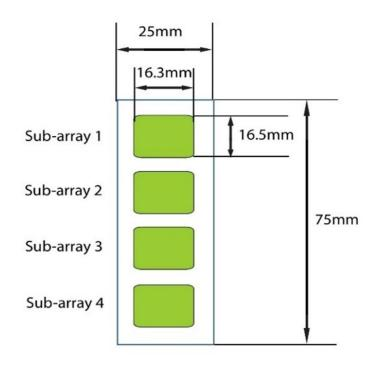
For optimal biotin labeling, it is necessary to determine the protein concentration in the cell/tissue lysate. We recommended using a BCA total protein assay (e.g., Pierce, Catalog # 23227).

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

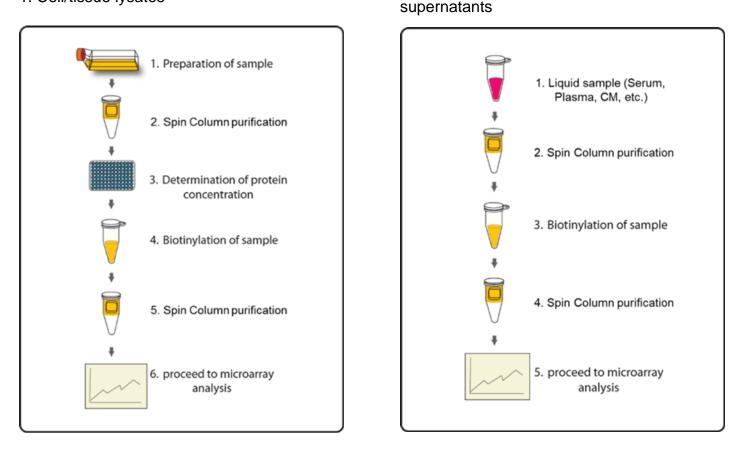
D. Incubations and Washes

- 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.

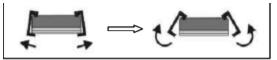
Note: Avoid exposure to light in Steps 19-25 by covering the Glass Slide Assembly with aluminum foil or incubate in a dark room.

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.

