

Proteomic Analysis of Tumor Biomarkers in Human Clinical Specimens Using the BioOdyssey™ Calligrapher™ MiniArrayer

Nere Navarro and Phillip E Schwartz,
Protein Biotechnologies, Inc., Ramona, CA 92065 USA

Introduction

Reverse-phase protein microarray (RPPA) technology is a powerful emerging analytic strategy for interrogating the proteomes of tissues and cells. As a high-throughput screening platform, RPPA permits rapid quantitative identification of novel cancer biomarkers associated with oncogenesis and disease progression. RPPA should be useful in accelerating the understanding of cellular differentiation, transformation, angiogenesis, tumorigenesis, and metastasis. This new technology has the ability to quickly elucidate alterations in protein expression levels, detect posttranslational modification and mRNA processing events, and dissect molecular networks associated with drug administration or exposure to environmental factors (for example, toxins, infectious agents, or radiation). With this new ability, cancer researchers, clinicians, and, more importantly, patients should realize a significant benefit in biomarker discovery, molecular diagnostics, drug development, and personalized medicine.

Innovations that increase the production, quality, and performance of protein microarrays are critical to realizing the full potential of this screening platform. Laboratories engaged in high-throughput protein biomarker identification and discovery require protein microarray products and methods that are easy to analyze, are easy to automate, and ensure consistent results. The amount of information obtained from current protein microarray experiments is limited to what one probe and slide can generate, and thus these experiments can consume an excessive amount of expensive resources, reagents, and clinical tissue samples. Realizing the full scientific and healthcare benefits of RPPA will require that sample throughput and sensitivity be maximized. Significant cost reductions can be achieved by using innovative multiplex detection systems to probe multiple samples with multiple analytes in a single assay.

In our laboratory, we have developed a line of ready-to-use protein microarrays for cancer research. Arrays are fabricated using lysates prepared from an extensive collection of human tumor and normal clinical specimens representing a variety of

organs, tissues, and cancer subtypes. Our approach is similar to that used to evaluate protein expression in the NCI-60 cancer cell lines (Nishizuka et al. 2003) and biopsy samples in clinical trial research (Gulmann et al. 2005). We are also developing sensitive detection systems that would permit automated multiplex analysis of multiple targets in a single assay. The development and manufacturing of our products require a robust and consistent arraying platform with exceptional spot-tracking features for quality control and run validation.

Detection sensitivity is an inherent challenge in the use of protein microarrays, especially for the detection of specific proteins in complex biological samples such as cell or tissue lysates. Spot sizes are often small (~100–300 μm), and the amount of target protein available for detection is correspondingly reduced. The choice of detection methods (chromogen, chemiluminescence, fluorescence, etc.) also influences the detection limits of the assay. Finally, data capture and image analysis hardware and software vary greatly in sensitivity, flexibility, and of course cost. Therefore, to maximize the data obtained from an experiment, it is important to maximize sensitivity.

We wished to evaluate parameters affecting the use of protein microarrays and the quality of the data obtained from experiments using them. To determine specificity and sensitivity criteria in our system, we prepared a special batch of microarray slides using lysates from lung, breast, and colon tissue. Slides were interrogated with individual antibodies specific for known biomarkers involved in oncogenesis, vascularization, the extracellular matrix, and cell maintenance. Target visualization was performed using a chromogen detection system, and microarray images were acquired and processed.

Methods

Six pairs of patient-matched normal and tumor tissue protein extracts from colon, breast, and lung were spotted in triplicate on specially prepared nitrocellulose-coated glass slides (Grace Bio-Labs, Inc.). Arrays were printed using a BioOdyssey Calligrapher miniarrayer with TeleChem Stealth SMP6 micro spotting pins, producing 160 μm spots, delivering ~2.0 nL/spot of a 1.0 mg/mL protein solution. Lysates were spotted in

