# RayBio<sup>®</sup> C-Series Human Protein Tyrosine Phosphorylation Antibody Array C3

For the semi-quantitative detection of 500 Tyrosine-phosphorylated human proteins in cell and tissue lysates

**Patent Pending Technology** 

User Manual (Revised Aug. 15<sup>th</sup>, 2022)

Cat# AAH-PTYR-3-2 (2 Sample Kit) Cat# AAH-PTYR-3-4 (4 Sample Kit)

Please read manual carefully before starting experiment



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



C-Series Antibody Arrays

# TABLE OF CONTENTS

I.	Introduction	2
II.	How It Works	3
III.	Components and Storage	4
IV.	Additional Materials Required	4
V.	Sample Tips and General Considerations	5
	A. Sample Collection, Preparation, and Storage	5
	B. Sample Types and Recommended Dilutions/Amounts	5
	C. Handling Membranes	5
	D. Incubations and Washes	6
VI.	Chemiluminescence Detection Tips	6
VII.	Component Preparation	7
VIII.	Protocol	8
	A. Blocking	8
	B. Sample Incubation	8
	C. First Wash	8
	D. Biotinylated Antibody Cocktail Incubation	8
	E. Second Wash	9
	F. HRP-Streptavidin Incubation	9
	G. Third Wash	9
	H. Chemiluminescent Detection	9
	I. Storage	10
IX.	Typical Results	10
Х.	Interpreting the Results	11
	A. Control Spots	11
	B. Data Extraction	11
	C. Data Analysis	11
XI.	Array Map	13
XII.	Troubleshooting Guide	14

# I. INTRODUCTION

Protein phosphorylation plays an unusually prominent role in cell signaling, development and growth. The RayBio<sup>®</sup> Human Protein Tyrosine Phosphorylation Antibody Array C3 is a very rapid, convenient and sensitive assay to simultaneous detect multiple protein phosphorylations and can be used to monitor activation or function of important biological pathways.

RayBiotech is committed to developing a series of phosphorylation antibody arrays. RayBio<sup>®</sup> Human Protein Tyrosine Phosphorylation Antibody Array C3 is specifically designed for simultaneously identifying the relative levels of Tyrosine phosphorylation of 500 different human proteins in cell lysates. By monitoring the changes in protein tyrosine phosphorylation in your experimental model system, you can verify pathway activation in your cell lines without spending excess time and effort in performing an analysis of immunoprecipitation and/or Western Blot.

By using RayBio<sup>®</sup> Human Protein Tyrosine Phosphorylation Antibody Array C3, treated or untreated cell lysate is added into antibody array membranes. The antibody array membranes are washed and biotinylated anti-phosphotyrosine antibody is used to detect phosphorylated tyrosines on target protein. After incubation with HRP-streptavidin, the signals are visualized by chemiluminescence.

RayBio<sup>®</sup> C-Series Antibody Arrays have several advantages over detection of cytokines using single-target ELISA kits:

- 1. <u>More Data, Same or Less Sample</u>: Antibody arrays provide high-content screening using about the same sample volume as traditional ELISA.
- 2. <u>Global View of Cytokine Expression</u>: Antibody array screening improves the chances for discovering key factors, disease mechanisms, or biomarkers related to cytokine signaling.
- 3. <u>Similar (sometimes better) Sensitivity</u>: As little as 4 pg/ml of MCP-1 can be detected using the C-Series array format. In contrast, our similar MCP-1 ELISA assay has a sensitivity of 40 pg/ml of MCP-1.
- 4. <u>Increased Range of Detection</u>: ELISA assays typically detect a concentration range of 100- to 1000-fold, however, RayBiotech arrays can detect IL-2 at concentrations of 25 to 250,000 pg/ml, a range of 10,000-fold.
- 5. <u>Better Precision</u>: As determined by densitometry, the inter-array Coefficient of Variation (CV) of spot signal intensities is 5-10%, comparing favorably with ELISA testing (CV = 10-15%).

# II. HOW IT WORKS

Array support	Samples		
* ****	Incubation of Sample with arrayed antibody supports	}	1-2 hrs
Biotin-Ab	Incubation with Biotinylated Ab	}	1-2 hrs
streptavidin	Incubation with labeled-Streptavidin	}	1 hrs
	Detection of signals		
	Data analysis and graph		

## III. COMPONENTS AND STORAGE

Store kit at ≤ -20 °C immediately upon arrival. Kit must used within the 6 month expiration date.