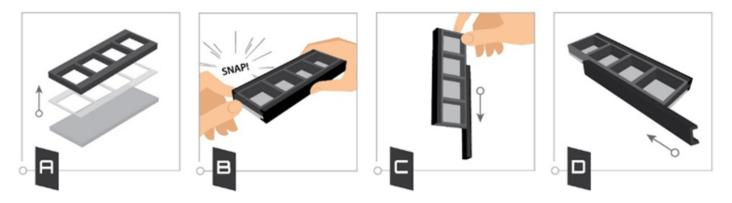
E. Fluorescence Detection

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	454	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	464	449	465	450	466	466	452	452	455	455	454	454	433	433	430	430	472	437
32	400	456	439	439	400	400	401	401	402	402	405	405	404	404	405	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	ACTR10	73	CDK5R1	145	FZD2	217	KCNJ6	289	NDUFV1	361	PSMA4	433	STX1A
2	ACTR1A	74	CHRM1	146	FZD3	218	KCNJ9	290	NDUFV2	362	PSMA5	434	TAF4
3	ACTR1B	75	CHRM3	147	FZD4	219	KIF5A	291	NEFH	363	PSMA6	435	TAF4B
4	ADAM10	76	CHRM5	148	FZD5	220	KIF5B	292	NEFL	364	PSMA7	436	TBP
5	ADAM17	77	CHRNA7	149	FZD6	221	KIF5C	293	NEFM	365	PSMA8	437	TFAM
6	ADCY1	78	CHUK	150	FZD7	222	KLC1	294	NFE2L2	366	PSMB1	438	TGM2
7	ADCY2	79	CLTB	151	GABRA1	223	KLC2	295	NFKB1	367	PSMB2	439	TH
8	ADCY3	80	CLTC	152	GABRA4	224	KLC3	296	NGF	368	PSMB3	440	TNFRSF1A
9	ADCY4	81	CLTCL1	153	GABRA5	225	KLC4	297	NOS1	369	PSMB4	441	TNFRSF1B
10	ADCY5	82	COX4I1	154	GABRB2	226	KRAS	298	NOS2	370	PSMB5	442	TOMM40
11 12	ADCY6 ADCY7	83	COX412 COX5A	155 156	GABRD GABRE	227	LPL LRP1	299 300	NOTCH1 NOX1	371 372	PSMB6 PSMB7	443 444	TOMM40L TP53
13	ADCY8	85	COX58	157	GABRG1	229	LRP1 LRP5	301	NOX1	372	PSMC2	444	TRAF2
14	ADCY9	86	COX6A1	158	GABRG2	230	LRP6	302	NRAS	374	PSMC3	446	TRPC1
15	ADORA2A	87	COX6A2	159	GABRQ	231	LRRK2	303	NRBF2	375	PSMC4	447	TUBA1A
16	ADRM1	88	COX6B1	160	GABRR2	232	MAOA	304	NRF1	376	PSMC5	448	TUBA1B
17	AKT1	89	COX7A1	161	GAPDH	233	MAOB	305	PIK3C3	377	PSMC6	449	TUBA1C
18	AKT2	90	COX7A2L	162	GLUL	234	MAP2K1	306	PIK3CA	378	PSMD1	450	TUBA3C
19	AKT3	91	COX7B	163	GNAI1	235	MAP2K2	307	PIK3CB	379	PSMD11	451	TUBA4A
20	AMBRA1	92	COX7B2	164	GNAI2	236	MAP2K3	308	PIK3CD	380	PSMD12	452	TUBA8
21	AP2A1	93	COX7C	165	GNAI3	237	MAP2K7	309	PIK3R1	381	PSMD13	453	TUBAL3
22	AP2A2	94	COX8A	166	GNAL	238	MAP3K10	310	PIK3R2	382	PSMD14	454	TUBB
23	AP2B1 AP2M1	95 96	CREB1 CREB3L1	167 168	GNAO1 GNB1	239 240	MAP3K5 MAPK1	311 312	PIK3R3 PIK3R4	383 384	PSMD2 PSMD3	455	TUBB1 TUBB2B
24	AP2IVI1 AP2S1	97	CREB3L1 CREB3L2	169	GNB1 GNB2	240	MAPK1 MAPK11	313	PINSR4 PINK1	385	PSMD3	450	TUBB3
26	APAF1	98	CSF1	170	GNB3	242	MAPK12	314	PLA2G4A	386	PSMD6	458	TUBB6
27	APBB1	99	CSNK1A1	171	GNB4	243	MAPK13	315	PLA2G4F	387	PSMD7	459	TXN
28	APH1A	100	CSNK1E	172	GNB5	244	MAPK14	316	PLCB1	388	PSMD8	460	TXN2
29	APOE	101	CSNK2A1	173	GNG11	245	MAPK3	317	PLCB2	389	PSMD9	461	UBA52
30	APP	102	CSNK2A2	174	GNG12	246	MAPK8	318	PLCB3	390	PTGS2	462	UBA7
31	ARAF	103	CTNNB1	175	GNG13	247	MAPK9	319	PLCB4	391	RAB5A	463	UBB
32	ATF4	104	CYC1	176	GNG4	248	MAPT	320	PLCG1	392	RAC1	464	UBC
33	ATF6	105	CYCS	177	GNG5	249	MFN1	321	POLR2A	393	RAF1	465	UBE2G1
34 35	ATG2A ATP2A1	106 107	DAXX DCTN1	178 179	GNG8 GNGT1	250 251	MFN2 MME	322 323	POLR2B POLR2C	394 395	RAP1A RAP1B	466	UBE2G2 UBE2J1
