

## Profiling of Human Angiogenesis Biomarkers in Sera of Cancer Patients Using the Bio-Plex<sup>®</sup> Suspension Array System

V. Gupta, L. Ma, J. Fedynyshyn, and W. Tan,  
Bio-Rad Laboratories, Inc., Hercules, CA 94547 USA

### Introduction

Angiogenesis is defined as the formation of new blood vessels from preexisting blood vessels. This process is considered natural during wound healing, pregnancy, and the female reproductive cycle. Angiogenesis becomes pathological during early tumor formation. Angiogenic transformation is a multistep process consisting of extracellular matrix remodeling, endothelial cell migration and proliferation, capillary differentiation, and abnormal production and interaction of various angiogenic factors. These processes subsequently lead to tumor vasculature and metastasis. A variety of angiogenic factors (soluble as well as membrane bound) involved in regulating these steps have been identified. The levels of these factors could be elevated or reduced in cancer patients. These angiogenic factors are potential candidate drug targets relevant for the development of cancer therapies. Development of cancer therapies often involves the measurement of levels of these angiogenic factors.

In this study, we evaluated the ability of the Bio-Plex suspension array system to detect the levels of nine biomarkers — angiopoietin-2 (Ang-2), VEGF, PDGF-BB, leptin, PECAM-1, IL-8, follistatin, HGF, and G-CSF in sera collected from four different types of cancer patients. These nine biomarkers were chosen for their potential clinical relevance to cancer development (Bach et al. 2007, Bauhofer et al. 2007, Grusch et al. 2006, Jackson et al. 2003, Schips et al. 2007, Shih et al. 2006, Somasundar et al. 2004, Vairaktaris et al. 2007, Wang et al. 2006, Yoshiji et al. 2005).

### Methods

A magnetic bead-based technology was used to determine profiles of the nine angiogenic biomarkers. This technology utilizes an 8  $\mu\text{m}$  magnetic bead coupled with two fluorophores. These classification dyes exhibit emission spectra at distinct wavelengths and absorption maxima at 635 nm. The reporter dye is a third fluorophore, phycoerythrin (PE), which shows emission at a third distinct wavelength and maximum absorption at 532 nm. PE was chosen for its high molar extinction coefficient, quantum yield, resistance to photobleaching, minimal self-quenching, and excellent stability. The detector unit consists of a flow cell that enables magnetic beads to travel in a single file (laminar flow) through a region illuminated by a pair of lasers. The beads emit light at three wavelengths, two from the classification dyes and one from the reporter dye.

Profiles of the nine angiogenic biomarkers collected from sera of four types of cancer patients, colorectal cancer ( $n = 10$ ), breast, lung, and prostate cancer ( $n = 15$ ), were evaluated using a magnetic bead and flow-based assay on a Bio-Plex suspension array system. This platform integrates a series of color-coded beads, each of which is coupled to a unique antibody specific for a biochemical marker. These capture antibody-coupled beads serve as solid phases for the capture of desired serum or plasma proteins, using a standard sandwich-based detection format, as schematically represented in Figure 1. The levels of biomarkers in cancer patients were compared to those in control sera, collected from healthy individuals ( $n = 117$ ).

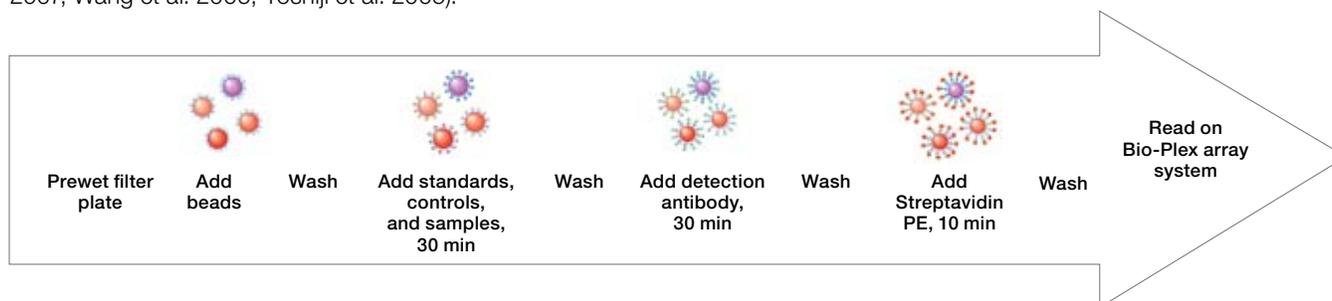


Fig. 1. A schematic representation of a multiplex bead-based sandwich immunoassay for the experimental workflow.