ITEM	COMPONENT	AAH-PTYR-3- 2	AAH-PTYR-3- 4	STORAGE TEMPERATURE AFTER THAWING**					
1	Antibody Arrays	2 membranes	4 membranes						
2	Blocking Buffer	3 vials (25ml/ea)	5 vials (25ml/ea)	≤-20 C					
3	Biotinylated Anti-Phosphotyrosine Antibody	1 vial	2 vials	2-8°C (for up to 3 days after dilution)					
4	1,000X HRP-Streptavidin Concentrate	1 vial (	50 µl)	2-8 °C					
5	20X Wash Buffer I Concentrate	1 vial (	30ml)						
6	20X Wash Buffer II Concentrate	1 vial (	30ml)						
7	2X Cell Lysis Buffer Concentrate	1 vial (1	16 ml)	2.8 °C					
8	Detection Buffer C	1 vial (10 ml)	2 vials (10 ml/ea)	270 C					
9	Detection Buffer D	1 vial (10 ml)	2 vials (10 ml/ea)						
10	Incubation Tray w/ Lid	2 trays	4 trays	Room Temperature					
11	Protease Inhibitor Cocktail 1 vial								
12	Phosphatase Inhibitor Cocktail II	1 v	ial	<u>2-20 C</u>					
Other Kit Components: Plastic Sheets, Array Map Template, User Manual									

\*Each package contains 2 or 4 membranes

\*\*For up to 3 months (unless stated otherwise) or until expiration date

# IV. ADDITIONAL MATERIALS REQUIRED

- Pipettors, pipet tips and other common lab consumables
- Orbital shaker or oscillating rocker
- Tissue paper, blotting paper or chromatography paper
- Adhesive tape or plastic wrap
- Distilled or de-ionized water
- A chemiluminescent blot documentation system:
  - o CCD Camera
  - X-Ray Film and a suitable film processor
  - o Gel documentation system
  - o Or another chemiluminescent detection system capable of imaging a western blot

# V. SAMPLE TIPS AND GENERAL CONSIDERATIONS

#### A. Sample Collection, Preparation, and Storage

**NOTE:** Optimal methods will need to be determined by each experimenter empirically based on researched literature and knowledge of the samples.

- If not using fresh samples, freeze samples as soon as possible after collection.
- Avoid multiple freeze-thaw cycles. If possible, sub-aliquot samples prior to initial storage.
- It is strongly recommended to add a protease inhibitor cocktail to cell and tissue lysate samples.
- Avoid sonication of 1 ml or less as this can quickly heat and denature proteins
- Most samples will not need to be concentrated. If concentration is required, a spin column concentrator with a chilled centrifuge is recommended.
- Always centrifuge the samples hard after thawing (~10,000 RPM for 2-5 minutes) in order to remove any particulates that could interfere with detection.
- The Cell Lysate can be prepared as follows:
  - For attached cells, remove supernatant from cell culture, wash cells twice with cold 1X PBS (for suspension cells, pellet the cells by spinning down the cells at 1500 rpm for 10 min) making sure to remove any remaining PBS before adding Lysis Buffer. Solubilize the cells at 2x10<sup>7</sup> cells/ml in 1X Lysis Buffer containing Protease Inhibitor Cocktail and Phosphatase Inhibitor Cocktail Set II (see preparation note shown on page 7 under Component Preparation Section). 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 14,000 x g for 10 min.

It is recommended that sample protein concentrations be determined using a total protein assay. For incubation with the Phosphorylation Antibody Array I, use at a protein concentration of 50-1000  $\mu$ g/ml for cell lysates.

Lysates should be used immediately or aliquot and stored at -70 °C. Thawed lysates should be kept on ice prior to use.

If you experience high background, you may further dilute your samples. If signals are too weak, the cell lysates can be pretreated by immunoprecipitations before incubation with array membranes. Immunoprecipitations can be done using anti-phosphotyrosine and protein A.

General tips for preparing lysate samples can be viewed on the online Resources page of the website.

#### B. Sample Types and Recommended Dilutions/Amounts

- **NOTE:** Optimal sample dilutions and amounts will need to be determined by each experimenter empirically but the below recommendations may be used as a starting point. Blocking Buffer (ITEM 2) should be used to dilute samples. Normalize by loading equal amounts of protein per sample.
- **Cell and Tissue Lysates**: load **50 to 1000 μg** of total protein (after at least a 5-fold dilution to minimize the effect of any detergent(s)). Therefore the original lysate concentration should be **250 μg to 5 mg/ml**.

#### C. Handling Membranes

- The antibody printed side of each membrane is marked by a dash (-) or number (#) in the upper left corner.
- Do not allow membranes to dry out during the experiment or they may become fragile and break OR high and/or uneven background may occur.
- Grasp membranes by the corners or edges only using forceps. DO NOT touch printed antibody spots.

