

# Chemiluminescent VEGF ELISA Using the SpectraMax L Microplate Luminometer

SPECTRAMAX APPLICATION NOTE #9



By Cathy Olsen, Ph.D., Molecular Devices, Inc., 1311 Orleans Drive, Sunnyvale, CA 94089.

## INTRODUCTION

Vascular endothelial growth factors (VEGFs) are a family of secreted polypeptides that have been implicated in mammalian vascular development and in disease processes involving abnormal blood vessel growth. VEGFs are expressed during embryogenesis, where deletion of even a single VEGF allele severely disrupts vasculogenesis and is embryonic lethal. VEGF<sub>165</sub> is the most abundant and biologically active isoform of VEGF found in mammals.<sup>1</sup> The QuantiGlo Chemiluminescent VEGF Immunoassay is a solid phase ELISA that measures VEGF<sub>165</sub> levels in cell culture supernatants, serum, plasma, saliva, and urine. It uses the quantitative sandwich enzyme immunoassay technique in which a monoclonal antibody specific for VEGF is coated onto a microplate and standards and samples are added to the wells.<sup>2</sup> Unbound material is washed away, and a horseradish peroxidase-linked polyclonal antibody-specific for VEGF is added to the wells. An additional series of washes removes unbound antibody-enzyme reagent, then an enhanced luminol/peroxide substrate is added to the wells. (See Figure 1.) Light produced in proportion to the amount of VEGF initially bound to the wells is then measured using a microplate luminometer.

Assay results were detected using the SpectraMax<sup>®</sup> L microplate luminometer. This instrument is suitable for both flash and glow applications and is compatible with 96- and 384-well microplate formats. Its dedicated luminescence optical design yields an extremely high signal-to-noise ratio and extremely low crosstalk, with a dynamic range of more than eight orders of magnitude. Similar results may be obtained using the luminescence mode of SpectraMax<sup>®</sup> M5/M5<sup>e</sup> and FlexStation<sup>®</sup> 3 multi-mode microplate readers. Data collection and calculations were performed using SoftMax<sup>®</sup> Pro software, which includes preconfigured protocols to simplify data acquisition and analysis.

## MATERIALS

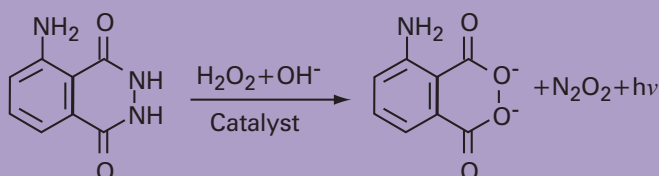
- QuantiGlo Human VEGF Immunoassay (R&D Systems P/N QVE00B). This kit contains all of the reagents and materials required to perform the assay, including assay standards and a microplate pre-coated with a monoclonal antibody specific for VEGF.
- SpectraMax L microplate luminometer. Note: this assay can also be detected using the luminescence mode of SpectraMax M5/M5<sup>e</sup> and FlexStation 3 multi-mode microplate readers.

## METHODS

### Preparation of reagents and standards

- Step 1. All reagents were brought to room temperature before use.
- Step 2. Wash Buffer Concentrate was diluted 10-fold with deionized water.
- Step 3. Working Glo Reagent was prepared 15 minutes to 4 hours before use by combining 1 part Glo Reagent A and 2 parts Glo Reagent B in a capped plastic tube, protected from light.
- Step 4. The VEGF standard was reconstituted with 0.5 mL deionized water and allowed to sit for a minimum of 15 minutes, with gentle agitation prior to making dilutions. Working standards were prepared by making a serial 1:10 dilution of the stock standard in

### Oxidation of Luminol (Figure 1)



An enhanced luminol/peroxide substrate is oxidized, resulting in emission of light.

Calibrator Diluent RD5L. Concentrations of working standards ranged from 20,000 to 1.3 pg/mL.

#### Assay procedure

- Step 1. 150  $\mu$ L of Assay Diluent RD1-8 was pipetted into each well.
- Step 2. 50  $\mu$ L of working standard or blank (Calibrator Diluent) was pipetted into triplicate wells of the supplied microplate. The microplate was incubated for 2 hours at room temperature with shaking (500  $\pm$  50 rpm recommended).
- Step 3. Wells were aspirated and washed 4 times with 400  $\mu$ L Wash Buffer using a multichannel pipettor. After the last wash, the remaining Wash Buffer was aspirated and the plate was blotted against clean paper towels.
- Step 4. 200  $\mu$ L of VEGF Conjugate was added to each well, and a new adhesive strip was applied. The microplate was incubated for 3 hours at room temperature, with shaking as in step 2.
- Step 5. Aspiration and washing steps were repeated as in step 3.
- Step 6. 100  $\mu$ L Working Glo Reagent was added to each well. The microplate was incubated for 5-20 minutes at room temperature, protected from light.
- Step 7. RLU values were determined using the SpectraMax L microplate luminometer. Instrument settings included the following: 1 second integration time, AutoRange PMT setting, and 470 nm target calibration wavelength.