## Results

### Assay Specificity

To study assay specificity, the nine recombinant antigens were independently evaluated with multiplexed capture and detection antibodies. This format determines the overall target specificity that is conferred by the capture antibody. Results are summarized in Table 1. Data showed <1% cross-reactivity for 8 of the 9 targets. The antibody pair for Ang-2 exhibited <3.4% cross-reactivity with IL-8, VEGF, PDGF-BB, and HGF antigens. Percentages of cross-reactivity were calculated using median fluorescence intensities (MFIs) of the multiplexed capture beads and detection antibodies in the presence of a single antigen.

### Assay Sensitivity and Limit of Detection

In a study designed to evaluate assay sensitivity and limits of detection, the standard curve for each target was tested in duplicate from eight independent assays for serum matrix and from five independent assays for RPMI matrix. The limit of detection (LOD) is defined as the concentration obtained at background MFI for the standard curve. Figure 2 depicts LOD values for both serum and RPMI matrices. Table 2 shows the range of concentrations in which the assay is both precise (intra-assay  $\leq 15\%$  CV and inter-assay  $\leq 25\%$  CV) and accurate (70–130% standard and spike recovery). These parameters

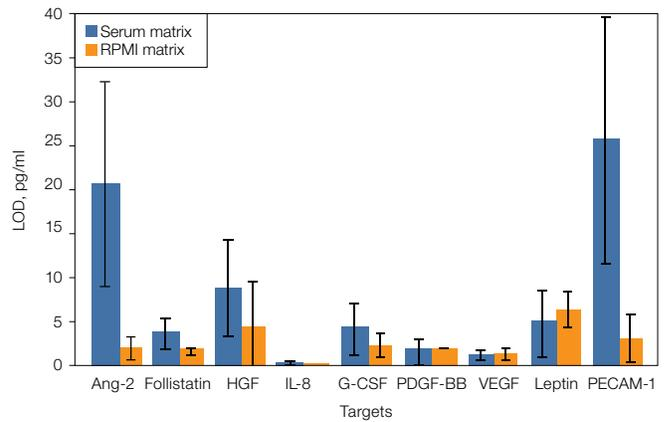


Fig. 2. Assay sensitivity of targets on serum and RPMI matrices.

define the upper and lower limits of quantitation for each target. The standard curves for each target are shown in Figure 3.

### Precision of Measurements

Values were determined from the variance of MFI of replicate wells for eight standard points. The intra-assay precision was measured from five replicate wells in one assay and the inter-assay precision was measured from duplicate wells in five independent assays. All targets recorded  $\leq 15\%$  CV.

Table 1. Percentage of cross-reactivity for different angiogenic factors.

| Target      | Ang-2 | Follistatin | HGF   | IL-8 | G-CSF | PDGF-BB | VEGF  | Leptin | PECAM-1 |
|-------------|-------|-------------|-------|------|-------|---------|-------|--------|---------|
| Ang-2       | 110.9 | 0.1         | 0.2   | 0.8  | 0.1   | 0.0     | 0.0   | 0.1    | 0.2     |
| Follistatin | 1.5   | 112.5       | 0.2   | 0.7  | 0.1   | 0.0     | 0.0   | 0.2    | 0.2     |
| HGF         | 3.4   | 0.4         | 104.5 | 0.9  | 0.1   | 0.1     | 0.1   | 0.3    | 0.2     |
| IL-8        | 2.4   | 0.2         | 0.2   | 98.8 | 0.0   | 0.0     | 0.0   | 0.5    | 0.3     |
| G-CSF       | 0.0   | 0.1         | 0.1   | 0.7  | 107.5 | 0.0     | 0.0   | 0.0    | 0.1     |
| PDGF-BB     | 3.3   | 0.3         | 1.1   | 0.9  | 0.2   | 106.1   | 0.3   | 1.4    | 0.9     |
| VEGF        | 1.2   | 0.2         | 0.3   | 0.8  | 0.1   | 0.0     | 106.5 | 0.2    | 0.2     |
| Leptin      | 0.3   | 0.0         | 0.0   | 0.7  | 0.0   | 0.0     | 0.0   | 116.4  | 0.0     |
| PECAM-1     | 0.2   | 0.1         | 0.0   | 0.7  | 0.0   | 0.0     | 0.0   | 0.0    | 117.0   |

Table 2. Assay range, as determined from the mean of five independent assays.

| Targets     | Assay Range, pg/ml |        |
|-------------|--------------------|--------|
|             | LLOQ*              | ULOQ** |
| Ang-2       | 59                 | 59,784 |
| Follistatin | 25                 | 25,584 |
| HGF         | 20                 | 20,201 |
| IL-8        | 1                  | 7,475  |
| G-CSF       | 11                 | 25,665 |
| PDGF-BB     | 3                  | 24,423 |
| VEGF        | 1                  | 16,906 |
| Leptin      | 8                  | 36,256 |
| PECAM-1     | 45                 | 32,691 |

\* Lower limit of quantitation.