#### D. Incubations and Washes

- Perform <u>ALL</u> incubation and wash steps under gentle rotation or rocking motion (~0.5 to 1 cycle/sec) using an orbital shaker or oscillating rocker to ensure complete and even reagent/sample coverage. Rocking/rotating too vigorously may cause foaming or bubbles to appear on the membrane surface which should be avoided.
- All washes and incubations should be performed in the Incubation Tray (ITEM 10) provided in the kit.
- Cover the Incubation Tray with the lid provided during all incubation steps to avoid evaporation and outside debris contamination.
- Ensure the membranes are completely covered with sufficient sample or reagent volume during each incubation.
- Avoid forceful pipetting directly onto the membrane; instead, gently pipette samples and reagents into a corner of each well.
- Aspirate samples and reagents completely after each step by suctioning off excess liquid with a pipette. Tilting the tray so the liquid moves to a corner and then pipetting is an effective method.
- Optional overnight incubations may be performed for the following steps to increase overall spot signal intensities:
  - Sample Incubation
  - Biotinylated Antibody Cocktail Incubation
  - HRP-Streptavidin Incubation
- **NOTE:** Overnight incubations should be performed at 4 °C (also with gentle rocking/shaking). Be aware that longer incubations can also increase the background response so complete liquid removal and washing is critical.

### VI. CHEMILUMINESCENCE DETECTION TIPS

- Beginning with adding the detection buffers and ending with exposing the membranes should take no more than 10-15 minutes as the chemiluminescent signals may start to fade at this point.
- Trying multiple exposure times is recommended to obtain optimum results.
- A few seconds to a few minutes is the recommended exposure time range, with 30 seconds to 1 minute being suitable for most samples.

## VII. COMPONENT PREPARATION

- **NOTE:** Thaw all reagents to room temperature immediately before use. If wash buffers contain visible crystals, warm to room temperature and mix gently until dissolved.
- **NOTE:** The Biotinylated Antibody Cocktail (ITEM 3) and the HRP-Streptavidin Concentrate (ITEM 4) vials should be briefly centrifuged (~1000 g) before opening to ensure maximum recovery and mixed well as precipitates may form during storage.

ITEM	COMPONENT	PREPARATION	EXAMPLE				
1	Antibody Arrays	No Proportion	NI/A				
2	Blocking Buffer	No Preparauori	N/A				
3	Biotinylated Antibody Cocktail*	Pipette 1 ml of Blocking Buffer into each vial. Mix gently with a pipette. Transfer the entire contents into a tube containing 11 ml of the Blocking Buffer.	N/A				
4	1,000X HRP-Streptavidin Concentrate	Dilute 1,000-fold with Blocking Buffer. Mix gently with a pipette.	10 μl of 1,000X concentrate + 9990 μl of Blocking Buffer = 10 ml of 1X working solution				
5	20X Wash Buffer I Concentrate	Dilute <b>each</b> 20-fold with distilled or deionized	10 ml of 20X concentrate + 190 ml of water				
6	20X Wash Buffer II Concentrate	water.	=200mlof1X working solution				
7	2X Cell Lysis Buffer Concentrate**	Dilute 2-fold with distilled or deionized water.	10mlof2Xconcentrate+10mlofwater= 20mlof1X working solution				
8	Detection Buffer C						
9	Detection Buffer D	No Preparation	N/A				
10	Incubation Tray w/ Lid						
11	Protease Inhibitor Cocktail	Pipette 60 µl of <b>1X</b> Cell Lysis Buffer into the vial to prepare 100X Protease Inhibitor Cocktail concentrate.					
12	Phosphatase Inhibitor Cocktail II	Add 180 μl of 1X Lysis Buffer into the vial to prepare 25X Phosphatase Inhibitor Cocktail Set II Concentrate. Dissolve the powder thoroughly by gentle mixing.					

 $^{*1}$  vial is enough to test 2 membranes

\*\*Only for use for preparing cell or tissue lysates. General tips for preparing lysates and other common sample types can be found on the online Resources Page

**Note**: Prior to preparing cell or tissue lysates: Add 20 μl Protease Inhibitor Cocktail Concentrate (100X) and 80 μl Phosphatase Inhibitor Cocktail Set II Concentrate (25X) into 1.9 ml 1X Lysis Buffer immediately before use. Mix well.

# VIII. PROTOCOL

- **NOTE:** Prepare all reagents and samples immediately prior to use. See Sections V and VII. <u>ALL</u> incubations and washes must be performed under gentle rotation/rocking (~0.5-1 cycle/sec)
- 1) Remove the kit from storage and allow the components to equilibrate to room temperature (RT).
- 2) Carefully remove the Antibody Arrays (ITEM 1) from the plastic packaging and place each membrane (printed side up) into a well of the Incubation Tray (ITEM 10). One membrane per well.

