

RayBio[®] Label-Based (L-Series)

Mouse Antibody Array L-2808 Membrane Kit

A combination of Mouse L-308, L-2, L-3, L-4, L-5, and L-6 arrays

Patent Pending Technology User Manual (February 1, 2024)

For the simultaneous detection of the relative expression of 2808 mouse proteins in serum, plasma, cell culture supernatants, cell/tissue lysates and other body fluids.

L-Series Mouse Antibody Array L-2808

Cat# AAM-BLM-2808-2 (2 Sample Kit)

Cat# AAM-BLM-2808-4 (4 Sample Kit)

**Please read manual carefully
before starting experiment**



Your Provider of Excellent Protein Array Systems and Services

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

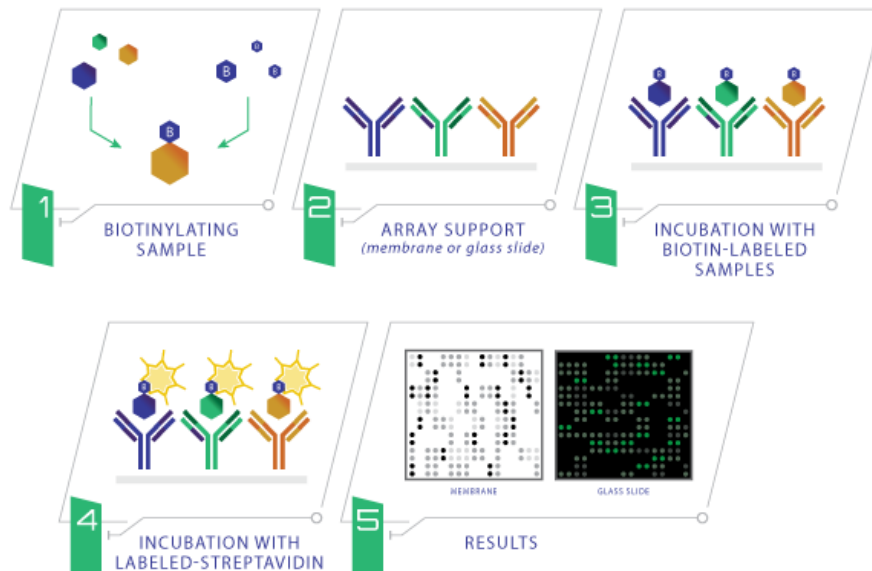
TABLE OF CONTENTS

I.	Introduction and How It Works.....	2
II.	Materials Provided.....	3
	A. Storage Recommendations.....	3
	B. Additional Materials Required.....	3
III.	Overview and General Considerations.....	4
	A. Preparation and Storage of Samples.....	4
	B. Handling the Array Membrane.....	6
	C. Incubation of Antibody Array.....	6
	D. Layout of Array Membrane.....	7
IV.	Protocol.....	8
	A. Sample Purification.....	8
	B. Biotin Labeling of Sample.....	9
	C. Blocking and Incubations.....	10
	D. Detection.....	12
V.	Antibody Array Map	13
VI.	Antibody Array Target Lists.....	15
VII.	Interpretation of Results.....	21
	A. Explanation of Controls Spots.....	21
	B. Typical Results.....	21
	C. Background Subtraction.....	22
	D. Normalization of Array Data.....	22
	E. Threshold of Significant Difference.....	23
	F. Pathway Analysis of the Array Proteins.....	23
VIII.	Troubleshooting Guide.....	24
IX.	Selected References.....	25

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 membrane arrays are then blocked, similar to a Western blot, and the biotin-labeled sample is added onto the membrane array which is pre-printed with capture antibodies and incubated to allow for interaction of target proteins. After incubation with HRP-Conjugated Streptavidin, the signals can be visualized by chemiluminescence.



II. Materials Provided

A. Storage Recommendations

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

ITEM	DESCRIPTION	2 SAMPLE KIT	4 SAMPLE KIT
B	Labeling Reagent	2 vials	4 vials
C	Labeling Buffer	1 bottle (30 ml)	1 bottle (30 ml)
D	Stop Solution	1 vial (50 µl)	1 vial (50 µl)
E	L-series Antibody Array Membranes	2 membranes each of Mouse L-308, L-2, L-3, L-4, L-5, and L-6	4 membranes each of Mouse L-308, L-2, L-3, L-4, L-5, and L-6
F	4X Blocking Buffer	2 bottles (30 ml)	4 bottles (30 ml)
I	500X HRP-Conjugated Streptavidin Concentrate	2 vials (100 µl)	4 vials (100 µl)
K	Detection Buffer C	3 bottles (10 ml)	6 bottles (10 ml)
L	Detection Buffer D	3 bottles (10 ml)	6 bottles (10 ml)
G	20X Wash Buffer 1 Concentrate	3 bottles (30 ml)	6 bottles (30 ml)
H	20X Wash Buffer 2 Concentrate	3 bottles (30 ml)	6 bottles (30 ml)
J-2	Spin Columns (10 ml)	4 columns	8 columns
N/A	Plastic Incubation Trays (w/lid)	12 trays	24 trays
N/A	2X Lysis Buffer	1 bottle (10 ml)	1 bottle (10 ml)
Other Kit Components: Plastic Sheets			

B. Additional Materials Required

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

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

III. Overview and General Considerations

A. Preparation and Storage of Samples

1) Preparation of Cell Culture Supernatants

1. Seed cells at a density of 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.**,†
4. To collect supernatants, centrifuge at 1,000 x g for 10 min and store as ≤ 1 ml aliquots at -80°C until needed.
5. If you want to use cell mass for inter-sample normalization, measure the total wet weight of cultured cells in the pellet and/or culture dish. You may then normalize between arrays by dividing densitometry signals by total cell mass (i.e., express results as the relative amount of protein expressed/mg total cell mass). Or you can normalize between arrays by determining the cell lysate concentration using a total protein assay (BCA Protein Assay Kit, Pierce, Prod #: 23227).

**The density of cells per dish used is dependent on the cell type. More or less cells may be required.*

***Optimal culture time may vary and will depend on the cell line, treatment conditions and other factors.*

†Bovine serum proteins produce detectable signals on the RayBio® L-Series Array in media containing serum concentrations as low as 0.2%. When testing serum-containing media, we strongly recommend testing an uncultured media blank for comparison with sample results.

2) Extracting Protein from Cells

1. Centrifuging Cells:

a. Adherent Cells:

- i. Remove supernatant from cell culture and wash cells gently twice with cold 1X PBS taking care not to disturb cell layer.
- ii. Add enough cold 1X PBS to cover cell layer and use cell scraper to detach cells.

b. Cells in Suspension: Pellet the cells by centrifuging using a microcentrifuge at 1500 rpm for 10 min.

2. Make sure to remove any remaining PBS before adding 1X Cell Lysis Buffer (2X Cell Lysis Buffer should be diluted 2-fold with ddH₂O). Solubilize the cells at 2×10^7 cells/ml in 1X Cell Lysis Buffer.
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 min at 2-8 °C.

Note: If the lysates appear to be cloudy, transfer the lysates to a clean tube, centrifuge again at 13,000 rpm for 20 minutes at 2-8°C. If the lysates are still not clear, store them at -20°C for 20 minutes. Remove from the freezer and immediately centrifuge at 13,000 rpm for 20 minutes at 2-8°C.

4. Transfer lysates to a clean tube. Determine cell lysate concentrations using a total protein assay (BCA Protein Assay Kit, Pierce, Prod# 23227). Aliquot the lysates and store at -80°C.

3) Extracting Protein from Crude Tissue

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

2. Homogenize the tissue according to homogenizer manufacturer instructions.
3. Transfer extracts to microcentrifuge tubes and centrifuge for 20 min at 13,000 rpm (4°C).

Note: If the supernatant appears to be cloudy, transfer the supernatants to a clean tube, centrifuge again at 13,000 rpm for 20 minutes at 2-8°C. If the supernatant is still not clear, store the lysate at -20°C for 20 minutes. Remove from the freezer, immediately centrifuge at 13,000 rpm for 20 minutes at 2-8°C.

4. Transfer supernatant to a clean tube and store at -80°C.

4) Determine the total protein concentration

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

B. Handling the Array Membranes

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

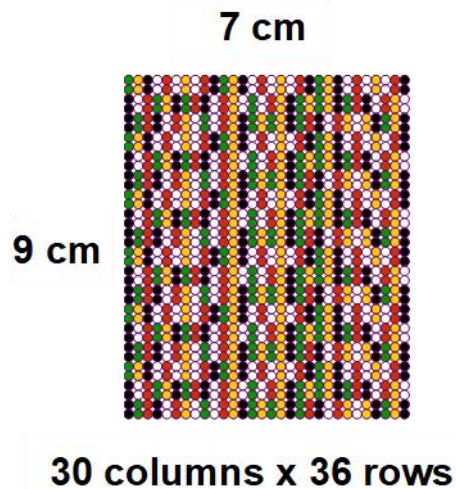
C. Incubations of Antibody Array

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

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

D. Layout of Array Membrane

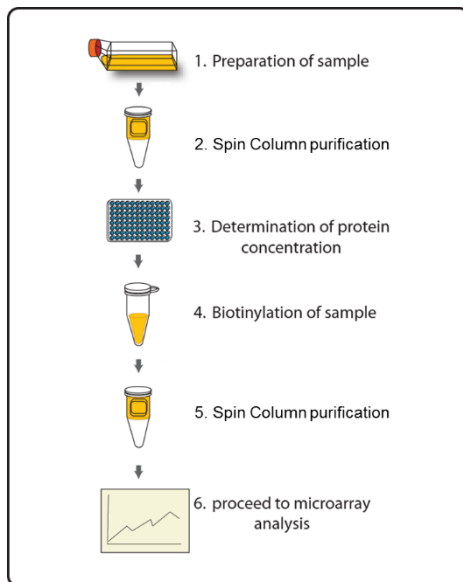
(The layout for AAM-BLM-308 is 30 x 30, while all the rest arrays are 30 x 36)



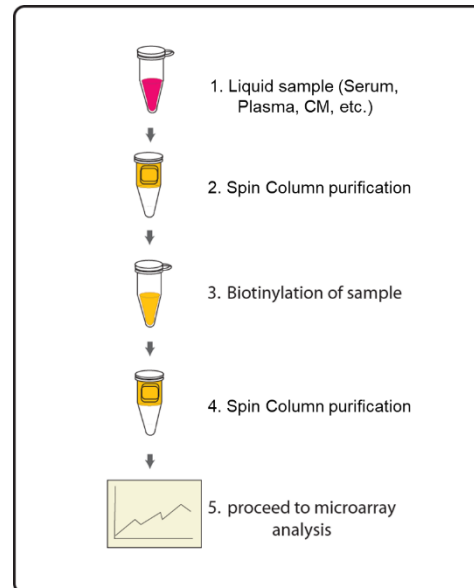
IV. Protocol

Assay Diagram

1. Cell/tissue lysates



2. Serum, plasma, body fluid, or Cell culture supernatants



A. Sample purification

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

1. Twist to remove the bottom closure of the Spin column and loosen the cap (Do not remove).
2. Place the Spin column into a 50 ml conical collection tube, centrifuge at 1,000 x g for 3 minutes to remove the storage buffer. Discard the flow-through.
3. Wash the column three times with 5 ml ultrapure water (ddH₂O) or 1xPBS (pH8.0) each, centrifuge 1,000 x g for 3 minutes to remove the

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

4. Apply sample on top of the resin within the next few minutes. Centrifuge at 1,000 x g for 3 minutes to collect the flow-through that contains sample. The recommended sample dilution as following:

- *Culture Media (CM): 3.5 ml neat supernatant*
- *Serum/Plasma: 50 µl serum/plasma in 3.5 ml labeling buffer*
- *Cell/tissue lysate: 100 µg lysate in 1 ml labeling buffer*

Note: Each labelled sample volume is enough for at least 6 membranes following the protocol below.