\*\* Upper limit of quantitation.

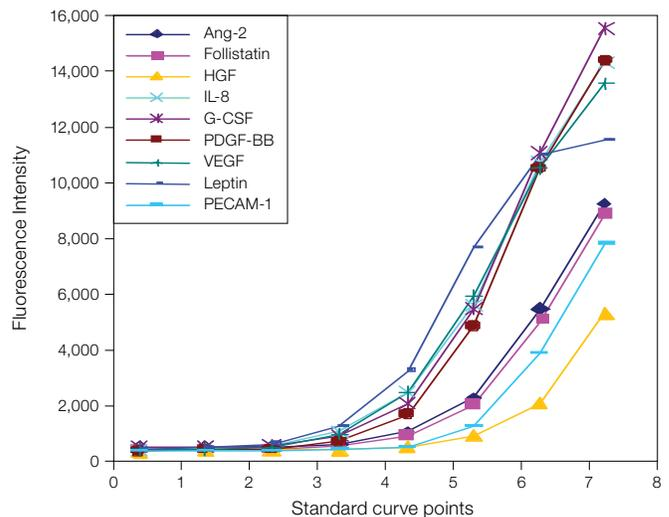
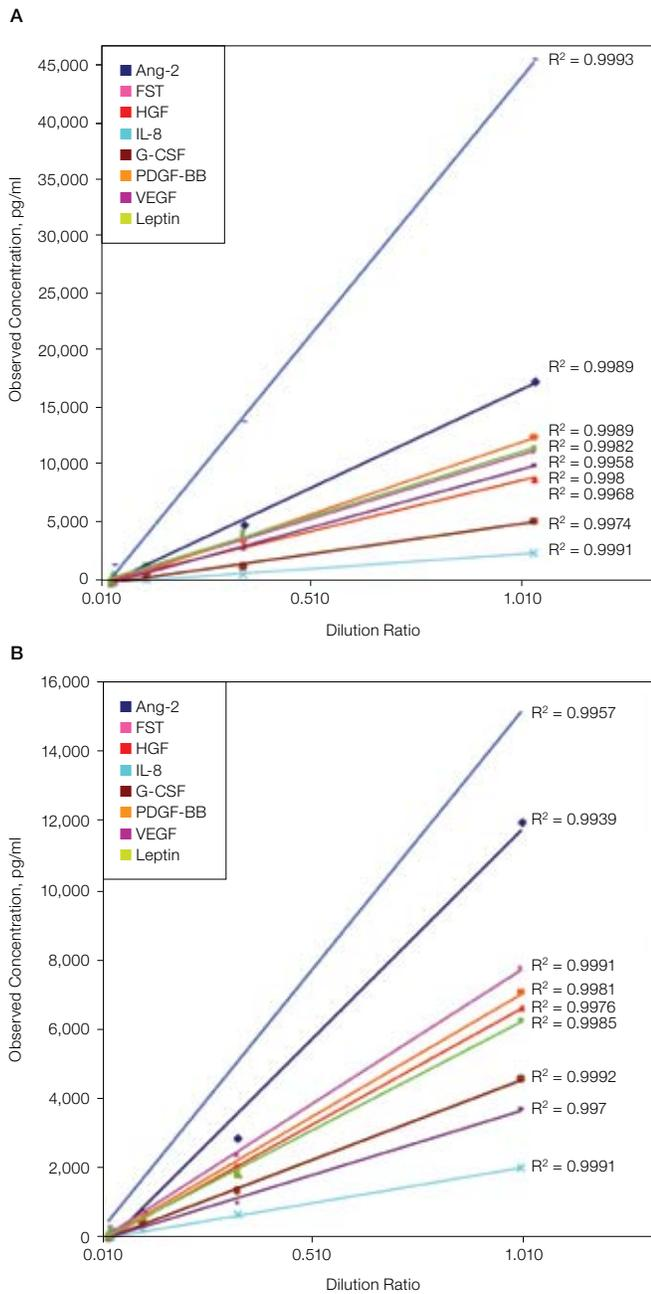