**NOTE:** The antibody printed side is marked by a dash (-) or number (#) in the upper left corner.

#### A. Blocking

- 3) Pipette 6 ml of Blocking Buffer (ITEM 2) into each well and incubate for 1 hour at RT.
- 4) Aspirate blocking buffer from each well with a pipette.

#### B. Sample Incubation

- 5) Pipette 6 ml of diluted or undiluted sample into each well and incubate for 1.5 to 5 hours at RT OR overnight at 4 °C.
- **NOTE:** Longer incubations can help maximize the spot signal intensities. However, doing so can also increase the background response so complete liquid removal and washing is critical.
- 6) Aspirate samples from each well with a pipette.

#### C. First Wash

- **NOTE:** The 20X Wash Buffer Concentrates I and II (ITEM 5 and 6) must be diluted 20-fold before use. See Section VII for details.
- 7) <u>Wash Buffer I Wash</u>: Pipette 20 ml of **1X** Wash Buffer I into each well and incubate for 5 minutes at RT. Repeat this 2 more times for a total of 3 washes using fresh buffer and aspirating out the buffer completely each time.
- 8) <u>Wash Buffer II Wash</u>: Pipette 20 ml of **1X** Wash Buffer II into each well and incubate for 5 minutes at RT. Repeat this 1 more time for a total of 2 washes using fresh buffer and aspirating out the buffer completely each time.

#### D. Biotinylated Antibody Cocktail Incubation

**NOTE:** The Biotinylated Antibody Cocktail (ITEM 3) must be prepared before use. See Section VII for details.

- 9) Pipette 6 ml of the **prepared** Biotinylated Antibody Cocktail into each well and incubate for 1.5 to 2 hours at RT OR overnight at 4°C.
- 10) Aspirate biotinylated antibody cocktail from each well.

#### E. Second Wash

11) Wash membranes as directed in Steps 7 and 8.

#### F. HRP-Streptavidin Incubation

- **NOTE:** The 1,000X HRP-Streptavidin Concentrate (ITEM 4) must be diluted before use. See Section VII for details.
- 12) Pipette 6 ml of **1X** HRP-Streptavidin into each well and incubate for 2 hours at RT OR overnight at 4°C.
- 13) Aspirate HRP-Streptavidin from each well.

#### G. Third Wash

14) Wash membranes as directed in Steps 7 and 8.

#### H. Chemiluminescence Detection

- **NOTE:** *Do not allow membranes to dry out during detection.*
- 15) Transfer the membranes, printed side up, onto a sheet of chromatography paper, tissue paper, or blotting paper lying on a flat surface (such as a benchtop).
- 16) Remove any excess wash buffer by blotting the membrane edges with another piece of paper.
- 17) Transfer and place the membranes, printed side up, onto a plastic sheet (provided) lying on a flat surface.
- **NOTE:** Multiple membranes can be placed next to each other and fit onto a single plastic sheet. Use additional plastics sheets if necessary.
- 18) Into a single clean tube, pipette equal volumes (1:1) of Detection Buffer C (ITEM 8) and Detection Buffer D (ITEM 9). Mix well with a pipette.
- **EXAMPLE**: 4.2 ml of Detection Buffer C + 4.2 ml of Detection Buffer D = 8.4 ml (enough for 2 membrane)
- 19) Gently pipette 4 ml of the Detection Buffer mixture onto each membrane and incubate for 2 minutes at RT (DO NOT ROCK OR SHAKE). <u>Immediately afterwards, proceed to Step 20</u>.
- **NOTE:** Exposure should ideally start within 5 minutes after finishing Step 19 and completed within 10-15 minutes as chemiluminescence signals will fade over time. If necessary, the signals can usually be restored by repeating washing, HRP-Streptavidin and Detection Buffers incubations (Steps 11-19)

- 20) Place another plastic sheet on top of the membranes by starting at one end and gently "rolling" the flexible plastic sheet across the surface to the opposite end to smooth out any air bubbles. The membranes should now be "sandwiched" between two plastic sheets.
- **NOTE:** Avoid "sliding" the top plastic sheet along the membranes' printed surface.
- 21) Transfer the sandwiched membranes to the chemiluminescence imaging system such as a CCD camera (recommended) and expose.
- **NOTE:** Optimal exposure times will vary so performing multiple exposure times is strongly recommended. See Section VI for additional details.

#### I. <u>Storage</u>

22) To store, without direct pressure, gently sandwich the membranes between 2 plastic sheets (if not already), tape the sheets together or use plastic wrap to secure them, and store at  $\leq$  -20 °C for future reference.