Note: The maximal sample volume is 4 ml for each Spin Column. Do not load over 4 ml of sample into a Spin Column.

B. Biotin-Labeling the Sample

Note: Amines (e.g., Tris, glycine) and azides quench the biotinylation reaction. Avoid contaminating samples with these chemicals prior to biotinylation.

5. Immediately before use, prepare Labeling Reagent. Briefly spin down the Labeling Reagent tube (Item B). Add 100 µl Labeling Buffer into the tube, then pipette up and down or vortex slightly to dissolve the lyophilized reagent.
6. Add Labeling Reagent to the sample tube. Incubate the reaction solution at RT with gentle rocking or shaking for 30 min. Mix the reaction solution by gently tapping the tube every 5 minutes.
 - a. For labeling cell culture supernatants: Add 80 µl of Labeling Reagent into the sample tube (for 3.5 ml supernatant).
 - b. For labeling serum or plasma: Add 80 µl of Labeling Reagent Solution into the sample tube (for 50 µl serum/plasma *in 3.5 ml labeling buffer*).

- c. For labeling cell or tissue lysates: Add 8 μ l of Labeling Reagent Solution into the sample tube (for 100 μ g lysate *in 1 ml labeling buffer*).
- d. For all other body fluid: Add 2 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 more or less amount sample is labelled, adjust this volume proportionally.

7. Add 5 μ l Stop Solution (Item D) into each reaction tube. Using a new spin column, repeat Steps 1-4 of section A. Sample Purification to remove the excess non-reacted biotin reagent from each sample.

Note: Biotinylated samples can be stored at -20°C or -80°C until you are ready to proceed with the assay.

C. Blocking and Incubations

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

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

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

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

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

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

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

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

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

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

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

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

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

16. Wash as directed in steps 11 through 13.

D. Detection

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

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

19. The signal can be detected directly from the membrane using a chemiluminescence imaging system or by exposing the array to x-ray film (we recommend using Kodak X-Omat™ AR film) with subsequent development. Expose the membranes for 40 seconds. Then re-expose the film according to the intensity of signals. If the signals are too strong (background too high), reduce the exposure time (e.g., 5–30 seconds). If the signals are too weak, increase the exposure time (e.g., 5–20 min or overnight) or re-incubate membranes overnight with 1X HRP-Conjugated Streptavidin, and repeat detection on the second day.

20. Save membranes at –20 °C to –80 °C for future reference.

V. Antibody Array Maps

A. RayBio® Mouse Antibody Array L-308 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	POS2	POS3	Blank	Blank	Blank	Blank	Blank	1	2	3	4	5	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	6	7	8	9	10	11	12
2	POS1	POS2	POS3	Blank	Blank	Blank	Blank	Blank	1	2	3	4	5	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	6	7	8	9	10	11	12
3	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	13	14	15	16	17	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	18	19	20	21	22	23	24
4	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	13	14	15	16	17	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	18	19	20	21	22	23	24
5	25	26	27	28	29	30	31	32	33	34	35	36	37	Blank	Blank	38	39	40	41	42	43	44	45	46	47	48	49	50	51	52	
6	25	26	27	28	29	30	31	32	33	34	35	36	37	Blank	Blank	38	39	40	41	42	43	44	45	46	47	48	49	50	51	52	
7	53	54	55	56	57	58	59	60	61	62	63	64	65	Blank	Blank	66	67	68	69	70	71	72	73	74	75	76	77	78	79	80	
8	53	54	55	56	57	58	59	60	61	62	63	64	65	Blank	Blank	66	67	68	69	70	71	72	73	74	75	76	77	78	79	80	
9	81	82	83	84	85	86	87	88	89	90	91	92	93	Blank	Blank	94	95	96	97	98	99	100	101	102	103	104	105	106	107	108	
10	81	82	83	84	85	86	87	88	89	90	91	92	93	Blank	Blank	94	95	96	97	98	99	100	101	102	103	104	105	106	107	108	
11	109	110	111	112	113	114	115	116	117	118	119	120	121	Blank	Blank	122	123	124	125	126	127	128	129	130	131	132	133	134	135	136	
12	109	110	111	112	113	114	115	116	117	118	119	120	121	Blank	Blank	122	123	124	125	126	127	128	129	130	131	132	133	134	135	136	
13	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	137	138	139	140	141	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	142	143	144	145	146	147	148
14	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	137	138	139	140	141	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	142	143	144	145	146	147	148
15	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	149	150	151	152	153	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	154	155	156	157	158	159	160
16	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	149	150	151	152	153	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	154	155	156	157	158	159	160
17	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	161	162	163	164	165	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	166	167	168	169	170	171	172
18	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	161	162	163	164	165	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	166	167	168	169	170	171	172
19	173	174	175	176	177	178	179	180	181	182	183	184	185	Blank	Blank	186	187	188	189	190	191	192	193	194	195	196	197	198	199	200	
20	173	174	175	176	177	178	179	180	181	182	183	184	185	Blank	Blank	186	187	188	189	190	191	192	193	194	195	196	197	198	199	200	
21	201	202	203	204	205	206	207	208	209	210	211	212	213	Blank	Blank	214	215	216	217	218	219	220	221	222	223	224	225	226	227	228	
22	201	202	203	204	205	206	207	208	209	210	211	212	213	Blank	Blank	214	215	216	217	218	219	220	221	222	223	224	225	226	227	228	
23	229	230	231	232	233	234	235	236	237	238	239	240	241	Blank	Blank	242	243	244	245	246	247	248	249	250	251	252	253	254	255	256	
24	229	230	231	232	233	234	235	236	237	238	239	240	241	Blank	Blank	242	243	244	245	246	247	248	249	250	251	252	253	254	255	256	
25	257	258	259	260	261	262	263	264	265	266	267	268	269	Blank	Blank	270	271	272	273	274	275	276	277	278	279	280	281	282	283	284	
26	257	258	259	260	261	262	263	264	265	266	267	268	269	Blank	Blank	270	271	272	273	274	275	276	277	278	279	280	281	282	283	284	
27	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	285	286	287	288	289	Blank	Blank	290	291	292	293	294	295	296	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	
28	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	285	286	287	288	289	Blank	Blank	290	291	292	293	294	295	296	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	
29	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	297	298	299	300	301	Blank	Blank	302	303	304	305	306	307	308	Blank	Blank	Blank	Blank	Blank	Blank	POS3	POS2	POS1
30	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	297	298	299	300	301	Blank	Blank	302	303	304	305	306	307	308	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	

B. RayBio® Mouse Antibody Array L-2, L-3, L-4, L-5, and L-6 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	POS2	POS3	Blank	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
2	POS1	POS2	POS3	Blank	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
3	Blank	Blank	Blank	27	28	29	30	31	32	33	34	35	36	37	38	39	40	41	42	43	44	45	46	47	48	49	50	51	52	53
4	Blank	Blank	Blank	27	28	29	30	31	32	33	34	35	36	37	38	39	40	41	42	43	44	45	46	47	48	49	50	51	52	53
5	54	55	56	57	58	59	60	61	62	63	64	65	66	67	68	69	70	71	72	73	74	75	76	77	78	79	80	81	82	83
6	54	55	56	57	58	59	60	61	62	63	64	65	66	67	68	69	70	71	72	73	74	75	76	77	78	79	80	81	82	83
7	84	85	86	87	88	89	90	91	92	93	94	95	96	97	98	99	100	101	102	103	104	105	106	107	108	109	110	111	112	113
8	84	85	86	87	88	89	90	91	92	93	94	95	96	97	98	99	100	101	102	103	104	105	106	107	108	109	110	111	112	113
9	114	115	116	117	118	119	120	121	122	123	124	125	126	127	128	129	130	131	132	133	134	135	136	137	138	139	140	141	142	143
10	114	115	116	117	118	119	120	121	122	123	124	125	126	127	128	129	130	131	132	133	134	135	136	137	138	139	140	141	142	143
11	144	145	146	147	148	149	150	151	152	153	154	155	156	157	158	159	160	161	162	163	164	165	166	167	168	169	170	171	172	173
12	144	145	146	147	148	149	150	151	152	153	154	155	156	157	158	159	160	161	162	163	164	165	166	167	168	169	170	171	172	173
13	174	175	176	177	178	179	180	181	182	183	184	185	186	187	188	189	190	191	192	193	194	195	196	197	198	199	200	201	202	203
14	174	175	176	177	178	179	180	181	182	183	184	185	186	187	188	189	190	191	192	193	194	195	196	197	198	199	200	201	202	203
15	204	205	206	207	208	209	210	211	212	213	214	215	216	217	218	219	220	221	222	223	224	225	226	227	228	229	230	231	232	233
16	204	205	206	207	208	209	210	211	212	213	214	215	216	217	218	219	220	221	222	223	224	225	226	227	228	229	230	231	232	233
17	234	235	236	237	238	239	240	241	242	243	244	245	246	247	248	249	250	251	252	253	254	255	256	257	258	259	260	261	262	263
18	234	235	236	237	238	239	240	241	242	243	244	245	246	247	248	249	250	251	252	253	254	255	256	257	258	259	260	261	262	263
19	264	265	266	267	268	269	270	271	272	273	274	275	276	277	278	279	280	281	282	283	284	285	286	287	288	289	290	291	292	293
20	264	265	266	267	268	269	270	271	272	273	274	275	276	277	278	279	280	281	282	283	284	285	286	287	288	289	290	291	292	293
21	294	295	296	297	298	299	300	301	302	303	304	305	306	307	308	309	310	311	312	313	314	315	316	317	318	319	320	321	322	323
22	294	295	296	297	298	299	300	301	302	303	304	305	306	307	308	309	310	311	312	313	314	315	316	317	318	319	320	321	322	323
23	324	325	326	327	328	329	330	331	332	333	334	335	336	337	338	339	340	341	342	343	344	345	346	347	348	349	350	351	352	353
24	324	325	326	327	328	329	330	331	332	333	334	335	336	337	338	339	340	341	342	343	344	345	346	347	348	349	350	351	352	353
25	354	355	356	357	358	359	360	361	362	363	364	365	366	367	368	369	370	371	372	373	374	375	376	377	378	379	380	381	382	383
26	354	355	356	357	358	359	360	361	362	363	364	365	366	367	368	369	370	371	372	373	374	375	376	377	378	379	380	381	382	383
27	384	385	386	387	388	389	390	391	392	393	394	395	396	397	398	399	400	401	402	403	404	405	406	407	408	409	410	411	412	413
28	384	385	386	387	388	389	390	391	392	393	394	395	396	397	398	399	400	401	402	403	404	405	406	407	408	409	410	411	412	413
29	414	415	416	417	418	419	420	421	422	423	424	425	426	427	428	429	430	431	432	433	434	435	436	437	438	439	440	441	442	443
30	414	415	416	417	418	419	420	421	422	423	424	425	426	427	428	429	430	431	432	433	434	435	436	437	438	439	440	441	442	443
31	444	445	446	447	448	449	450	451	452	453	454	455	456	457	458	459	460	461	462	463	464	465	466	467	468	469	470	471	472	473
32	444	445	446	447	448	449	450	451	452	453	454	455	456	457	458	459	460	461	462	463	464	465	466	467	468	469	470	471	472	473
33	474	475	476	477	478	479	480	481	482	483	484	485	486	487	488	489	490	491	492	493	494	495	496	497	498	499	500	Blank	Blank	Blank
34	474	475	476	477	478	479	480	481	482	483	484	485	486	487	488	489	490	491	492	493	494	495	496	497	498	499	500	Blank	Blank	Blank
35	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank
36	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank	Blank