#### RESULTS

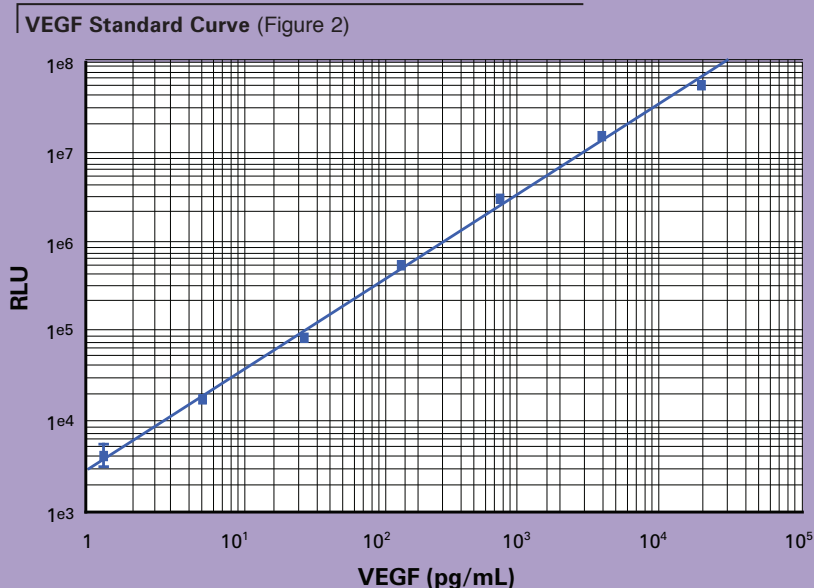
Figure 1 shows the VEGF standard curve obtained with the SpectraMax L microplate luminometer using the settings outlined above. The dynamic range for this assay spans greater than four orders of magnitude, with a calculated sensitivity of about 1.5 pg/mL, based on a concentration giving a signal 3 times the standard deviation of the background. This correlated well with R&D Systems' claim of a mean minimum detectable dose of 3.3 pg/mL for the kit. The SpectraMax M5 and FlexStation 3 multi-mode microplate readers yielded similar results (data not shown). Results were analyzed and the standard curve was plotted using SoftMax Pro software.

#### SUMMARY

The Quantiglo Human VEGF Immunoassay is a highly sensitive method for assaying VEGF in a variety of sample types. The SpectraMax L microplate luminometer with SoftMax Pro software provides excellent sensitivity and dynamic range for this assay as well as simplified data acquisition and analysis.

#### REFERENCES

1. Holmes, David IR and Zachary, Ian (2005). The vascular endothelial growth factor (VEGF) family: angiogenic factors in health and disease. *Genome Biology* 6: 209.
2. Quantiglo Human VEGF Immunoassay (P/N QVE00B) package insert.



Standards ranging from 20,000 to 1.3 pg/mL were assayed as described in the assay procedure. The microplate was read on a SpectraMax L microplate luminometer with a 1-second integration time and target calibration wavelength of 470 nm. Results were plotted using SoftMax Pro software.

#### SALES OFFICES

**United States & Canada**  
Molecular Devices  
Tel. +1-800-635-5577  
Fax +1-408-747-3601

**Brazil**  
Molecular Devices Brazil  
Tel. +55-11-3616-6607  
Fax +55-11-3616-6607

**China**  
Molecular Devices Shanghai  
Tel. +86-21-6887-8820  
Fax +86-21-6887-8890

**Germany**  
Molecular Devices GmbH  
Tel. +49-89/96-05-88-0  
Fax +49-89/9-62-02-34-5

**Japan**  
Molecular Devices Japan, Osaka  
Tel. +81-6-6399-8211  
Fax +81-6-6399-8212

Molecular Devices Japan, Tokyo  
Tel. +81-3-5282-5261  
Fax +81-3-5282-5262

**South Korea**  
Molecular Devices Korea, LLC  
Tel. +82-2-3471-9531  
Fax +82-2-3471-9532

**United Kingdom**  
Molecular Devices Ltd.  
Tel. +44-118-944-8000  
Fax +44-118-944-8001

[www.moleculardevices.com](http://www.moleculardevices.com)

FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

The trademarks used herein are the property of Molecular Devices, Inc. or their respective owners.

©2010 Molecular Devices, Inc. Printed in U.S.A. 6/10 #0120-14628