### IX. TYPICAL RESULTS



Typical results obtained with RayBio<sup>®</sup> C-Series Antibody Arrays

The preceding figures present typical images obtained with RayBio<sup>®</sup> C-Series Antibody Arrays. These membranes were probed with conditioned media from two different cell lines. Membranes were exposed with Kodak X-Omat<sup>®</sup> film at room temperature for 1 minute.

Note the strong signals of the Positive Control spots in the upper left and lower right corners. (See below for further details on the control spots.)

The signal intensity for each antigen-specific antibody spot is proportional to the relative concentration of the antigen in that sample. Comparison of signal intensities for individual antigen-specific antibody spots between and among array images can be used to determine relative differences in expression levels of each analyte sample-to-sample or group-to-group.

# X. INTERPRETING THE RESULTS

#### A. Control Spots

<u>Positive Control Spots (POS)</u> – controlled amount of biotinylated antibody printed onto the array. Used for normalization and to orientate the arrays.

<u>Blank Spots (BLANK)</u> – nothing is printed here. Used to measure the background response.

#### B. Data Extraction

Visual comparison of array images may be sufficient to see differences in relative protein expression. However, most researchers will want to perform numerical comparisons of the signal intensities (or more precisely, signal *densities*), using 2-D densitometry. Gel/Blot documentation systems and other chemiluminescent or phosphorescent detection systems are usually sold as a package with compatible densitometry software.

Any densitometry software should be sufficient to obtain spot signal densities from your scanned images. One such software program, ImageJ, is available for free from the NIH website along with an array plug-in.

We suggest using the following guidelines when extracting densitometry data from our array images:

- For each array membrane, identify a single exposure that the exhibits a high signal to noise ratio (strong spot signals and low background response). Strong Positive Control Spot signals but not too strong that that they are "bleeding" into one another is ideal. The <u>exposure time does not need to be identical for</u> <u>each array</u>, but Positive Control signals on each array image should have similar intensities.
- Measure the density of each spot using a circle that is roughly the size of one of the largest spots. Be sure to use the <u>same extraction circle</u> dimensions (area, size, and shape) for measuring the signal densities on every array for which you wish to compare the results.
- For each spot, use the <u>summed signal density</u> across the entire circle (ie, total signal density per unit area)

#### C. Data Analysis

**NOTE:** RayBiotech offers Microsoft<sup>®</sup> Excel-based Analysis Software Tools for each array kit for automatic analysis. Please visit the website at <u>www.raybiotech.com</u> or contact us for ordering information.

Once the raw numerical densitometry data is extracted, the background must be subtracted and the data normalized to the Positive Control signals to analyze.

<u>Background Subtraction</u>: Select values which you believe best represent the background. If the background is fairly even throughout the membrane, the Negative Control Spots (NEG) and/or Blank Spots (BLANK) should be similar and are accurate for this purpose.

<u>Positive Control Normalization</u>: The amount of biotinylated antibody printed for each Positive Control Spot is consistent from array to array. As such, the intensity of these Positive Control signals can be used to

normalize signal responses for comparison of results across multiple arrays, much like housekeeping genes and proteins are used to normalize results of PCR gels and Western Blots, respectively.

To normalize array data, one array is defined as "Reference Array" to which the other arrays are normalized to. The choice of the Reference Array is arbitrary.

**NOTE:** The RayBio<sup>®</sup> Analysis Software Tools always designate Array 1/Sample 1 as the Reference Array.

Next, the simple algorithm below can be used to calculate and determine the signal fold expression between like analytes.

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

Where:

P1 = mean signal density of Positive Control spots on reference array P(y) = mean signal density of Positive Control spots on Array "y" X(y) = mean signal density for spot "X" on Array for sample "y" X(Ny)= normalized signal intensity for spot "X" on Array "y"

For example:

Let's determine the relative expression for IL-6 on two different arrays (Arrays 1 and 2). Let's assume that the duplicate signals for the IL-6 spots on each array are identical (or that the signal intensity used in the following calculation is the mean of the two duplicates spots). Also assume the following:

P1 = 2500 P2 = 2700 IL-6 (1) = 300 IL-6 (2) = 455

Then IL-6(N2) = 455 \*2500/2700 = 421.30

The fold increase of IL-6(N2) vs IL-6(1) = 421.3/300 = 1.40-fold increase or a 40% increase in the signal intensity of IL-6 in Array 2 vs. Array 1.