VI. Antibody Array Target Lists

A. RayBio® Mouse Antibody Array L-308 Target List

Number	Name	Number	Name	Number	Name	Number	Name	Number	Name
1	6CKine	63	DPPIV	125	IGFBP-1	187	IL-28B	249	SCF R
2	Activin A	64	DR3	126	IGFBP-2	188	IL-31	250	SDF-1
3	Activin C	65	Dtk	127	IGFBP-3	189	IL-31 RA	251	SAA1
4	Activin R1B	66	EDAR	128	IGFBP-5	190	Insulin	252	Shh-N
5	Adiponectin	67	EGFR	129	IGFBP-6	191	Integrin beta-2	253	SIGIRR
6	AgRP	68	EG-VEGF	130	IGFBP-L1	192	I-TAC	254	SLPI
7	ALCAM	69	Endocan	131	IGF-1	193	GRO alpha	255	Soggy-1
8	ANGPTL2	70	Endoglin	132	IGF-2	194	Kremen-1	256	SPARC
9	ANGPTL3	71	Endostatin	133	IL-1 alpha	195	Kremen-2	257	Spinesin
10	Amphiregulin	72	Eotaxin-1	134	IL-1 beta	196	Lefty-1	258	TACI
11	Artemin	73	Eotaxin-2	135	IL-1 R4	197	Leptin R	259	TARC
12	Axl	74	Epigen	136	IL-1 R6	198	LEPTIN	260	TCA-3
13	bFGF	75	Epiregulin	137	IL-1 R9	199	LIF	261	IL-27 R alpha
14	B7-1	76	Erythropoietin	138	IL-1 R1	200	LIGHT	262	TECK
15	BAFF R	77	E-Selectin	139	IL-1 R2	201	LIX	263	TFPI
16	BCMA	78	FADD	140	IL-2	202	LRP-6	264	TGF beta 1
17	beta-Catenin	79	FAM3B	141	IL-2 R alpha	203	L-Selectin	265	TGF beta 2
18	BLC	80	Fas	142	IL-2 R beta	204	Lungkine	266	TGF beta 3
19	Betacellulin	81	Fas Ligand	143	IL-3	205	Lymphotactin	267	TGF beta R1
20	Cardiotrophin-1	82	Fc gamma RIIB	144	IL-3 R alpha	206	LTBR	268	TGF beta R2
21	IL-1ra	83	FGF R3	145	IL-3 R beta	207	MAdCAM-1	269	TSP-1
22	CCL28	84	FGF R4	146	IL-4	208	MCP-1	270	CXCL7
23	MIP-1 beta	85	FGF R5 beta	147	IL-4 R	209	MCP-5	271	Tie-2
24	MCP-3	86	FGF-21	148	IL-5	210	M-CSF	272	TIMP-1
25	MCP-2	87	Flt-3 Ligand	149	IL-5 R alpha	211	MDC	273	TIMP-2
26	CCR10	88	FLRG	150	IL-6	212	MFG-E8	274	TIMP-4
27	CCR3	89	Follistatin-like 1	151	IL-6 R	213	MFRP	275	TL1A
28	CCR4	90	Fractalkine	152	IL-7	214	MIG	276	TLR1
29	CCR6	91	Frizzled-1	153	IL-7 R alpha	215	MIP-1 alpha	277	TLR2
30	CCR7	92	Frizzled-6	154	IL-9	216	MIP-1 gamma	278	TLR3
31	CCR9	93	Frizzled-7	155	IL-9 R	217	MIP-2	279	TLR4
32	CD11b	94	Galectin-3	156	IL-10	218	MIP-3 alpha	280	TMEFF1
33	CD14	95	GCSF	157	IL-10 R alpha	219	MIP-3 beta	281	TNF RI
34	CRP	96	GDF-1	158	IL-11	220	MMP-2	282	TNF RII
35	CD27	97	GDF-3	159	IL-12 p40	221	MMP-3	283	TNF alpha
36	CD27 Ligand	98	GDF-5	160	IL-12 p70	222	MMP-9	284	TNF beta
37	CD30 Ligand	99	GDF-8	161	IL-12 R beta 1	223	MMP-12	285	Thrombopoietin
38	CD30	100	GDF-9	162	IL-13	224	MMP-14	286	TRAIL
39	CD40	101	GFR alpha-2	163	IL-13 R alpha 2	225	MMP-24	287	TRAIL R2
40	CD40 Ligand	102	GFR alpha-3	164	IL-15	226	NRG3	288	TRANCE
41	Cerberus 1	103	GFR alpha-4	165	IL-15 R alpha	227	Neurturin	289	TREM-1
42	Chordin-Like 2	104	GITR	166	IL-16	228	NGFR	290	TROY
43	F3	105	GITR Ligand	167	IL-17A	229	NOV	291	TSLP
44	IL-2 R gamma	106	Glut2	168	IL-17 RB	230	Osteoactivin	292	TSLP R
45	IP-10	107	GM-CSF	169	IL-17C	231	Osteopontin	293	TWEAK
46	Cripto-1	108	Granzyme B	170	IL-17D	232	Osteoprotegerin	294	TWEAK R
47	Crossveinless-2	109	Granzyme D	171	IL-17E	233	OX40 Ligand	295	Ubiquitin+1
48	Cryptic	110	Granzyme G	172	IL-17F	234	PDGF-C	296	uPAR
49	CSK	111	Gremlin-1	173	IL-17 RA	235	PDGF R alpha	297	Urokinase
50	CTACK	112	GHR	174	IL-17 RC	236	PDGF R beta	298	VCAM-1
51	CTLA-4	113	HGFR	175	IL-17 RD	237	Pentraxin-3	299	VE-Cadherin
52	CXCL14	114	HGF	176	IL-18 R alpha	238	PF4	300	VEGF-A
53	CXCL16	115	HVEM	177	IL-20	239	PIGF-2	301	VEGFR1
54	CXCR2	116	ICAM-1	178	IL-20 R alpha	240	Progranulin	302	VEGFR2
55	CXCR3	117	ICAM-2	179	IL-21	241	Prolactin	303	VEGFR3
56	CXCR4	118	ICAM-5	180	IL-21 R	242	P-Selectin	304	VEGF-B
57	CXCR6	119	ICK	181	IL-22	243	RAGE	305	VEGF-C
58	EGF	120	IFN-alpha/beta R1	182	IL-22BP	244	RANTES	306	VEGF-D
59	Decorin	121	IFN-alpha/beta R2	183	IL-23	245	RELN beta	307	WIF-1
60	DKK-1	122	IFN-beta	184	IL-23 R	246	Resistin	308	WISP-1
61	Dkk-3	123	IFN-gamma	185	IL-24	247	S100A10		
62	Dkk-4	124	IFN-gamma R1	186	IL-27	248	SCF		

