

Pathway oriented phospho-proteomics analysis with reverse phase protein array (RPPA) technology in cell growth and cancer therapy

Takeomi Inoue, Naoko Iwata, Eiji Nishiwaki, Yasuyuki Kirii
Carna Biosciences, Inc., Kobe, Japan

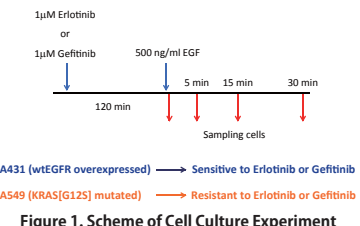
Abstract

Reverse Phase Protein Array (RPPA) is a powerful technology that facilitates simultaneously measuring the level of a single protein in multiple samples in a high-throughput format. We have developed a signal transduction oriented RPPA system that focuses on phospho-protein profiles in samples. Thus far, our system utilizes 180 phospho-protein specific antibodies and can profile the phosphorylation status of more than 100 Gene Ontology (GO) processes covering more than 10 components in each process. In this study, we show the phospho-protein profiles of two cancer cell lines, A431 and A549 treated with Erlotinib and Gefitinib. Our RPPA system could detect dynamic changes in phosphorylation profiles of both cell lines as well as several differences between them with high sensitivity and reproducibility.

On the other hand, the results of our RPPA experiments were in accordance with those of the *in vitro* kinase assays we have obtained. We are now expanding the panel of cell lines and drugs to devise a method to guide strategies for effective targeting in drug discovery by integrating kinase-inhibition-potency data and cellular-signaling-profile data.

Methods

Cell Culture: A431 (human epidermoid carcinoma) and A549 (human non-small cell lung carcinoma) cells were incubated in serum-free DMEM or F-12 medium, respectively overnight prior to treatment with TKIs. Cells were pre-incubated with 1μM Erlotinib or Gefitinib for 2 hours before stimulating with 500 ng/ml rhEGF. Treated cells were collected at 0 (immediately after adding EGF), 5, 15, and 30 min (Figure 1).



RPPA: 1-4 x 10⁶ cells were homogenized in our proprietary lysis buffer. Lysates were printed onto glass slides with an arrayer equipped with 32 pins. Serially diluted lysates were also spotted in order to place the expression level of samples in a dynamic range for signal detection (Figure 2). Each slide was probed with an anti-phospho antibody and signals were generated by employing a catalyzed signal amplification (CSA) system (Dako) and Alexa-647-conjugated streptavidin. The targets of anti-phospho antibodies utilized in this study span a wide variety of Gene Ontology categories and pathways (Table 1 & 2).

Data Analysis: Stained slides were scanned on a desktop scanner (InnoScan 710AL, INNOPSYS, Carbonne, France). SuperCurve (<http://bioinformatics.mdanderson.org/main/OOMPA:Overview>) algorithms were used to estimate a single value of logarithmic concentration from the four serial dilutions for each sample. Two sample t-test was used to compare values between two groups. All analyses were performed using the R statistical Programming Environment (v.2.15.1).

Pathway Mapping: UniProt ID was ascribed to each target of an anti-phospho antibody based on the antibody information, while the records in UniProt were linked to the information in KEGG pathway and Gene Ontology. This combination (Figure 3) resulted in classification of our targets into GO groups (Table 1) and mapping onto KEGG pathways (Table 2, Figure 4).

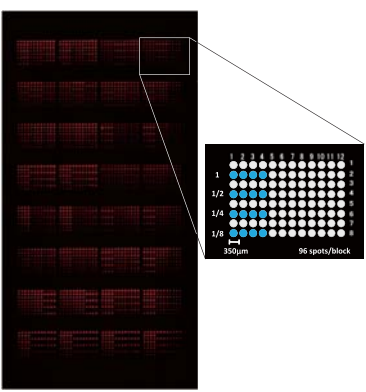


Figure 2. Stained Slide Image and Array Design
A representative stained glass slide image and array design (inset) are shown. Four two-fold serial dilutions of lysates were printed in quadruplicate. Neighboring blocks have the same layout to reduce the effects of local background variation. Each glass slide can accommodate at most 96 samples that are arrayed as two blocks of 192 lysate spots that each contains 6 samples.

10	14	71	40	GO:0000001	protein kinase C activity	P
11	14	41	40	GO:0001003	stress-activated MAPK cascade	P
12	14	41	40	GO:0006267	protein kinase receptor signaling pathway	P
13	14	40	40	GO:0004036	regulation of blood vessel endothelial cell migration	P
14	14	40	40	GO:0001008	MAPK stress kinase activity	P
15	14	37	40	GO:0006691	JAK-STAT signaling involved in growth hormone signaling pathway	P
16	14	37	40	GO:0001363	transforming-factor-mediated signaling pathway	P
17	14	32	40	GO:0000007	cytokine-cytokine receptor signaling pathway	P
18	14	32	40	GO:0003066	TGF-beta type II receptor signaling pathway	P
19	14	31	40	GO:0001003	cell-to-cell signaling by receptor activity	P
20	14	31	40	GO:0001010	the receptor	P
21	14	30	40	GO:0001034	cell-to-cell receptor 2 signaling pathway	P
22	14	29	40	GO:0001010	the receptor	P
23	14	28	40	GO:0000075	glycoprotein tyrosine kinase receptor signaling pathway	P
24	14	28	40	GO:0001010	the receptor	P
25	14	27	40	GO:0000063	tyrosine phosphorylation	P
26	14	27	40	GO:0000063	trans signaling	P
27	14	27	40	GO:0001042	cell-to-cell receptor 1 signaling pathway	P
28	14	26	40	GO:0001010	the receptor	P
29	14	26	40	GO:0001038	positive regulation of tyrosine autophosphorylation	P
30	14	26	40	GO:0004030	regulation of fatty acid oxidation	P
31	14	25	40	GO:0001010	signal transduction	P
32	14	25	40	GO:0005983	cell receptor signaling pathway	P
33	14	25	40	GO:0004005	transforming-factor-mediated signaling pathway	P
34	14	24	40	GO:0001032	positive regulation of protein kinase C activity	P
35	14	24	40	GO:0004983	protein kinase of intracellular cell cycle	P
36	14	23	40	GO:0004034	flattened growth factor receptor signaling pathway	P
37	14	21	40	GO:0000086	insulin receptor signaling pathway	P
38	14	21	40	GO:0000073	RBM protein (L1 protein) receptor specific DNA binding	P
39	14	21	40	GO:0000173	epinephrine growth factor receptor signaling pathway	P
40	14	20	40	GO:0004941	protein kinase B signaling cascade	P
41	14	20	40	GO:0001036	activation of MAPK activity	P
42	14	20	40	GO:0004977	protein autophosphorylation	P