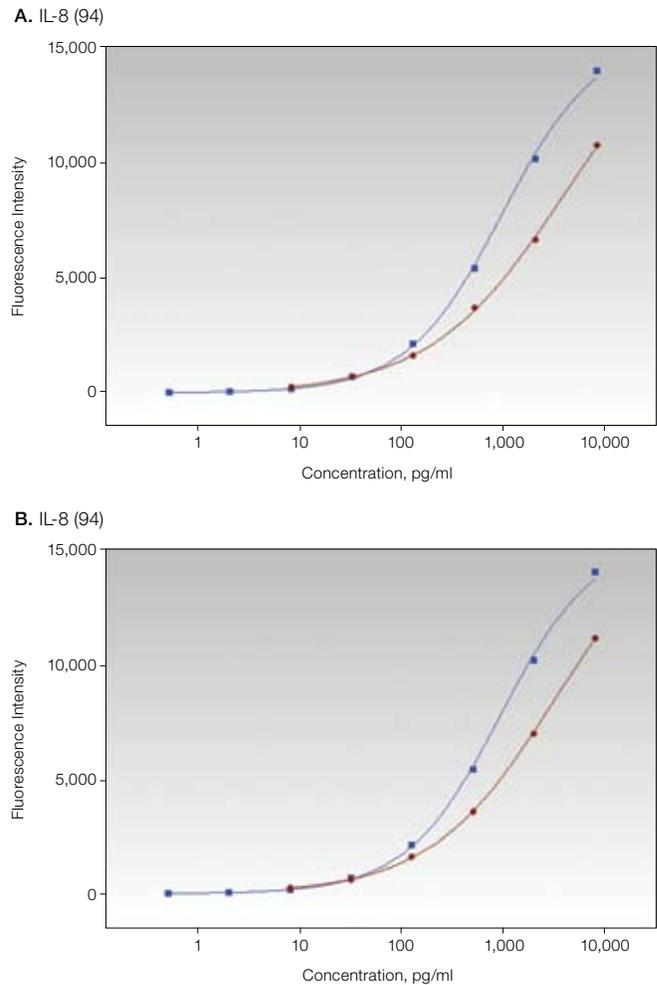
Fig. 3. Plot of 8-point standard curves for each target.



**Fig. 4. Linearity of dilution in different targets.** A, concentration of samples measured from serum matrix; B, concentration of samples measured from a plasma matrix.

#### Dilutional Linearity and Parallelism

Dilutional linearity ensures analyte present in concentrations above the ULOQ can be diluted linearly and measured accurately within the assay range. Linearity was demonstrated in both serum and plasma samples with  $R^2 > 0.99$  for all nine targets. Figure 4 shows one representative serum and plasma sample tested by means of spiked sampling.

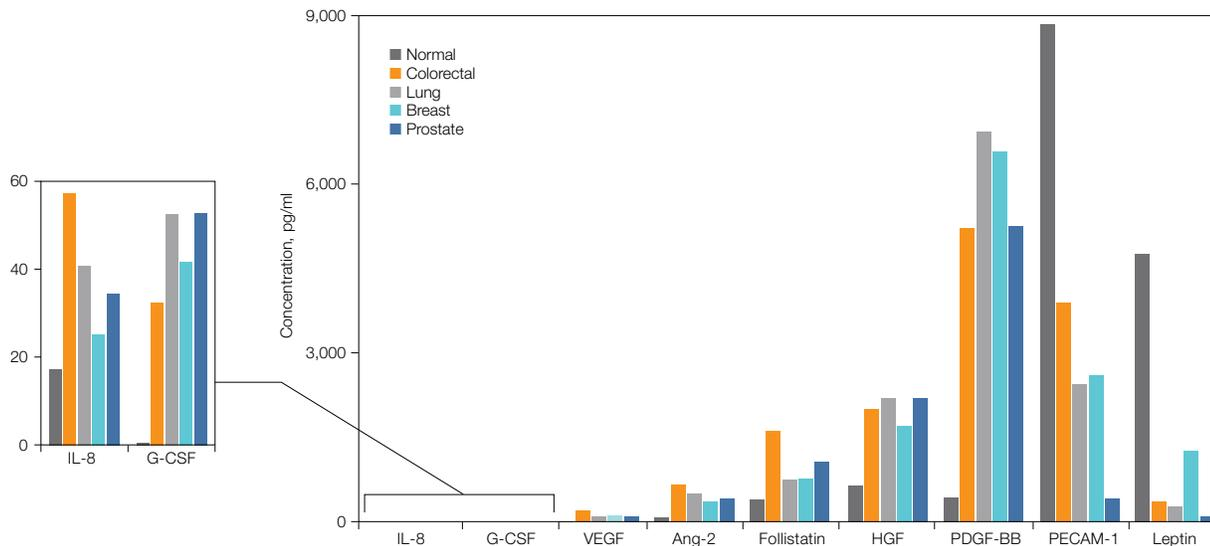


**Fig. 5. Examples of parallelism in serum and plasma matrices using an 8-point standard (■) and 6-point spiked serum or plasma sample (●) dilutions with 4-PL curve fitting.**