B. RayBio® Mouse Antibody Array L-2 Target List

Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name
1	14-3-3 beta	73	ASGR2	145	CD21	217	D4	289	Fodrin alpha	361	hnRNP A2B1	433	Lubricin
2	14-3-3 zeta	74	ASH2L	146	CD39L4	218	DAN	290	Frizzled 8	362	hnRNP C1+C2	434	LUZP1
3	53BP1	75	ASL	147	CD41	219	DARS2	291	FRY	363	hnRNP G	435	LYZL1
4	AMY1	76	AspAT	148	CD42b	220	DBH	292	FSH-B	364	hnRNP L	436	MAGI2
5	AAT1	77	DNPEP	149	CD48	221	DCXR	293	FTL1	365	hnRNP M	437	MAN1
6	ABAT	78	ASXL1	150	CD5L	222	DDAH1	294	FUCA2	366	hnRNP U	438	MAN1A1
7	ABCF1	79	ATP5A1	151	CD98	223	DDT	295	FUS	367	Hornerin	439	Mannosidase II
8	ABI3BP	80	ATPB	152	CDA	224	DDX3Y	296	G3BP1	368	Hoxb3	440	MAP1A
9	ACAA1	81	B3GNT2	153	CDK2	225	DEFA6	297	G6PD	369	HOXD11	441	MAPRE1
10	ACAA2	82	B4GalT1	154	CED-6	226	Desmocollin 1	298	GALNT2	370	HP1BP3	442	MARCKS
11	ACAA	83	B7-H2	155	CENPF	227	Desmocollin-2	299	GANAB	371	HPD	443	MASP3
12	ACLY	84	BAD	156	CEP57	228	Desmocollin-3	300	GAPDH	372	HRPT1	444	MBD2
13	ACO1	85	BASP1	157	CE51	229	Desmoglein-1	301	GARNL1	373	HRG	445	MBP
14	ACTBL2	86	Bassoon	158	Cezanne	230	Desmoglein-2	302	GART	374	HRP12	446	MCAM
15	ACTC1	87	Bcl2L2	159	CFB	231	Desmoplakin 3	303	Gastrokine 1	375	HSPA1A	447	Mcl-1
16	ACTG1	88	BCOR	160	CFHR1	232	DGK-theta	304	GATM	376	HTRA1	448	MCM
17	ACTG2	89	beta I Spectrin	161	CFI	233	DISC 1	305	GBE1	377	HUWE1	449	MDH1
18	ACTN1	90	beta I Tubulin	162	CFVII	234	DMRN9	306	GCDPF 15	378	IDH1	450	MEP1A
19	ADA	91	beta III Tubulin	163	Chitobiase	235	DOT1L	307	GCLC	379	IFRD1	451	MT-2
20	ADAMDEC1	92	BID	164	Chitotriosidase	236	DPP3	308	GCSH	380	IGF2BP2	452	Metavinculin
21	ADAS	93	BIN2	165	Cholinesterase	237	DRIL1	309	GDA	381	IGFBP7	453	MFAP4
22	ADGRF5	94	Biotinidase	166	CHORDC1	238	DSCAM	310	GDF7	382	IGSF4B	454	MF12
23	ADGRL4	95	BIRC6	167	CHREBP	239	DSPG3	311	GDI1	383	ILK	455	mGLUR5
24	ADH1	96	BMP-1	168	Chromogranin B	240	ECHS1	312	GDI2	384	Inhibin beta	456	Mimcan
25	ADH1C	97	BPGM	169	CKB	241	ECI1	313	Gephyrin	385	Integrin b1	457	MLCK
26	ADH4	98	BPIFB1	170	CLIC1	242	ECM1	314	GFAP	386	Integrin beta 6	458	MMR
27	ADH5	99	BPIFB2	171	CLUP1	243	EEF1G	315	GGCT	387	Integrin a6	459	MN1
28	ADM	100	Brevican	172	CL-P1	244	EEF2	316	GGH	388	IQGAP2	460	Moesin
29	Advillin	101	BRG1	173	CLTA	245	EFEMP2	317	GIP	389	IRE1	461	MP1
30	AEBP1	102	BRSK1	174	CNOT1	246	EFTUD2	318	GLUPR2	390	IRS2	462	MPCA
31	AFG3L2	103	C1QA	175	CO4A2	247	EH03	319	GLUD1	391	ISOC2	463	MPO
32	AGA	104	C1QB	176	Colfin-1	248	Eif4a1	320	Glycoprotein V	392	ITGB4BP	464	MRP 1
33	Aggrecan	105	C1QR	177	COG4	249	ELAVL1	321	GMZA	393	ITIH2	465	MSH6
34	Agrin	106	C1RL	178	COL19A1	250	EMSY	322	GMF beta	394	ITIH3	466	Mtor
35	AGXT	107	C1s	179	COL4A3	251	EN2	323	GNB1	395	ITIH4	467	Multimerin 2
36	Ahsp	108	C4BPA	180	Col6A2	252	Endorepellin	324	GNPTG	396	JAM-A	468	MyBPC3
37	AIFM1	109	C6	181	COL9A3	253	ENO3	325	GOLM4	397	JPT1	469	MYH2
38	AKAP9	110	C8A	182	COLEC10	254	ENSA	326	GOLM1	398	KDM4B	470	MYH6
39	AKR1B1	111	C8G	183	Collagen I a1	255	EPB41	327	GPD1	399	Keratin 36	471	MYH7
40	AKR7A2	112	C9orf40	184	Collagen III	256	EPCR	328	GPLD1	400	KIAA0319L	472	MYHC 2x
41	ALAD	113	CA1	185	Collagen IVa6	257	Ephrin B1	329	GRHRP	401	KIAA1468	473	MYL12B
42	ALDH16A1	114	CA150	186	Collagen IX	258	Eps 15	330	GRP170	402	KLKB1	474	MYOSA
43	ALDH1A1	115	CACNB4	187	Collagen V	259	ERAB	331	GSS	403	KMT2D	475	Myoferlin
44	ALDH9A1	116	Cadherin 22	188	Collagen X	260	ERp29	332	GSTM1	404	KRT31	476	Myosin 18B
45	alpha Actinin 4	117	Cadherin-6	189	Collagen XV	261	Erp57	333	GSTO1	405	KRT33B	477	Myosin9
46	alpha Synuclein	118	CALD1	190	COMP	262	Erp72	334	GSTP1	406	KRT73	478	NABC1
47	alpha Tubulin 4	119	Calpain S1	191	Corneodesmosin	263	ESD	335	Guanylin	407	KRT82	479	NAGLU
48	ALPL	120	Calpastatin	192	Cortactin	264	ESR1	336	GZMM	408	KRT85	480	NAP1L1
49	ALS	121	Calponin-2	193	COTLL1	265	Ezrin	337	H6PD	409	KSR1	481	NAPRT1
50	Alsin	122	Calretinin	194	CPB2	266	FABP5	338	HABP2	410	LAF4	482	NASP
51	Aminoacylase 1	123	Calumenin	195	CPE	267	Factor IX	339	HBB	411	LAIR1	483	NCAM2
52	Aminopeptidase A	124	CAP1	196	CPEB3	268	Factor V	340	HDGF	412	LAMB1	484	Nebulin
53	Androgen Receptor	125	CAPZA1	197	CPM	269	Factor XI	341	Hemoglobin	413	LMNA	485	Nectin-1
54	ANGPTL6	126	CA2	198	CPNE3	270	Factor XII	342	Hemoglobin A1c	414	LMNB2	486	Nectin-3
55	ANGPTL8	127	CA3	199	CRHBP	271	Factor XIII	343	HEXB	415	LAMA2	487	Neogenin
56	Ankrd26	128	Caspase-14	200	Crkl	272	FAH	344	HGFA	416	LAMB2	488	Nesprin2
57	Annexin A1	129	Catalase	201	CRMP2	273	FAM20C	345	HIBADH	417	LAMC1	489	Neurofibromin
58	Annexin A2	130	Cathelicidin	202	CRTAC1	274	FAM3C	346	HINT1	418	LAMP1	490	Neurogranin
59	Annexin A5	131	Cathepsin A	203	CRY2	275	FASN	347	HIP1R	419	LASP1	491	Neuropeptide B
60	Annexin A6	132	Cathepsin G	204	Cyclophilin A	276	FASTKDS5	348	Histone H1.2	420	LCAT	492	Neuropilin-1
61	ANP	133	Cathepsin H	205	Cyclophilin B	277	FBP 38	349	Histone H1.4	421	LCMT2	493	Neurotrimin
62	ANP32A	134	Cathepsin Z	206	Cystatin	278	FDP5	350	Histone H2A	422	LDH-H	494	NF-M
63	Antithrombin III	135	CBS	207	CYTL1	279	FGG	351	Histone H2A.Z	423	LEDGF	495	NIF3L1
64	APLP1	136	CCAR2	208	Cytochrome b5	280	Fibrillin 1	352	Histone H2B K	424	Umbin	496	NME3
65	AQR	137	CCDC126	209	Cytochrome c	281	Fibrinogen-like 2	353	Histone H3.3	425	LIMS1	497	nNOS1
66	ARFGF3	138	CCDC25	210	Cytokeratin 1	282	Fibrinopeptide B	354	Histone H4	426	LMW-PTP	498	Notch-2
67	Arp3	139	CCS	211	Cytokeratin 10	283	Fibulin 3	355	HMGB1	427	LOK	499	NPAS3
68	ARPC2	140	CD109	212	Cytokeratin 13	284	Ficolin 2	356	HMGB2	428	LOX	500	NPM1
69	ARPC3	141	CD133	213	Cytokeratin 14	285	Filamin C	357	HMGB3	429	LOXL1		
70	ARPP19	142	CD148	214	Cytokeratin 15	286	FKBP1A	358	HMGN2	430	LPA		
71	ART3	143	CD155	215	Cytokeratin 20	287	FKBP25	359	HNF-3 alpha	431	LSAMP		
72	ARTS1	144	CD157	216	Cytokeratin 9	288	FKBP51	360	hnRNP A1	432	LTBP4		