36	ATP2A1	107	DCTN1 DCTN2	1/9	GNGT2	252	NAE1	323	POLR2D	396	RB1CC1	468	UBE2J2
37	ATP5D	100	DCTN2 DCTN3	181	GPR37	253	NCAM1	325	POLR2E	397	RCOR1	469	UBE2L6
38	AXIN1	110	DCTN4	182	GPR83	254	NCSTN	326	POLR2F	398	RELA	470	UCHL1
39	AXIN2	111	DCTN5	183	GPX1	255	NDUFA1	327	POLR2G	399	REST	471	UCP1
40	BACE1	112	DCTN6	184	GPX2	256	NDUFA10	328	POLR2H	400	RPS27A	472	ULK2
41	BACE2	113	DDC	185	GPX3	257	NDUFA12	329	POLR21	401	RPS6KA1	473	UQCRB
42	BAD	114	DDIT3	186	GPX5	258	NDUFA13	330	POLR2J	402	RPS6KA2	474	UQCRC1
43	BAX	115	DERL1	187	GPX7	259	NDUFA2	331	POLR2J2	403	RPS6KA3	475	UQCRC2
44	BCL2	116	DKK1	188	GRIA1	260	NDUFA3	332	POLR2K	404	RPS6KA6	476	UQCRFS1
45	BCL2L1 BDNF	117 118	DKK2 DKK4	189 190	GRIA2 GRIN1	261	NDUFA4 NDUFA4L2	333 334	POLR2L PPARG	405	RTN3 RTN4	477	UQCRH UQCRQ
40	BECN1	119	DLG4	190	GRIN1	263	NDUFA5	335	PPARGC1A	400	SDHB	479	VDAC1
48	BID	120	DNAH17	192	GRIN2B	264	NDUFA6	336	PPID	408	SDHC	480	VDAC2
49	BRAF	121	DNAH8	193	GRIN2C	265	NDUFA7	337	PPIF	409	SIN3A	481	WIPI1
50	C1QA	122	DNAI1	194	GRIN2D	266	NDUFA8	338	PPP1CA	410	SLC17A6	482	WIPI2
51	C1QC	123	DNAI2	195	GRM1	267	NDUFA9	339	PPP1CC	411	SLC17A7	483	WNT1
52	C5	124	DNAL1	196	GRM5	268	NDUFAB1	340	PPP2CA	412	SLC18A1	484	WNT10A
53	C6	125	DNAL4	197	GSK3B	269	NDUFB1	341	PPP2CB	413	SLC18A2	485	WNT10B
54	C7	126	DNALI1	198	HAP1	270	NDUFB10	342	PPP2R1A	414	SLC1A1	486	WNT11
55	C8B	127	DRD1	199	HDAC1	271	NDUFB11	343	PPP2R1B	415	SLC1A2	487	WNT16
56	CACNA1C CACNA1S	128 129	DRD2 DUSP1	200	HDAC2 HIP1	272	NDUFB2 NDUFB3	344 345	PPP3CA PPP3CB	416	SLC1A3 SLC25A31	488 489	WNT2 WNT2B
58	CALM15	129	DUSP1 DVL1	201	HRAS	275	NDUFB5 NDUFB4	345	PPP3CB PPP3R1	417	SLC25A51 SLC25A4	489	WNT26
59	CALIVII CALIVII	130	DVL1 DVL2	202	HSPA1A	275	NDUFB5	340	PPP3R1 PPP3R2	418	SLC25A5	490	WNT4
60	CALML3	132	DVL3	204	HSPA5	276	NDUFB6	348	PRKACA	420	SLC25A6	492	WNT5A
61	CALML5	133	EGR1	205	HTRA2	277	NDUFB7	349	PRKACB	421	SLC32A1	493	WNT5B
62	CAMK2A	134	EIF2AK2	206	IDE	278	NDUFB8	350	PRKACG	422	SLC38A1	494	WNT6
63	CAMK2B	135	EIF2AK3	207	IFT57	279	NDUFB9	351	PRKCA	423	SLC38A2	495	WNT7A
64	CAMK2D	136	EIF2S1	208	IKBKB	280	NDUFC1	352	PRKCB	424	SLC38A3	496	WNT7B
	CAMK4	137	ELK1	209	IL1A	281	NDUFS1	353	PRKCG	425	SLC6A1	497	WNT8B
65		138	EP300	210	IL1B	282	NDUFS2	354	PRPH	426	SLC6A3	498	WNT9A
66	CASP1		E D L L L		IL6	283	NDUFS3	355	PSEN1	427	SNCA	499	WNT9B
66 67	CASP3	139	ERN1	211		0.0.1	NIDUCES	055	DOCTOR	4.5.5	0.010.010		VET
66 67 68	CASP3 CASP7	139 140	FADD	212	INSR	284	NDUFS4	356	PSEN2	428	SNCAIP SOD1	500	XBP1
66 67 68 69	CASP3 CASP7 CASP8	139 140 141	FADD FAS	212 213	INSR IRS1	285	NDUFS5	357	PSENEN	429	SOD1	500	XBP1
66 67 68	CASP3 CASP7	139 140	FADD	212	INSR							500	XBP1

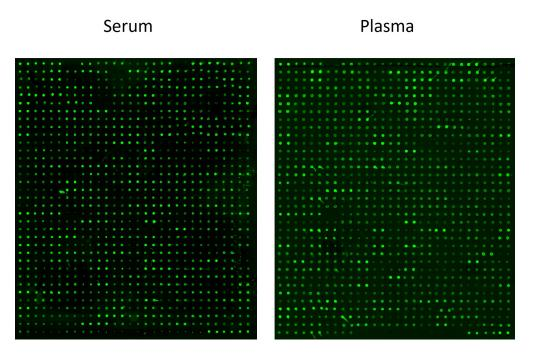
VII. Interpretation of Results

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