RayBio[®] Label-Based (L-Series) Human DNA Damage Screening Array

Patent Pending Technology User Manual (Apr 14, 2023)

For the simultaneous detection of the relative expression of 308 Human proteins in serum, plasma, cell culture supernatants, cell/tissue lysates or other body fluids.

AAH-BLG-DNA-4 (4 Sample Kit) AAH-BLG-DNA-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

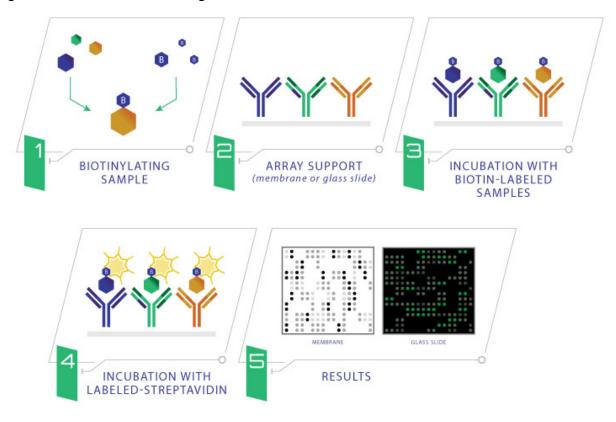
Table of Contents

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

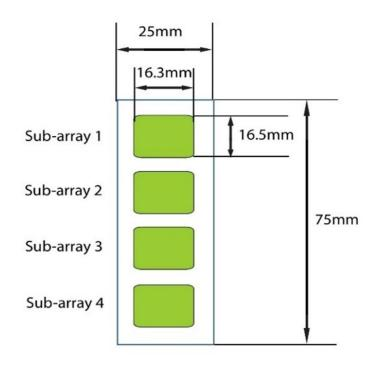
For optimal biotin labeling, it is necessary to determine the protein concentration in the cell/tissue lysate. We recommended using a BCA total protein assay (e.g., Pierce, Catalog # 23227).

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

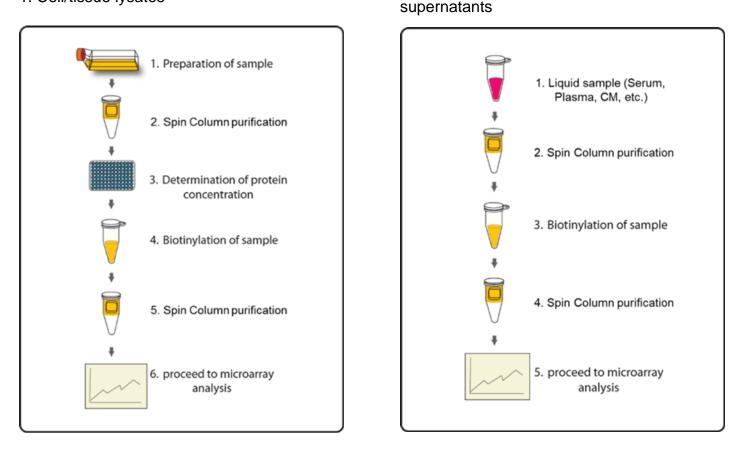
D. Incubations and Washes

- 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.

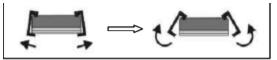
Note: Avoid exposure to light in Steps 19-25 by covering the Glass Slide Assembly with aluminum foil or incubate in a dark room.

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.

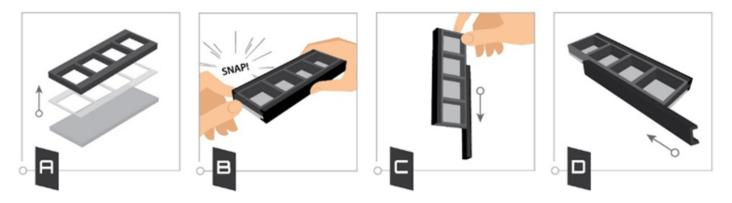
E. Fluorescence Detection

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	282	282	283	283	284	284	285	285	286	286	287	287	288	288	289	289	290	290	291	291	292	292	293	293	294	294	295	295	296	296
21	297	297	298	298	299	299	300	300	301	301	302	302	303	303	304	304	305	305	306	306	307	307	308	308	POS3	POS3	POS2	POS2	POS1	POS1