C. RayBio® Mouse Antibody Array L-3 Target List

Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name
1	AARE	73	Filaggrin	145	PABP1	217	PREP	289	RPL22	361	SIM2	433	TRAP1
2	ACAT1	74	FITM1	146	PACS1	218	PRG2	290	RPL23A	362	SIRPB1	434	TRAP220
3	ACOT2	75	GARS	147	PNLIP	219	PrP	291	RPL3	363	Six3	435	TRF 2
4	ADAM28	76	GCC2	148	PARVB	220	Profilin 1	292	RPL32	364	SLC4A1	436	TRIM14
5	AHCY	77	GLI-2	149	PCAP	221	Prolargin	293	RPL4	365	SLTRK1	437	Tropomyosin 3
6	AK1	78	GLOD4	150	PCBP1	222	Prosaposin	294	RPL7	366	SLURP1	438	TRP-1
7	AKR1A1	79	GLUL	151	PCBP2	223	PTGDS	295	RPL7A	367	SMAD6	439	TRPS1
8	ALDH2	80	GMPR1	152	PCCA	224	PSMD2	296	RPLP0	368	SMC4	440	Trypsinogen-2
9	DEFA5	81	GOLGA3	153	PCDH12	225	Protein C	297	RPLP2	369	SMPD4	441	TSR2
10	ANKRD9	82	GP2	154	PCDH8	226	Protein Z	298	RPS10	370	SNRPD1	442	TTC3
11	ANXA3	83	gp340	155	PCK2	227	PRR4	299	RPS11	371	SOD1	443	TTF1
12	AP180	84	GTF2F1	156	PCMT1	228	PRRC2A	300	RPS12	372	SOD2	444	TUBA6
13	AP3S2	85	HA1	157	PCNA	229	PRSS23	301	RPS13	373	SOD-3	445	TWF2
14	APLP2	86	HARS	158	PCPE-1	230	PRSS3	302	RPS14	374	Somatoliberin	446	TXNDC15
15	ApoAV	87	HIC1	159	PCSK9	231	PRTN3	303	RPS15A	375	Somatostatin	447	TXNDC4
16	ASPM	88	HIP55	160	PDAP1	232	PSMA1	304	RPS16	376	SORD	448	TXNDC5
17	ASS1	89	H1F0	161	PDE1B	233	PSMA2	305	RPS18	377	SorLA	449	TXNRD2
18	ATOX1	90	HIST1H1B	162	PDIA6	234	PSMA4	306	RPS19	378	SOX4	450	UBA1
19	ATPG	91	HIVEP2	163	PDLIM1	235	PSMA5	307	RPS2	379	SOX5	451	UBE2D3
20	AUTS2	92	hnRNP K	164	PDLIM3	236	PSMA6	308	RPS20	380	SP-D	452	Ube2L3
21	BAI2	93	hnRNP R	165	PDZD2	237	PSMB1	309	RPS23	381	Spectrin	453	UBE2N
22	BarX1	94	HNRNPUL2	166	PEBP1	238	PSMB2	310	RPS25	382	SPEN	454	UCH-L1
23	BBS1	95	HNRPA3	167	PEBP4	239	PSMB3	311	RPS3	383	SPG48	455	UFM 1
24	UBC9	96	HP1 gamma	168	PENK	240	PSMB4	312	RPS3A	384	SPINK5	456	UGGT
25	BLM	97	Importin 7	169	PEPD	241	PSMB5	313	RPS4X	385	SPS2L	457	CMPK1
26	BOLA2	98	Involucrin	170	perilipin-3	242	PSMB6	314	RPS5	386	SPTBN2	458	UNC13D
27	C10orf58	99	ISLR	171	Perilipin-1	243	PSMB7	315	RPS8	387	SPTLC1	459	UNC45A
28	CACNA1H	100	ITPR2	172	Periostin	244	PSMC3	316	RPS9	388	Src	460	UNC5H4
29	Calpain-2	101	ITPR3	173	Periplakin	245	PSMD1	317	RREB1	389	SSC5D	461	UPB1
30	CaMK2	102	KCNAB3	174	Peroxiredoxin-2	246	PSMD5	318	RSF1	390	STAT3	462	UQCRRB
31	CaMK2D	103	LAMA5	175	Peroxiredoxin-3	247	PSMD9	319	RSU1	391	Stathmin 1	463	UQCRH
32	CBL	104	LDB3	176	Peroxiredoxin-1	248	PSME1	320	RUSC1	392	ST11	464	URB
33	CBR1	105	LHPP	177	PFAS	249	PSME2	321	Septin 7	393	STOM	465	URB2
34	CCDC58	106	LIPG	178	PFDN6	250	PTBP1	322	S100A1	394	STXBP2	466	UROC1
35	CCT6A	107	MAP4K4	179	PFKL	251	PTEN	323	S100A11	395	SUCLG1	467	UROD
36	CHCHD3	108	MICALL2	180	PGAM1	252	PTGR1	324	S100A7	396	SUMO3	468	Uroguanylin
37	Cingulin	109	MON2	181	PGAM2	253	PTK7	325	S100A9	397	SVEP1	469	URP2
38	CIT	110	MPST	182	PGK-1	254	PTMA	326	SDC4	398	Symplekin	470	USP14
39	CMG1	111	MRC2	183	PGLS	255	PTPRG	327	SAA4	399	SynCAM	471	USP2
40	CNBP	112	MSH3	184	PG-M	256	PTPRK	328	SBP-1	400	Synemin	472	USP5
41	CNPY2	113	MTA2	185	PGM1	257	PTPRM	329	SC35	401	SYNPO2L	473	Utroglobin
42	Coilin	114	MTHFD1	186	PGRPL	258	PTPRZ	330	SCG	402	Syntaxin 7	474	Utrophin
43	COL8A2	115	MUC5B	187	PHGDH	259	PZP	331	SCN3A	403	TAB182	475	VARS
44	COLEC11	116	MVD	188	Piccolo	260	QARS	332	SCP2	404	Talin1	476	VAP-1
45	COPG2	117	Myosin IIB	189	pIgr	261	QDPR	333	SDNSF	405	TARS	477	VAP-A
46	CORO1B	118	NACA1	190	PIK3C2B	262	QPRT	334	SDPR	406	TAX1BP3	478	VCP
47	CPA3	119	NAGPA	191	PIN	263	Quiescin Q6	335	SECISBP2	407	TBCA	479	VDAC1
48	CPI17 alpha	120	NAV2	192	PIP5K2 alpha	264	Rab1A	336	Secretogranin V	408	TCEB2	480	VILIP3
49	CrkRS	121	NFATC4	193	PISD	265	Rab7a	337	Semaphorin 6B	409	Tcf20	481	Vimentin
50	CRLF3	122	NNT	194	PLA2G1B	266	Ran	338	Semaphorin 7A	410	TCP1 delta	482	VNN1
51	CSRP3	123	NPEP3	195	Plastin 3	267	RanBP1	339	SERBP1	411	TCP1 eta	483	VPS4B
52	CTNNA1	124	NQO2	196	Plastin L	268	RanGAP1	340	Serpin A11	412	TCP1 theta	484	VSIG4
53	CTNND1	125	NSFL1C	197	PLBD2	269	RAP1B	341	Serpin A7	413	TC1P	485	WDR1
54	Cyclophilin F	126	NUCB1	198	PLD4	270	Rbm15	342	Serpin B3D	414	TDIF2	486	WDR44
55	Cystatin C	127	NUP214	199	Plectin	271	RCL	343	Serpin B6	415	Tenascin C	487	WISP2
56	DCAMK1L	128	OAF	200	Plexin B1	272	RECQ4	344	Serpin B8	416	Tenascin XB	488	WNK2
57	Dematin	129	OIT3	201	Plexin B2	273	Reg3A	345	Serpin F2	417	TFF2	489	XPG
58	DIAPH1	130	OPCML	202	PLOD1	274	REV3L	346	Serpin H1	418	TGM3	490	YB1
59	DKC1	131	ORM2	203	PLOD2	275	RHOC	347	Serpin A10	419	Thioredoxin-1	491	SYN1
60	DLST	132	OSBP1	204	Plxdc2	276	RHOG	348	SERPINB1	420	THOP1	492	YY1
61	DMRT1	133	OSCAR	205	PMCA	277	RNASE1	349	SerpinB4	421	TIF1 alpha	493	ZAK
62	Dystrophin	134	OSM R beta	206	PNP	278	RNASET2A	350	SerpinE2	422	TMEM103	494	zbtb11
63	Ebf4	135	Osteoadherin	207	POLD2	279	RLF	351	SerRS	423	TOB2	495	ZBTB4
64	EBP50	136	OTC	208	POLR2A	280	RNASE4	352	SET	424	TOMM70A	496	ZC3H18
65	ECHDC1	137	OTUB1	209	POR	281	Rnose2	353	SEZ6L2	425	TOP2B	497	ZC3H4
66	EHHADH	138	OTUD7A	210	PPOX	282	RP1	354	SF20	426	TPD52L2	498	ZC3H8
67	EIF3D	139	OT-NPI	211	PPP1CC	283	RPL10	355	SHANK1	427	TPM4	499	ZNF295
68	eIF4A2	140	p16 ARC	212	PPP1R9A	284	RPL10A	356	SHC1	428	TPP1	500	Zyxin
69	eIF4GII	141	p23	213	PPP2R1B	285	RPL11	357	SHMT1	429	TPPP3		
70	ENDOD1	142	p39	214	PPP2R4	286	RPL12	358	SHOX	430	TPR		
71	EYA2	143	P4HB	215	PRCP	287	RPL14	359	SHP-1	431	TALDO1		
72	F8	144	p73	216	PRDM13	288	RPL17	360	Siglec-1	432	Transthyretin		

