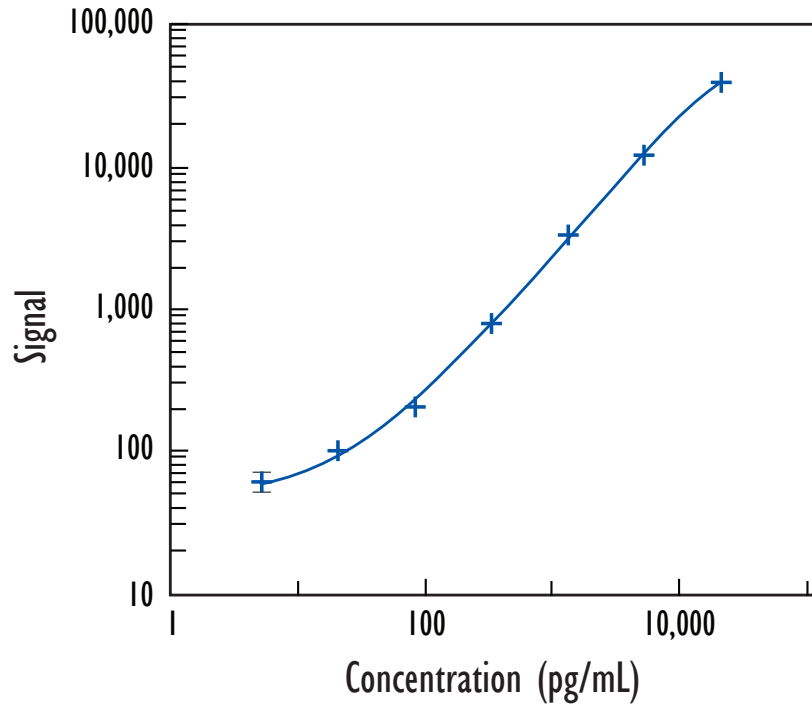


# MULTI-ARRAY<sup>®</sup> Human VEGF Tissue Culture Assay

## Detection of Human Vascular Endothelial Growth Factor (VEGF) in Tissue Culture

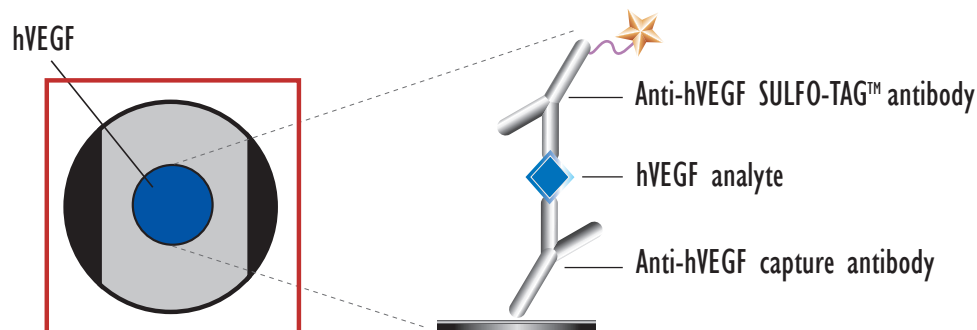


| Concentration (pg/mL) | hVEGF  |     |
|-----------------------|--------|-----|
|                       | Mean   | %CV |
| 0                     | 44     | 9   |
| 6                     | 56     | 17  |
| 24                    | 92     | 4   |
| 98                    | 192    | 4   |
| 391                   | 766    | 7   |
| 1,563                 | 3,324  | 2   |
| 6,250                 | 13,370 | 10  |
| 25,000                | 41,523 | 2   |

Above is representative calibration curve data.

|              | hVEGF    |
|--------------|----------|
| Average LLOD | 19 pg/mL |

LLOD is defined as 2.5x stdev above the background.



| Kit Size         | Catalog Number |
|------------------|----------------|
| 1 plate          | K151BMB-1      |
| 5 plates         | K151BMB-2      |
| 20 plates        | K151BMB-3      |
| 20 plates (Base) | K151BMA-3      |