# XI. 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	204	205	200	207	200	209	270	271	272	275	274	275	270	277	270	279	200	201	202	203	204	205	200	207	200	209	290	291	292	295
20	204	205	200	207	200	209	270	2/1	202	2/3	2/4	2/5	270	207	2/0	2/9	200	201	202	205	204	205	200	207	200	209	230	291	292	295
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
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	POS3	POS2	POS1																										
36	Blank	POS3	POS2	POS1																										

POS = Positive Control Spot BLANK = Blank Spot

**NOTE:** Protein alternative names, accession numbers, and official symbols can be accessed on <u>www.raybiotech.com</u> via the Resources Page.

# XII. ARRAY TARGET LIST

Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name	Number	Name
1	14-3-3 beta	73	Antithrombin III	145	C4BPA	217	CHREBP	289	Cytokeratin 9	361	EVC2	433	Glyoxalasell
2	14.2.2 opcilon	74	ADA	146	CED 0	210	Chromograpin P	200	D4 GDI	262	Errin	124	GM2A
2	14.3.2 sta	74		140	0.00 0	210	Chromogramm B	200	DAK	262	E21111	435	CME hate
3	14-3-3 eta	75	APLP-1	147	C6	219	Chromogranin C	291	DAK	363	FII	435	GIVIF Deta
4	14-3-3 gamma	76	APINI2	148	C8G	220	CIP29	292	Contactin-4	364	FABP5	436	GNB1
5	14-3-3 sigma	77	Apo (a)	149	C9orf40	221	СКВ	293	DARS2	365	Factor IX	437	GNPTG
6	14-3-3 theta	78	APOA1BP	150	CA1	222	CLIC1	294	DCI	366	Factor V	438	GOLPH2
7	14-3-3 zeta	79	ApoF	151	CA150	223	CLIC4	295	DCXR	367	Factor XII	439	GOLPH4
8	53BP1	80	ApoL1	152	CA2	224	CLIP170	296	DDAH1	368	Factor XIII	440	GOT2
9	67LR	81	ApoL2	153	CA3	225	CL-P1	297	DDT	369	FAM20C	441	GPR116
10	ABAT	82	ARFBP1	154	CACNB4	226	CLPS	298	DDX3Y	370	FAM3C	442	GPLD1
11	ABCF1	83	AREGEE3	155	CAD	227		299	DEEA6	371	Fascin	443	GRHL1
12	ABI3BD	8/		156	Cadherin 22	228	CNN2	300	DEP-1	372	EASN	444	Granzyme M
12	A01301	07	AraBE	150	Codhorin C	220	CNOT1	201	DNER	372	fact cholotal Mussin	444	CRUPP
15	ACAAI	65	ArgKS	157	Caulierin-6	229	CNUTI	301	DINER	3/3		445	GRHPR
14	ACAA2	86	ARP19	158	CALD1	230	CO4A2	302	Dermcidin	374	FASTKD5	446	GRP
15	ACACA	87	Arp2	159	CALML5	231	COG4	303	Desmocollin 1	375	FBP38	447	GSTM1
16	ACAA	88	ARP2/3	160	Calmodulin	232	COL19A1	304	Desmocollin-2	376	FBP2	448	GSTP1
17	ACLP	89	Arp3	161	Calpain 1	233	COL4A3	305	Desmocollin-3	377	FBPase 1	449	Guanylin
18	ACLY	90	ARPC2	162	Calpain S1	234	Col6A2	306	Desmoglein-1	378	FCGBP	450	GULP1
19	Aconitase 1	91	ARPC3	163	Calpastatin	235	COL9A3	307	Desmoglein-2	379	FDPS	451	H6PD
20	ACTBL2	92	ART3	164	Calretinin	236	COLEC10	308	Desmoplakin	380	FH	452	HABP2
21	ACTC1	93	ARTS1	165	Calumenin	237	Collagen La1	309	Desmuslin	381	Fibrillin 1	453	HB7
22	Actinin alpha 1	94	ARX	166	CAP1	238	Collagen III	310	Destrin	382	FGG	454	HCEC1
22		05		167	Cane	220	Collagon Waf	311	DCV	392	Fibringen like 2	154	HDGE
23	ADAIVIDECT	30	ASITZE	107	CAPZAA	2.59	Callana	242	DISCA	202	Fibringen-title 2	455	
24	ADAS	96	ASGR2	108	CAPZA1	240	Collagen IX	312	DISC 1	384	FIDTINOPEPTIDE B	456	HEG1
25	ADH1B	97	ASK1	169	CPB2	241	Collagen V	313	DMGDH	385	Fibulin 3	457	Hemoglobin
26	ADH1C	98	AST	170	CARHSP1	242	Collagen VI	314	DMRN9	386	Ficolin-2	458	Hemoglobin A1c
27	ADH4	99	DNPEP	171	Caspase-14	243	Collagen X	315	DBH	387	Filamin A	459	HBB
28	ADH5	100	ASXL1	172	Catalase	244	COL15A1	316	DOT1L	388	Filamin B	460	HBD
29	ADM	101	ATBF1	173	Cathelicidin	245	COMP	317	DPEP2	389	Filamin C	461	HBG2
30	Advillin	102	ATP5A	174	Cathepsin A	246	CFB	318	DPP3	390	FKBP12	462	HEXB
31	AFG3L2	103	ATP50	175	Cathensin G	247	Contactin-3	319	DPPI	391	FKBP25	463	HGFA