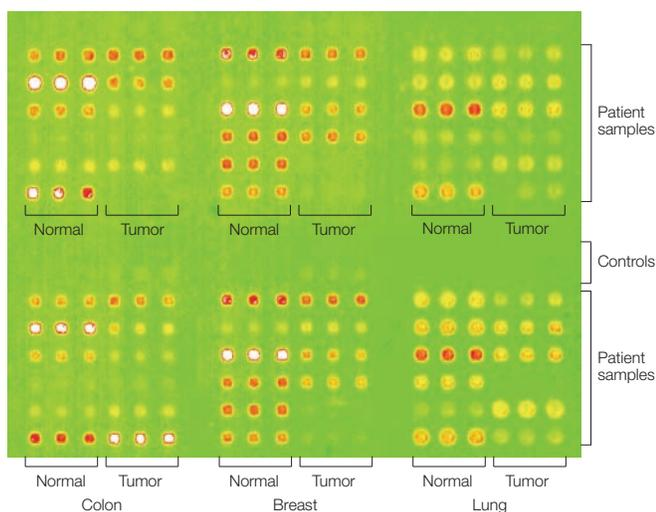


Fig. 1. False-color image of protein levels in a microarray of human clinical specimens. Normal and tumor lysates from six patients were arrayed and probed with an antibody to a circulating protein thought to correlate with differences in vascularization. Each row in a grid represents samples from a single patient. The first three spots in a row are from normal cells and the last three are from tumor cells. Upper and lower grids are duplicates of the same patient samples. The top two rows of the bottom grid consist of negative control spots to assess background noise of the assay.

modified RIPA protein extraction buffer according to the manufacturer's recommended protocols. Positive (purified IgGs) and negative (buffer only and BSA) controls were also included in the arrays. Sample tracking and spot placement were validated using the BackTracker™ file system, which is built into the miniarrayer software.

Slides were treated using protocols similar to those for western blotting of proteins transferred to nitrocellulose membranes (Krajewski et al. 1996). Monoclonal and polyclonal primary antibodies raised against β -actin, tubulin (Imgenex Corp.), glyceraldehyde-3-phosphate dehydrogenase and laminin-1 (Proteus BioSciences, Inc.), human IgG, and rabbit IgG (Jackson ImmunoResearch Laboratories, Inc.) were used according to suppliers' recommendations and detected using HRP-conjugated secondary antibodies (Protein Biotechnologies, Inc.). Color development was achieved by incubating the slides in tetramethylbenzidine (TMB) peroxidase substrate (Sigma-Aldrich Co.), and the slides were subsequently scanned using an ArrayIt SpotWare colorimetric microarray scanning system (TeleChem International, Inc.).

Results and Discussion

Protein arrays were used to evaluate several criteria associated with the manufacture and use of this platform as a tool for cancer research. After slides were spotted, labeled, and imaged, false-color images were generated to reveal differential protein target content within each matched pair of samples. Figure 1 shows the differential abundance of a circulating protein marker in normal and tumor samples derived from multiple tissues and multiple patients. For example, in breast tumors, all normal samples displayed higher levels of the protein than the tumor samples. Spot size and deposition were within acceptable limits (10% CV) as was signal-to-noise ratio for colorimetric detection of protein biomarkers using immunological detection methods. Sensitivity using TMB was in the low picogram range.

In our laboratory, the BackTracker software feature that comes with the BioOdyssey Calligrapher miniarrayer adds a unique advantage over those offered in other array systems we have tried. Because we have to create our own source plates for the different protein array products we manufacture, the ability to easily predefine the robot printing program for each unique array layout requirement, which the BackTracker file system provides, is especially valuable.

This study demonstrates the successful use of the BioOdyssey Calligrapher miniarrayer for accurately and reproducibly printing protein arrays of complex biological samples derived from human tumors and normal tissues. The BackTracker software feature enhances our ability to rapidly redefine protein array templates and has significantly improved our product manufacturing capabilities. Rapid screening of multiple clinical specimens for protein biomarkers in a single assay in an easy-to-use platform not only is possible but also is a valuable first step in determining further analytical strategies.

References

- Gulmann C et al., Proteomic analysis of apoptotic pathways reveals prognostic factors in follicular lymphoma, *Clin Cancer Res* 11, 5847–5855 (2005)
- Krajewski S et al., Detection of multiple antigens on western blots, *Anal Biochem* 236, 221–228 (1996)
- Nishizuka S et al., Proteomic profiling of the NCI-60 cancer cell lines using new high-density reverse-phase lysate microarrays, *Proc Natl Acad Sci USA* 100, 14229–14234 (2003)

ArrayIt, SpotWare, and Stealth are trademarks of TeleChem International, Inc.

Information in this tech note was current as of the date of writing (2006) and not necessarily the date this version (rev A, 2006) was published.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 4BIORAD Australia 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400
Canada 905 712 2771 China 86 21 6426 0808 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65
Germany 089 318 84 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 455 8800 India 91 124 4029300/5013478 Israel 03 963 6050
Italy 39 02 216091 Japan 03 5811 6270 Korea 82 2 3473 4460 Mexico 55 5200 05 20 The Netherlands 0318 540666 New Zealand 64 9415 2280
Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 095 721 14 04 Singapore 65 6415 3188 South Africa 27 0861 246 723
Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189/2578 7241 United Kingdom 020 8328 2000