D. RayBio® Mouse Antibody Array L-4 Target List

Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name
1	2B4	73	Cathepsin B	145	Contactin-6	217	FGF-1	289	LAMP2	361	Nope	433	SerpinA6
2	4-1BB	74	Cathepsin C	146	CPA1	218	FGF-10	290	Latexin	362	Notch-1	434	SerpinB10
3	4-1BB Ligand	75	Cathepsin D	147	CPA2	219	FGF-23	291	Layilin	363	Notch-3	435	SerpinB6b
4	A33	76	Cathepsin E	148	CPB1	220	FGF-4	292	LBP	364	NTB-A	436	SerpinD1
5	ACE	77	Cathepsin L	149	CRACC	221	FGF-6	293	LDHA	365	NTRK1	437	SerpinF1
6	ACE-2	78	CCL6	150	CREG	222	FGF-8	294	LDLR	366	NXPH1	438	sFRP-2
7	ACPP	79	CD160	151	CRELD1	223	FGFR2	295	Legumain	367	Olfactomedin-1	439	sFRP-3
8	ACVR2B	80	CD164	152	CRELD2	224	FLRT2	296	LIFR	368	Oncostatin M	440	sFRP-4
9	ADAM15	81	CD180	153	CRISP-4	225	Flt-3	297	LILRC1	369	OX40	441	SGSH
10	ADAM9	82	CD1d1	154	CSNK2A1	226	Follistatin	298	Limitin	370	PAI-1	442	Siglec-2
11	ADAMTS1	83	CD2	155	CSRP1	227	FOLR1	299	LIMPII	371	P-Cadherin	443	Siglec-3
12	AFP	84	CD200	156	CST7	228	Frizzled-2	300	Lipocalin-2	372	PD-1	444	Siglec-E
13	AGER	85	CD200 R1	157	CTSS	229	Frizzled-9	301	LOX-1	373	PDGF-AA	445	Siglec-F
14	AGT	86	CD200R4	158	CXADR	230	Galectin-1	302	LRIG1	374	PD-L1	446	SIGNR1
15	AIMP1	87	CD229	159	CXCL17	231	Galectin-2	303	LRPAP	375	PD-L2	447	SIRPA
16	AKT3	88	CD28	160	Cystatin B	232	Galectin-4	304	LRRC32	376	PECAM-1	448	SIRPB
17	ALK-1	89	CD2F-10	161	DCC	233	Galectin-7	305	LTA4H	377	Persephin	449	SLAM
18	Ameloblastin	90	CD300b	162	DcTRAIL R1	234	Galectin-9	306	LTF	378	PGRP-S	450	Slit2
19	AMICA	91	CD320	163	DcTRAIL R2	235	Gas 1	307	Lumican	379	PILRA	451	Smad5
20	ANG-1	92	CD34	164	DDC	236	Gas 6	308	LYVE-1	380	PILRB	452	SMOC-1
21	ANG-2	93	CD36	165	DDR2	237	G-CSF R	309	MANF	381	PIR-B	453	SorCS2
22	ANG-3	94	CD38	166	Dectin-1	238	GDF-11	310	MAP2K1	382	PLA2G2A	454	SorCS3
23	Angiogenin	95	CD39	167	Dectin-2	239	GFR1	311	Marapsin	383	PLA2G7	455	SPARCL1
24	ANGPTL7	96	CD39L3	168	Dkk-2	240	GLA	312	MARCO	384	Plexin A1	456	SR-AI
25	ANPEP	97	CD4	169	DLL1	241	Glypican 2	313	Matrilin-2	385	Plexin C1	457	STGALNAC2
26	APCS	98	CD44	170	DLL4	242	Glypican 5	314	Matrilin-3	386	PLTP	458	STC-2
27	ApoH	99	CD45	171	DNAM-1	243	gp130	315	Matrilin-4	387	Podocalyxin	459	Syndecan-1
28	ARSA	100	CD47	172	DNER	244	GPVI	316	MBL-1	388	Podoplanin	460	Syndecan-2
29	ARSG	101	CD5	173	DPP7	245	H60	317	MBL-2	389	PPM1A	461	Syndecan-3
30	ART4	102	CD53	174	E-Cadherin	246	HAI-1	318	Mcpt1	390	PRDC	462	TCAM-1
31	ASAH2	103	CD55	175	EFNA3	247	HAI-2	319	Mcpt6	391	PRDX5	463	TCN2
32	ASAH1	104	CD59a	176	EFNA5	248	HDAC8	320	Mcpt7	392	Pref-1	464	Testican 3
33	ASAM	105	CD6	177	EGFL6	249	HIN-1	321	M-CSF R	393	Prostasin	465	TF
34	B2M	106	CD69	178	ELA2	250	Hip	322	MD-1	394	Prss34	466	TFPI-2
35	B7-2	107	CD7	179	EMMPRIN	251	HPGD	323	MDL-1	395	PTK6	467	TFRC
36	B7-H3	108	CD73	180	Endoglycan	252	HPX	324	MEP1B	396	RANK	468	TGF-beta RIII
37	B7-H4	109	CD74	181	Endothelin-1	253	HS6ST3	325	MEPE	397	RBP4	469	TGM2
38	B7-H5	110	CD79B	182	ENPP2	254	ICOS	326	Mer	398	Reg2	470	Thrombomodulin
39	BACE-1	111	CD8 alpha	183	ENPP-7	255	IDS	327	MESDC2	399	Reg3B	471	THSD1
40	BACE-2	112	CD83	184	EpCAM	256	IFNa	328	METAP2	400	Relaxin-1	472	TIGIT
41	BAFF	113	CD90	185	EphA1	257	IFNA2	329	Meteorin	401	Renin 1	473	TIM-1
42	BAMBI	114	CD96	186	EphA2	258	IFNA5	330	METRNL	402	Ret	474	TIM-2
43	BCL2L1	115	CD97	187	EphA3	259	IFNGR2	331	MGL1	403	RGM-A	475	TIM-3
44	BDNF	116	CD99	188	EphA4	260	IGSF8	332	MGL2	404	RGM-B	476	TIM-4
45	beta IG-H3	117	CD99-L2	189	EphA5	261	IL-1 R7	333	MIF	405	RGM-C	477	TLR6
46	beta-NGF	118	CDC37	190	EphA6	262	IL10RB	334	MMP-10	406	ROBO3	478	TNFRH3
47	BLAME	119	CDCP1	191	EphA7	263	IL11RA1	335	MMP-7	407	ROBO4	479	Tpo R
48	BLMH	120	CDNF	192	EphA8	264	IL12RB2	336	MMP-8	408	R-Spondin 1	480	TREM-2
49	BMP-10	121	CEACAM-1	193	EphB3	265	IL-17B	337	MOG	409	R-Spondin 4	481	TremL1
50	BMP-2	122	CES2	194	EphB4	266	IL-18 BpC	338	MSP	410	Ryk	482	TrkB
51	BMP-4	123	CESSA	195	EphB6	267	IL-1F6	339	MSP R	411	S100A13	483	TrkC
52	BMP-5	124	CFH	196	Ephrin-A1	268	IL-1F8	340	NAALADL1	412	S100A3	484	TROP-2
53	BMP-6	125	Chemerin	197	Ephrin-A2	269	IL20RB	341	NCAM-1	413	S100A4	485	Tryptase-5
54	BMP-7	126	CHI3L1	198	Ephrin-A4	270	IL22RA1	342	NCSTN	414	S100A5	486	TSC22D1
55	BMPR-II	127	CHL-1	199	Ephrin-B2	271	IL-28A	343	Nectin-2	415	S100A6	487	TSG
56	BSSP-4	128	Chordin	200	Epimorphin	272	IL-33	344	NELL1	416	SBDS	488	TSG-6
57	BTLA	129	CHST3	201	EPOR	273	IL-34	345	Nephrin	417	SCARA5	489	TSP-2
58	BTN1A1	130	CHST4	202	ErbB2	274	IL36G	346	Nephronectin	418	SCGF	490	TSP-4
59	C1QBp	131	CHST7	203	ErbB3	275	JAM-B	347	Neprilysin	419	Sclerostin	491	TSPAN7
60	C1ra	132	CLEC3B	204	ESAM	276	JAM-C	348	Netrin-1	420	SELPLG	492	Vitronectin
61	C1S1	133	CLEC4F	205	EXTL2	277	JNK1	349	Netrin-4	421	SEMA3C	493	VLDL R
62	C4.4a	134	CLEC9a	206	F2	278	Kirrel2	350	Netrin-G1a	422	SEMA3F	494	Wnt-10b
63	C5a	135	Clusterin	207	FABP4	279	Kirrel3	351	Netrin-G2a	423	SEMA4C	495	Wnt-2b
64	CA10	136	CMG-2	208	FCER2	280	KLB	352	Neurocan	424	SEMA4D	496	Wnt-3a
65	CA12	137	CNDP1	209	FCGR1	281	KLK1	353	Neuroglycan C	425	SEMA4F	497	Wnt-4
66	CA14	138	CNPY3	210	FCGR3	282	KLK7	354	Neuroplastin	426	SEMA4G	498	Wnt-8a
67	CA4	139	CNTF	211	FCGR4	283	Klrb1a	355	Neuropoietin	427	SEMA6A	499	XPNPEP2
68	CA8	140	CNTN1	212	FCRL1	284	KLRC1	356	NgR	428	SEMA6C	500	YM1
69	CA9	141	CNTN2	213	FCRL5	285	KNG1	357	Nidogen-2	429	SerpinA1c		
70	Cadherin-4	142	CNTNAP2	214	FCRN	286	KYNU	358	NKG2D	430	SerpinA1D		
71	Calprotectin	143	COCO	215	FETUB	287	LAG-3	359	Nkp46	431	SerpinA3c		
72	CAMK4	144	Contactin-4	216	Fetuin A	288	LAMA4	360	Noggin	432	SerpinA3N		