VI. Antibody Array Target List

Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name
1	AAAS	46	DBI	91	FABP4	136	MUS81	181	POLR2J2	226	PSMD3	271	SPCS2
2	ACADL	47	DCLRE1C	92	FABP5	137	NBN	182	POLR2K	227	PSMD4	272	SRP72
3	ACIN1	48	DCPS	93	FABP6	138	NR1H3	183	POLR2L	228	PSMD6	273	SRRM1
4	ACOX1	49	DDB1	94	FABP7	139	NTHL1	184	POLR3B	229	PSMD7	274	SSBP1
5	ACOX3	50	DDB2	95	FANCA	140	NUP107	185	POLR3E	230	PSMD8	275	SUM01
6	ACSBG1	51	DDX46	96	FANCC	141	NUP188	186	POLR3H	231	PSME1	276	SUMO2
7	ACSL1	52	DDX5	97	FANCD2	142	NUP50	187	POP4	232	PSME2	277	SUMO3
8	ACSL4	53	DHX15	98	FANCG	143	NUP62	188	PPARA	233	PSME3	278	SYCP3
9	ACSL5	54	DIS3L	99	FANCL	144	NUP85	189	PPARD	234	PSMF1	279	TACC3
10	ADIPOQ	55	DNTT	100	FXR1	145	NXF1	190	PPARG	235	RAD23A	280	TAF4
11	ADRM1	56	EEF1A2	101	GTF2A1L	146	NXF3	191	PPIE	236	RAD50	281	TAF4B
12	ANGPTL4	57	EIF1AY	102	GTF2F2	147	NXF5	192	PPIL1	237	RAD51	282	TAF5L
13	APEX1	58	EIF2B1	103	HBS1L	148	NXT2	193	PPP1CA	238	RAD54B	283	TAF7L
14	APOA1	59	EIF2B3	104	HES1	149	OLR1	194	PPP1CC	239	RBM17	284	TBP
15	APOA2	60	EIF2B4	105	HMGB1	150	OXA1L	195	PPP2CA	240	REV1	285	THOC1
16	APOC3	61	EIF2B5	106	HMGCS1	151	PABPC1	196	PPP2CB	241	RFC1	286	THOC3
17	ATM	62	EIF2S1	107	HNRNPA1	152	PABPC3	197	PPP2R1A	242	RFC3	287	THOC7
18	ATR	63	EIF3B	108	HNRNPC	153	PABPC4	198	PPP2R2A	243	RFC4	288	TOB1
19	ATRIP	64	EIF3D	109	HSPA1A	154	PAPOLA	199	PPP2R2B	244	RNASEH1	289	TOPBP1
20	BCAS2	65	EIF3F	110	HSPA1L	155	PARP1	200	PPP2R2C	245	RNASEH2B	290	TXNL4A
21	BRCA1	66	EIF3G	111	HSPA2	156	PARP4	201	PPP2R5C	246	RNGTT	291	U2AF1
22	BRCA2	67	EIF3I	112	HSPA5	157	PCBP1	202	PPP2R5E	247	RNMT	292	UBC
23	BRIP1	68	EIF3J	113	HSPA6	158	PCF11	203	PQBP1	248	RPA2	293	UBE2I
24	BTG2	69	EIF4A1	114	HSPA8	159	PCK1	204	PRKDC	249	RPA3	294	UBE2T
25	BTG3	70	EIF4A3	115	HSPA9	160	PCNA	205	PRMT5	250	RPP14	295	UCP1
26	BTG4	71	EIF4B	116	HSPD1	161	PDPK1	206	PRPF18	251	RXRA	296	UNG
27	BUD31	72	EIF4E	117	IFNG	162	PELO	207	PRPF19	252	RXRB	297	UPF1
28	C1D	73	EIF4EBP1	118	ILK	163	PFKM	208	PSMA1	253	RXRG	298	UPF3B
29	CD36	74	EIF4G1	119	KPNB1	164	PFKP	209	PSMA2	254	SCD	299	WDR33
30	CDC5L	75	EIF4G2	120	LIG1	165	PLTP	210	PSMA8	255	SEC11C	300	WDR61
31	CDK7	76	EIF4G3	121	LIG3	166	PNPT1	211	PSMB10	256	SEC13	301	XPA
32	CETN2	77	EIF5	122	LPL	167	POLB	212	PSMB3	257	SF3B4	302	XPC
33	CHERP	78	ELAC2	123	LSM1	168	POLD4	213	PSMB6	258	SLC27A1	303	XPO1
34	CLNS1A	79	ENO2	124	LSM3	169	POLH	214	PSMB7	259	SLC27A2	304	XRCC1
35	CNOT6	80	ERCC2	125	LSM7	170	POLR1B	215	PSMB8	260	SLC27A4	305	XRCC2
36	CPSF2	81	ERCC3	126	MAGOH	171	POLR2A	216	PSMB9	261	SLC27A5	306	XRCC3
37	CPSF4	82	ERCC5	127	MAGOHB	172	POLR2B	217	PSMC2	262	SLC27A6	307	XRCC5
38	CPSF6	83	ERCC8	128	MBD4	173	POLR2C	218	PSMC4	263	SLU7	308	XRCC6
39	CPT1A	84	EXOSC10	129	MCM2	174	POLR2D	219	PSMC5	264	SMG1		
40	CPT1B	85	EXOSC2	130	MCM7	175	POLR2E	220	PSMC6	265	SMN1	6	
41	CPT2	86	EXOSC4	131	MMP1	176	POLR2F	221	PSMD11	266	SNRPB		
42	CSTF3	87	EXOSC7	132	MSH2	177	POLR2G	222	PSMD12	267	SNRPD1	6	
43	CUL4B	88	FABP1	133	MSH6	178	POLR2H	223	PSMD13	268	SNRPD2		
44	CYFIP2	89	FABP2	134	MSI1	179	POLR2I	224	PSMD14	269	SNRPD3		
45	CYP7A1	90	FABP3	135	MSI2	180	POLR2J	225	PSMD2	270	SNW1		

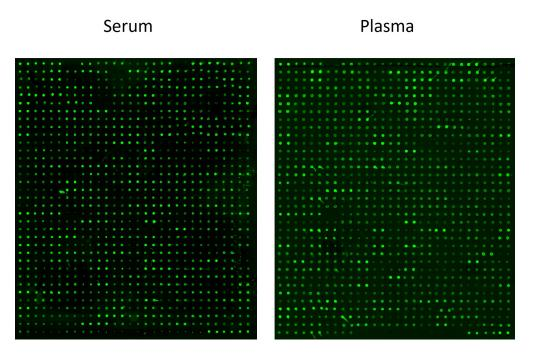
VII. Interpretation of Results

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