32	AGA	104	ATPR	176	Cathensin H	2/18	COPS8	320	DRU1	302	EKBD51	464	hGH
22	Aggreeen	105	DOCNTO	177	Cathopsin 7	240	Corneedeemeein	221	DECAM	202	FIC2	404	huppoh
33	Aggrecan	105	B3GINT2	177	Catriepsin 2	249	Corneodesmosti	321	DSCAIVI	393	FLG2	405	IIHR250
34	AGXI	106	B4Gall1	1/8	CBS	250	Coronin 3	322	DSPG3	394	FOLR3	466	HIBADH
35	AHNAK	107	B7-H2	179	CCDC126	251	Cortactin	323	Dystroglycan	395	Frizzled 8	467	HINT1
36	Ansp	108	B7-H3	180	CCDC25	252	COIL1	324	UBA1	396	FRY	468	HIP1R
37	AIF	109	BAD	181	CCT3	253	CPE	325	ECHS1	397	FSH	469	Histone H1.2
38	AK2	110	Band 3	182	CD109	254	CPEB3	326	ECM-1	398	Azurocidin	470	Histone H1.3
39	AKAP9	111	BASP1	183	CD133	255	CPM	327	EEF1G	399	FUCA1	471	Histone H2A
40	AKR1B1	112	Bassoon	184	CD155	256	CPN1	328	EEF2	400	FUCA2	472	Histone H2A.Z
41	AKR1C3	113	BAZ2B	185	CD157	257	CPNE3	329	EFEMP2	401	FAH	473	Histone H2B K
42	AKR7A2	114	BCHE	186	CD16	258	CPS1	330	EFTUD2	402	G0/G1switch 2	474	Histone H3.3
43	ALAD	115	Bcl-w	187	CD21	259	СКММ	331	FHD1	403	G3BP	475	Histone H4
44	ΔΙΤ	116	BCOR	188	CD32	260	CRE21	332	EHD3	404	GALNT2	476	HLA-C
45		117	hota 1 Sportrin	190	CD2E	260	CRUPP	222	ELESS	405	gamma Catonin	477	HMGR1
45	AD/1	110	CDVDD1	100	CD3014	201	Criti	224	elF4A1	405	CARDU	470	HIMODI
40	AUXI	110	CRIBBI	190	CD39L4	202	CIKL	334	err4A1	400	GAPUH	478	HIVIGB2
47	ALDH16A1	119	beta 1 lubulin	191	CD41	263	CRMP2	335	elf5A	407	GARNL1	479	HMGB3
48	ALDH1A1	120	CUBB3	192	CD42b	264	CRTAC1	336	ELAVL1	408	GART	480	HMGN2
49	ALDH9A1	121	BID	193	CD48	265	CS	337	EMILIN1	409	Gastrokine 1	481	HN1
50	ALKP	122	BIN2	194	CD5L	266	Ctip2	338	EMSY	410	GATM	482	FoxA1
51	ALP	123	BIRC6	195	CD9	267	Cux2	339	EN2	411	GBE1	483	hnRNP A1
52	MAN1A1	124	BLMH	196	CD98	268	Cyclophilin A	340	Endorepellin	412	GCDFP 15	484	hnRNP A2B1
53	alpha Actinin 4	125	BLVRB	197	CDA	269	Cyclophilin B	341	ENO1	413	GCLC	485	hnRNP C1+C2
54	Alpha Fodrin	126	BMP-1	198	CDC5L	270	Cystatin D	342	ENO1+ENO2+ENO3	414	GCSH	486	hnRNP G
55	alpha Glucosidase II	127	BPGM	199	CDK2	271	Cystatin E	343	ENSA	415	GDA	487	hnRNP L
56	alpha-Synuclein	128	BPIFB1	200	CEACAM-8	272	Cystatin S	344	Envoplakin	416	GDF7	488	hnRNP M1-M4
57	alpha Tubulin	120	RDI 1	201	CECP1	272	Cystatin CN	3/5	EDN	A17	GDI1	490	hnRND II
50		120	PPCAD	201	CENIDE	273	CCDD1	240	EDD 44	410	GDI1	400	Hornerin
58	CRYAA	130	DRCA Z	202	CENPF	2/4	CSKP1	340	EP 641	418	GUI2	490	nomerin
59	ALS	131	BRD2	203	CEP57	275	CYTL1	347	EPCR	419	Gepnyrin	491	Hoxb3
60	Als2	132	Brevican	204	CES1	276	Cytochrome b5	348	Ephrin B1	420	GFAP	492	HOXD11
61	ALS2CR1	133	Brg1	205	CETP	277	Cytochrome c (n)	349	Ephrin B2	421	GHRF	493	HP1BP3
62	Aminoacylase	134	BRSK1	206	Cezanne	278	Cytokeratin 1	350	EPHX2	422	GIP	494	HPD
63	Androgen Receptor	135	BTD	207	CFHR1	279	Cytokeratin 10	351	EPPK1	423	GLIPR2	495	HPR
64	ANGPTL6	136	BTF3	208	CFHR4	280	Cytokeratin 13	352	Eps15	424	GLRX1	496	HPRT
65	ANGPTL8	137	C1q	209	CFHR5	281	Cytokeratin 14	353	ERAB	425	G6PD	497	HRG
66	ANK	138	C1nA	210	CFI	282	Cytokeratin 15	354	ERAP2	426	PRKCSH	498	HRSP12
67	Ankrd26	139	C1oB	211	CFI 1	283	Cytokeratin 16	355	ERp29	427	GLUD1	499	HSC70
60	Annevin A1	1/0	C1cP1	212	CEVII	200	Cytokeratin 17	355	ERp57	120	0.001	500	HSP47
60	Annovin A2	1.41	C101	212	CI (17	204	Cutokoratin 20	257	EPr 70	420	65704	500	115647
70	Annexin Az	141	CIKL	213	Chitada	285	Cytokeratin 20	357	ERp/2	429	03101		h
/0	Annexiñ Ab	142	U1S	214	Chitobiase	286	Cytokeratin 3	358	ESD	430	655		
71	Annexin V	143	ELP6	215	Cnitotriosidase	287	Cytokeratin 4	359	ESR1	431	GPD1		
72	ANP	144	C4.4A	216	CHORDC1	288	Cytokeratin 5	360	ETL	432	Glycoprotein V		i