E. RayBio® Mouse Antibody Array L-5 Target List

Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name
1	A1BG	73	BTNL2	145	Cited-2	217	EBF-3	289	HSP90	361	Neuropilin-2	433	SEMA3B
2	A2B5	74	Bub3	146	c-Jun	218	EBI3	290	HSPA8	362	NFIL3	434	SEMA3E
3	A2M	75	C1qTNF10	147	CLEC12A	219	EDG-1	291	HSPB8	363	NGK2C	435	SEMA4A
4	ACTA1	76	C3	148	CLEC1B	220	EEA1	292	ID1	364	NKX3.1	436	SerpinA5
5	ACTA2	77	C3AR1	149	CLEC2d	221	EED	293	IFNA4	365	NLRP3	437	Serpin1
6	ACTN2	78	C5AR1	150	CLEC2i	222	EGLN1	294	IFNK	366	Nodal	438	SEZ6L
7	ADAM10	79	Cadherin-17	151	CLEC4A	223	EHMT1	295	IGF1R	367	NPTXR	439	SHIP
8	ADAM19	80	Calreticulin	152	CLEC4B	224	EHMT2	296	IGFBP-rP10	368	NR1F3	440	SHIP2
9	ADAM33	81	CAMK2A	153	c-Myc	225	eIF4E	297	IGSF4C	369	NR4A2	441	SHP-2
10	ADAMTS15	82	CARD11	154	COL13A1	226	eIF5A	298	IGSF9	370	NRAGE	442	Siglec-G
11	ADCYAP1	83	Carm1	155	Collagen II	227	ELF3	299	ihh	371	NXPH-3	443	SIRP delta
12	ADGRE1	84	CART	156	Collectrin	228	EMAP-II	300	IkB-epsilon	372	Oct-3/4	444	SLC39A4
13	Adipsin	85	Caspase-3	157	Complexin-1	229	EMX2	301	IKK beta	373	OLFM-L3	445	Smad9
14	ADNP	86	Caspase-4	158	Contactin-3	230	Endomucin	302	IL-17RE	374	Olig1	446	SOX2
15	ADRB2	87	Caspase-7	159	COX-1	231	Enolase 2	303	IL-36Ra	375	Opticin	447	Spi-1
16	Adropin	88	Cathepsin 3	160	COX4	232	ENTPD2	304	IL-38	376	p38 alpha	448	Spi-B
17	Afamin	89	Cathepsin 6	161	CPXM1	233	ENTPD6	305	IL-3R beta	377	p38 gamma	449	SSEA-3
18	AGP	90	Cathepsin 7	162	CRADD	234	EOMES	306	IMPAD1	378	Park7	450	STAT5b
19	AHR	91	Cathepsin F	163	CREB	235	EPAS1	307	Intelectin-2	379	PARP	451	STAT6
20	AIRE	92	Caveolin-1	164	CRISP-1	236	ERK1	308	IRF1	380	Patched 1	452	STK3
21	AJAP1	93	Caveolin-2	165	Crk	237	ESM-1	309	IRF3	381	Patched 2	453	SUMF1
22	Akt1	94	CBFB	166	CRY1	238	Ets-1	310	IRF5	382	PCK1	454	SUMF2
23	Akt2	95	CBP	167	CRYAA	239	ETV2	311	IRF6	383	PD-ECGF	455	Syntaxin 12
24	AMH	96	CCDC134	168	CRYAB	240	EXTL1	312	ISLR-2	384	PDGFRL	456	Syntaxin 1A
25	AMIGO2	97	CCK-AR	169	CSE1L	241	EZH2	313	ITGA10	385	PEAR1	457	TAGLN
26	AMIGO3	98	CCN2	170	CSNK1A1	242	FABP1	314	ITGA11	386	PGCP	458	TARDBP
27	Amnionless	99	CCR1	171	CSNK1D	243	FABP2	315	ITGA2	387	PI16	459	TDRD1
28	AMPH	100	CCR2	172	CSNK1E	244	FABP3	316	ITGA3	388	Pin1	460	Tex11
29	ANGPTL4	101	CCR5	173	CSNK1G1	245	FAM3D	317	ITGA5	389	PKC gamma	461	TEX19.1
30	Annexin A11	102	CCR8	174	CSNK2B	246	FANCD2	318	ITGA7	390	PKN1	462	TGN38
31	Annexin A13	103	CCR2L2	175	CTHRC1	247	FAP	319	ITGA9	391	Plasminogen	463	Thioredoxin-2
32	Annexin A4	104	CD101	176	CTRP3	248	FBP1	320	ITGAE	392	PLET-1	464	THRSP
33	Annexin A7	105	CD116	177	Cubilin	249	FCMR	321	ITGB3	393	Plexin A4	465	TICAM2
34	APBA2	106	CD11c	178	CX3CR1	250	FCRLA	322	ITGB4	394	Plfr	466	TLE3
35	APBB1IP	107	CD151	179	CXCL3	251	FCRLB	323	ITGB7	395	PLUNC	467	TLR11
36	APE	108	CD177	180	CXCR1	252	FGF-15	324	JAB1	396	PLZF	468	TLR12
37	ApoA4	109	CD200R1L	181	CXCR5	253	FGF-16	325	Jagged 2	397	Podocan	469	TLR9
38	ARC	110	CD205	182	CXCR7	254	FGL1	326	Kallikrein 5	398	PON3	470	TM4SF1
39	ARG1	111	CD209d	183	Cyclin A2	255	Fgr	327	Keap1	399	PPP1CA	471	Tmp21
40	ARHGAP1	112	CD209e	184	Cyclin B2	256	FMOD	328	Klre-1	400	PPP1R2	472	TORC1
41	ASCL1	113	CD24	185	Cyclin D1	257	FOLR4	329	Laminin-1	401	PPP2CA	473	TORC2
42	ATF1	114	CD248	186	Cystatin D	258	FosB	330	LAR	402	PRAT4B	474	TORC3
43	ATG3	115	CD2AP	187	DACH2	259	FoxC2	331	LC3B	403	PRDM1	475	TREML2
44	ATG4A	116	CD300a	188	DACT3	260	FoxD3	332	Lefty-2	404	PRDM14	476	TRHDE
45	ATG5	117	CD300f	189	DAPP1	261	FoxN1	333	Lgr5	405	PRKACB	477	TRIM5
46	ATG7	118	CD300g	190	DARC	262	FoxO3	334	Lipin 3	406	PRKCI	478	Troponin T
47	ATN1	119	CD302	191	DCBLD2	263	Frizzled-3	335	Lipocalin-13	407	Properdin	479	TSP50
48	Axin-1	120	CD3D	192	DCTN1	264	GAD1	336	LPHN3	408	PROS1	480	TSPAN8
49	B4GAT1	121	CD3E	193	DCTN2	265	GATA-1	337	LY6G	409	PSD-95	481	Tsukushi
50	BACH1	122	CD59b	194	DCX	266	GATA-2	338	MC3R	410	PVRIG	482	UBE2K
51	BAG1	123	CD63	195	DDX1	267	GATA-3	339	M-Cadherin	411	Pygopus-1	483	ULBP-1
52	BAG4	124	CD68	196	DDX5	268	GBL	340	MCSP	412	Rad17	484	Urotensin-II R
53	BAG6	125	CD79A	197	Desmin	269	GDF-15	341	MEA-1	413	RAE1	485	UTF1
54	BAK	126	CD81	198	DGCR2	270	GFRA1	342	MEK2	414	Rae-1 epsilon	486	VDBP
55	Bax	127	CD9	199	DGKE	271	GH	343	Metadherin	415	RAMP2	487	VG5Q
56	BCAM	128	CD94	200	DGKG	272	Ghrelin	344	MFAP2	416	Rap1A	488	VGf
57	Bcl-10	129	CDC14	201	DGKI	273	GLG1	345	MGMT	417	RARRES1	489	VHR
58	Bcl-2	130	CDC25A	202	DHFR	274	GLRX1	346	MINA	418	Ras-GAP	490	VSIG2
59	Bcl-6	131	CDC25C	203	Dhh	275	Glypican 1	347	MIXL1	419	RBBP4	491	VWC2
60	BEX3	132	CDC73	204	DIXDC1	276	GPR64	348	MKI67	420	Reg3D	492	WARP
61	Blk	133	CDK4	205	DMP-1	277	GPR77	349	MKK6	421	Renin R	493	Wnt-5a
62	BLVRB	134	CDO	206	DNMT3A	278	GRIN1	350	MR1	422	RIPK4	494	Wnt-5b
63	BMP-8b	135	CEACAM-2	207	DOC2A	279	GRIN2A	351	MS4A4B	423	ROCK2	495	WWOX
64	BMPR-IB	136	CELF1	208	DOK7	280	GRK2	352	MUCDHL	424	R-Spondin 2	496	XPB
65	BNIP3L	137	CELSR2	209	DPPA3	281	HAND2	353	MUP-1	425	RTVP-1	497	XPD
66	BOC	138	CES3	210	DPPA4	282	Haptoglobin	354	Myoglobin	426	RUNX3	498	ZAG
67	BOK	139	CHD1	211	DPPA5	283	H8-EGF	355	NAC1	427	S100A16	499	ZNF281
68	B-Raf	140	Chk1	212	DYRK2	284	Hepcidin	356	NAIP	428	SIP5	500	ZNRF3
69	BRCA1	141	Chk2	213	DYRK3	285	HMOX2	357	N-Cadherin	429	S5a		
70	BST-2	142	CHMP2B	214	E2A	286	HRP-1	358	NEDD4	430	SCD-1		
71	BTF3	143	ciAP-2	215	EBF-1	287	HSF1	359	NELL2	431	SCGB3A2		
72	BTN2A2	144	CIB1	216	EBF-2	288	HSF4	360	Neurofascin	432	SEC13		