Parallelism ensures that the standards and samples have the same binding characteristics. This is demonstrated by comparing a spiked sample dilution-response curve with the standard concentration-response curve. Figure 5 shows parallelism between an 8-point standard and a 6-point spiked dilution using a representative serum or plasma sample with 4-PL curve fitting. With the exception of PECAM-1, the slope difference was  $<20\%$  for the remaining 8 targets in the serum and plasma samples tested.

#### Physiological Levels of Angiogenesis Biomarkers

The physiological levels of these markers were determined in 117 serum samples collected from apparently healthy donors. The mean and median values were compared to serum levels of samples collected from four cancer types. Seven of the nine targets showed marked elevation across all four cancer types. In contrast, the levels of leptin and PECAM-1 were reduced (Figure 6).



**Fig. 6. Changes in biomarker levels in sera of different cancer groups.** Concentration levels represent median values.

### Conclusion

Initial screenings of cancer biomarkers are both costly and labor-intensive. In many cases, studies generated from different test formats often yield conflicting, and therefore, inconclusive results. The Bio-Plex system is capable of measuring the levels of multiple targets simultaneously in less than 2 hours, using as little as 12.5 µl of serum, plasma, and other matrices. This feature significantly reduces the time and cost spent on preliminary screening of serum samples for biomarker profiling. Overall, the assays are specific, sensitive, and reproducible. The system also provides a broad range for measurements, which can be utilized depending on specific sensitivity requirements.

Preliminary analysis of different cancer samples using the Bio-Plex suspension array system showed marked elevation of 7 of the 9 biomarkers studied, suggesting the importance of these markers as key regulators of angiogenesis. The reduction of the levels of leptin and PECAM-1 also provide interesting insight into the modulation of these markers in the associated cancers.

### References

- Bach F, et al. Angiotensins in malignancy. *EJSO* (2006) doi: 10.1016: 1–9.
- Bauhofer A, et al. Perioperative prophylaxis with granulocyte colony-stimulating factor (G-CSF) in high-risk colorectal cancer patients for an improved recovery: A randomized, controlled trial. *Surgery* (2007) 501–502.
- Grusch M, et al. Deregulation of the activin/follistatin system in hepatocarcinogenesis. *J. Hepatology* (2006) 45: 673–680.
- Jackson D E, et al. The unfolding tale of PECAM-1. *FEBS Letters* (2003) 540: 7–14.
- Shih A H, et al. Platelet-derived growth factor (PDGF) and glial tumorigenesis. *Cancer Letters* (2006) 232: 139–147.
- Schips L, et al. Serum levels of vascular endothelial growth factor (VEGF) and endostatin in renal cell carcinoma patients compared to a control group. *European Urology* (2007) 51: 168–174.
- Somasundar P, et al. Leptin is a growth factor in cancer. *J. Surgical Research* (2004) 116: 337–349.
- Vairaktaris E, et al. The interleukin-8 (-251A/T) polymorphism is associated with increased risk for oral squamous cell carcinoma. *EJSO* (2007) 33: 504–507.
- Wang P, et al. Bladder cancer cell invasion is enhanced by cross-talk with fibroblasts through hepatocyte growth factor. *Urology* (2007) 69: 780–783.
- Yoshiji H, et al. Angiotensin 2 displays a vascular endothelial growth factor dependent synergistic effect in hepatocellular carcinoma development in mice. *Gut* (2005) 54(12): 1768–1775.

The Bio-Plex suspension array system includes fluorescently labeled microspheres and instrumentation licensed to Bio-Rad Laboratories, Inc. by the Luminex Corporation.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA 800 4BIORAD Australia 61 02 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 21 3237 9400  
Canada 905 364 3435 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 Israel 03 963 6050  
Italy 39 02 216091 Japan 03 6361 7000 Korea 82 2 3473 4460 Mexico 52 555 488 7670 The Netherlands 0318 540666 New Zealand 0508 805 500  
Norway 23 38 41 30 Poland 48 22 331 99 99 Portugal 351 21 472 7700 Russia 7 495 721 14 04 Singapore 65 6415 3188 South Africa 27 861 246 723  
Spain 34 91 590 5200 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan 886 2 2578 7189 United Kingdom 020 8328 2000