# XIII. TROUBLESHOOTING GUIDE

PROBLEM	CAUSE	RECOMMENDATION						
	Chemiluminescent imager is not working properly	Contact image manufacturer						
No signals	Too Short Exposure	Expose the membranes longer						
(not even the positive controls spots)	Degradation of components due to improper storage	Store entire kit at ≤ - 20°C. Do not use kit after expiration date. See storage guidelines.						
	Improper preparation or dilution of the HRP-Streptavidin	Centrifuge vial briefly before use, mix well, and do not dilute more than 1000-fold						
	Waiting too long before exposing	The entire detection process should be completed in 10-15 minutes						
Positive controls spots	Low sample protein levels	Decrease sample dilution, concentrate samples, or load more protein initially						
signals visible but no other	Skinned Sample Incubation Sten	Samples must be loaded after the blocking step						
spots	Too Short of Incubations	Ensure the incubations are performed for the appropriate time or try the optional overnight incubation(s)						
	Bubbles present on or below membrane	Don't rock/rotate the tray too vigorously or pipette the sample or reagent with excessive force						
Uneven signals and/or	Insufficient sample or reagent volume	Load enough sample and reagent to completely cover the membrane						
background	Insufficient mixing of reagents	Gently mix all reagents before loading onto the membrane, especially the HRP-Streptavidin and Biotin Antibody Cocktail						
	Rocking/Rotating on an uneven surface while incubating	Rock/rotate on a flat surface or the sample or reagent can "pool" to one side						
	Too much HRP-Streptavidin or Biotinylated Antibody Cocktail	Prepare these signal enhancing components precisely as instructed						
	Membranes dried out	Do not let the membranes dry out during the experiment. Cover the incubation tray with the lid to minimize evaporation						
High background signals or all spots visible	Too High of Sample Protein Concentration	Increase dilution of the sample or load less protein						
• • • • •	Exposed Too Long	Decrease exposure time						
	Insufficient Washing	Ensure all the wash steps are carried out and the wash buffer is removed completely after each wash step						
	Non-specific binding	Ensure the blocking buffer is stored and used properly.						

RayBio<sup>®</sup> is the trademark of RayBiotech, Inc.

The RayBio<sup>®</sup> Cytokine Antibody Array C-Series is 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 6 months from the date of purchase 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.

Kodak X-Omat is a registered trademark of Eastman Kodak Company. Microsoft Excel is a registered trademark of Microsoft Corporation. Note:

This product is for research use only.



©2010 RayBiotech, Inc.