F. RayBio® Mouse Antibody Array L-6 Target List

Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name
1	11 beta-HSD1	73	FKBP12.6	145	HSP27	217	MBD4	289	PAK6	361	Rad23	433	SYT1
2	14-3-3 epsilon	74	FKBP13	146	HSP32	218	MCHR1	290	Pannexin-1	362	Ra1A	434	TAB1
3	14-3-3 eta	75	FKBP38	147	HSP60	219	MCM2	291	PARL	363	Rap2A	435	TAF5
4	14-3-3 gamma	76	FLIP	148	HSP90AA1	220	MCPIP1	292	Pax3	364	RARB	436	TANK
5	14-3-3 sigma	77	FNDC5	149	HSP90B1	221	MDM2	293	Pax5	365	RASSF2	437	TBX5
6	14-3-3 theta	78	FoxA2	150	HSPB7	222	MEF2C	294	Pax7	366	RBFOX3	438	TC-PTP
7	5T4	79	FoxF1	151	HSPH1	223	MEKK2	295	Paxillin	367	Reelin	439	Tec
8	AARS	80	FoxJ3	152	HTRA2	224	MIB1	296	PBR	368	Rel1	440	Tenascin R
9	ABCG2	81	FoxP2	153	IFITM1	225	Midkine	297	PDCD5	369	RelA	441	Teneurin-2
10	ACHE	82	FRG1	154	IFITM3	226	MIOS	298	PDCD6	370	Relaxin R1	442	Teneurin-4
11	ACO2	83	Frk	155	IFNA14	227	Mitofusin 2	299	PDGF-BB	371	Rheb	443	TER-119
12	ADAMTS10	84	FRMPD4	156	IGF2R	228	MKK3	300	PDGFD	372	RND3	444	TH
13	ADAMTS4	85	FRS2	157	IGFBP4	229	MLK4 alpha	301	PDHX	373	RNF168	445	THAP11
14	AGO2	86	FSTL4	158	IGLL1	230	MNX1	302	PDX-1	374	RNF2	446	Themis
15	ALDH3A2	87	GABAB R1	159	IGSF5	231	MPP5	303	PDZK1	375	ROCK1	447	Tie1
16	alpha-Interneixin	88	GABAB R2	160	Ikb-alpha	232	MSH2	304	PEA-15	376	RP56	448	TIM-5
17	ANGEL1	89	GABRA1	161	IKK epsilon	233	MSK2	305	PFKM	377	RPTOR	449	TIM-6
18	APOF	90	GABRA4	162	IKK gamma	234	MSMB	306	PGPEP-1	378	RSK1	450	TLR7
19	ART1	91	GABRA5	163	IL13Ra1	235	MSX1	307	PHOX2B	379	R-Spondin 3	451	TMSB10
20	ASGR1	92	GAD2	164	IL-19	236	Musashi-1	308	PIM2	380	SALL4	452	TNFRSF21
21	ATP1A3	93	GALNT10	165	ILKAP	237	MYB	309	PIWIL1	381	SAM68	453	TNNC1
22	B4GALT6	94	GATE-16	166	IMPA1	238	MyD88	310	PKLR	382	SAMSN1	454	TRAF-2
23	BCAP	95	GCK	167	IMPDH1	239	MYF-5	311	PKM1	383	SAP	455	TRAF-3
24	BMP-9	96	GDF-6	168	IMPDH2	240	MYL2	312	PKN2	384	SATB1	456	TREM-3
25	BMPR1A	97	GDPD2	169	ING1	241	Myocardin	313	PLA2G12B	385	SCAMP3	457	TREML4
26	Brachyury	98	GGT1	170	INOS	242	Myogenin	314	PLC-beta 1	386	SDF2	458	TRIM32
27	BVES	99	GIT1	171	Intelectin-1	243	NADK	315	PLC-beta 4	387	Secretagogin	459	TRIM63
28	Cadherin-13	100	GLI-1	172	IRF2	244	Nanog	316	PLC-delta3	388	SECTM1A	460	TRP14
29	Calcineurin A	101	GLI-3	173	ITGA1	245	NBS1	317	PLC-gamma 1	389	SEMA5A	461	TRRAP
30	CAMK1	102	GLIS1	174	ITGA8	246	NCOA3	318	PLC-gamma 2	390	SEN8	462	TRXR1
31	CCBL1	103	GLRX3	175	ITGAV	247	NCOR1	319	Plexin A2	391	SerpinB3c	463	TSC2
32	CD19	104	Glucagon	176	ITGB5	248	NDFP1	320	Plexin A3	392	SEZ6	464	UBASH3B
33	CD300E	105	Glyoxalase I	177	Jagged 1	249	NEDD8	321	Plexin B3	393	SFRP1	465	UBR5
34	CD84	106	Glyoxalase II	178	Jak1	250	NEGR1	322	PNPLA2	394	SFXN5	466	UCH-L3
35	CEBPB	107	Glypican 3	179	Jak2	251	NEK3	323	PNUTS	395	SH2B1	467	UCP1
36	CEL	108	Glypican 6	180	JARID2	252	Nesfatin-1	324	PON1	396	SH2D2A	468	UCP2
37	CHCHD4	109	GPX1	181	JIP1	253	Nestin	325	POU5F1	397	SH3GL2	469	UNC13A
38	CHODL	110	GPX2	182	JMJD6	254	NETO2	326	PPA1	398	Shisa-4	470	Uromodulin
39	CIDEB	111	GPX3	183	JNK2	255	NeuroD2	327	PPM1B	399	SIN3A	471	USP7
40	CLEC14A	112	GRAP2	184	Junb	256	Neurolysin	328	PPM1L	400	SLAIN1	472	VAMP-1
41	CLEC4N	113	GRB2	185	KDM4A	257	NF2	329	PPP3R1	401	SLC1A3	473	VAMP-2
42	CLOCK	114	GRB7	186	Keil	258	NFKB1	330	PRDM16	402	SLC27A3	474	VANGL2
43	CLTC	115	GRK1	187	Kir2.1	259	NF-L	331	PRDX4	403	SLC6A3	475	Vav-1
44	CNDP2	116	GRK5	188	KIR3DL1	260	Nicalin	332	PRDX6	404	Slit3	476	VDR
45	CoREST	117	GRK7	189	KIRREL1	261	Ninjurin-2	333	PRKAA1	405	Smad2	477	Visfatin
46	COX-2	118	GRM1	190	KITLG	262	NIPP1	334	PRKAA2	406	Smad7	478	VPREB1
47	CRABP2	119	GRP78	191	KLF4	263	NKX1-2	335	PRKACA	407	SMURF2	479	VSIG1
48	c-Rel	120	GSAP	192	KLF5	264	NKX2.2	336	PRKAR1A	408	SNAP25	480	VSIG3
49	DARPP-32	121	GSK3A	193	KLK6	265	NLRP10	337	PRKAR2A	409	Snapin	481	WASF1
50	DRAK2	122	GSK3B	194	Klotho	266	NM23-H1	338	PRKCA	410	SOCS-3	482	WASF3
51	DZIP1	123	GSN	195	KLRG1	267	Nogo-B	339	PRKCB	411	SOCS-4	483	WDR5
52	E2f6	124	GYPA	196	KMO	268	Norrin	340	PRKCQ	412	SOSTDC1	484	WNK1
53	ECD	125	HAO-1	197	KRT17	269	NPC1	341	PRL-3	413	SOX1	485	Wnt-1
54	EFNB3	126	Hck	198	L1CAM	270	NPR1	342	PRLR	414	SPOCK2	486	Wnt-10a
55	EGFL7	127	HCLS1	199	LAMP3	271	NRARP	343	PRMT1	415	SPRY2	487	Wnt-11
56	EphB2	128	HDAC2	200	LHX5	272	NrCAM	344	Prohibitin 2	416	SQSTM1	488	WNT16
57	ERAS	129	HDAC4	201	LILRB4	273	Nrf1	345	Proliferin	417	SR140	489	Wnt7a
58	ERBB4	130	HFE	202	LIPA	274	Nrf2	346	Protogenin	418	SSEA-1	490	Wnt-8b
59	ERK4	131	HHEX	203	Lipin 2	275	NSA2	347	PRPH	419	SSEA-4	491	Wnt-9b
60	ERK5	132	HIF1A	204	LMO4	276	NUAK1	348	Prss21	420	ST3GAL2	492	XIAP
61	ESGP	133	HIF-1AN	205	LPL	277	Nucleostemin	349	PTGDR2	421	STAT4	493	YAP1
62	ESRRB	134	Histamine H3 R	206	LPP	278	NUP85	350	PTGFRN	422	STAT5a	494	Yes
63	EVI2B	135	Histone H2AY	207	LRP1	279	N-WASP	351	PTN	423	STIM1	495	ZAP70
64	FABP6	136	HMGAI8	208	LRP2	280	Olig2	352	PTP1B	424	STMN2	496	ZBTB7A
65	FAK	137	HMGGA2	209	LRRC4	281	Olig3	353	PTP-MEG2	425	STX6	497	ZFP42
66	Fascin	138	HNFA4G	210	LRRN1	282	Omgp	354	PUM1	426	Substance P	498	ZFP57
67	FCAMR	139	HOMER1	211	LRRTM2	283	Orexin A	355	PVALB	427	SULF2	499	ZIC1
68	FGF-18	140	Hoxb8	212	LRRTM3	284	Osteocrin	356	PXN	428	SUZ12	500	ZIC3
69	FGF-7	141	HSSST4	213	Ly6	285	OTOR	357	Rab11A	429	SWAP70		
70	FGF-9	142	HSP2	214	MAEA	286	p55PIK	358	Rab27a	430	Syntaxin 1B		
71	FGFR1	143	HSP10	215	MAFK	287	PAK3	359	RAB9A	431	Syntaxin 8		
72	FIZZ1	144	HSP20	216	MAPK1	288	PAK4	360	RACK1	432	Syntaxin-BP1		

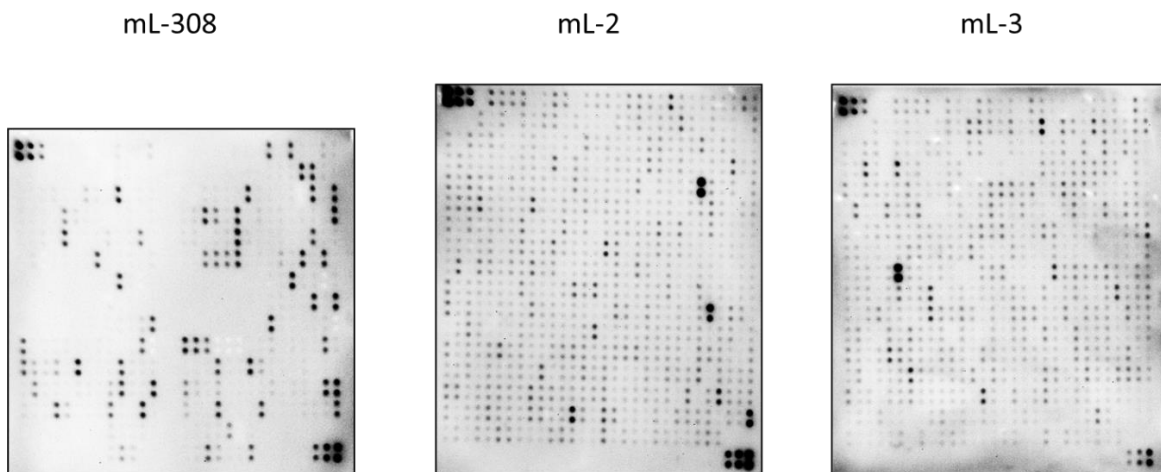
VII. Interpretation of Results:

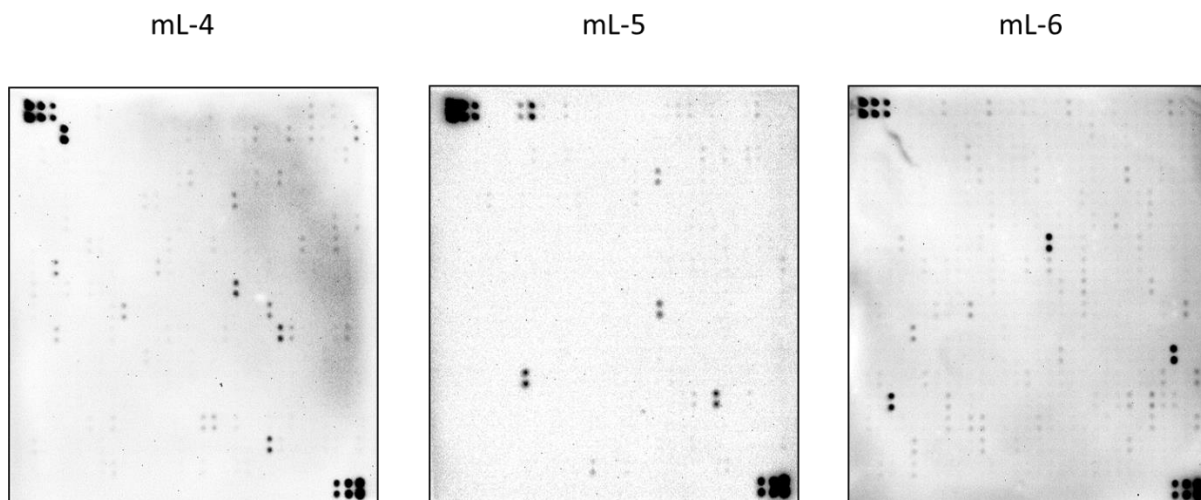
A. Explanation of Controls Spots

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

B. Typical Results

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





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

C. Background Subtraction

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

D. Normalization of Array Data

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

You can calculate the normalized values as follows:

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

Where:

P1 = mean signal intensity of POS spots on reference array

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

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

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

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

E. Threshold of Significant Difference

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

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

F. Pathway Analysis of the Array Proteins

Mouse antibody array L-2808 detects 2808 unique mouse proteins, including most analyzed cytokines, chemokines, adipokines, extracellular matrix proteins, growth factors, angiogenic factors, proteases, enzymes, soluble and transmembrane receptors and transport proteins, adhesion molecules and other proteins. All the array proteins are provided with their Uniprot number and GeneID, which are essential for further data mining. Raybiotech offers affordable biostatistics and bioinformatics service, including data clean-up,

differential expression analysis, cluster analysis, biomarker selection, pathway analysis and experimental design. See more details on the website: <https://www.raybiotech.com/biostatistics-and-bioinformatics-services>

VIII. Troubleshooting Guide

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

IX. Selected References

Christina Scheel et al., *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 hepatocellular 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 al., *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 and LV 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 endometriosis. 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.

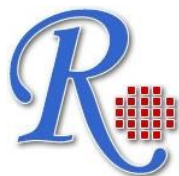
This product is intended for research only and is not to be used for clinical diagnosis. Our products may not be resold, modified for resale, or used to manufacture commercial products without written approval by RayBiotech, Inc.

Under no circumstances shall RayBiotech be liable for any damages arising out of the use of the materials.

Products are guaranteed for six months from the date of shipment when handled and stored properly. In the event of any defect in quality or merchantability, RayBiotech's liability to buyer for any claim relating to products shall be limited to replacement or refund of the purchase price.

RayBio® is a registered trademark of RayBiotech, Inc.

This product is for research use only.



©2024 RayBiotech, Inc.