

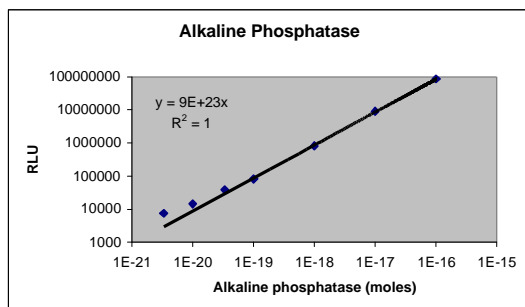
# A Veritas™ Microplate Luminometer Method for Alkaline Phosphatase



## 1. INTRODUCTION

The Turner BioSystems' Veritas™ Microplate Luminometer has a large dynamic range that is highly suited for alkaline phosphatase assays. Alkaline phosphatase is a stable enzyme often conjugated to secondary antibodies for immunoassays. Alkaline phosphatase is also a useful genetic reporter and a helpful tool in DNA cloning.

The superior sensitivity of the permits detection of very low levels of alkaline phosphatase activity. The Veritas can detect as little as  $3 \times 10^{-21}$  moles alkaline phosphatase enzyme. Measurements are linear over four orders of magnitude (Figure 1). All tests were conducted using Lumigen's APS-5 (Lumigen, Southfield, MI) and purified calf intestinal mucosa alkaline phosphatase (Biozyme, San Diego, CA).



*Figure 1: Detection of alkaline phosphatase with the Veritas Microplate Luminometer. 90  $\mu$ L APS-5 was added to 10  $\mu$ L alkaline phosphatase diluted in dH<sub>2</sub>O. Samples were incubated at room temperature for 5 minutes before measurement.*

## 2. MATERIALS REQUIRED

- Veritas™ Microplate Luminometer (P/N 9100-000)
- 96-well plates, white (E&K Scientific EK-25075)
- Lumigen APS-5
- Alkaline phosphatase (Biozyme ALP112G)
- p200 pipette and pipette tips
- p20 pipette and pipette tips

## 3. EXPERIMENT PROTOCOL

### 3.1 Reagent Preparation

**Lumigen APS-5:** Use as supplied. Store at 2–8°C, where it is stable for up to 12 months.

**Alkaline phosphatase:** Store at 4°C.

### 3.2 Instrument Setup

3.2.1 Double-click on the Veritas icon to start the software.

3.2.2 Click on "Create New Protocol" from the "Welcome to Veritas" dialog box.

3.2.3 Using the new protocol wizard, set up a protocol with 0 injectors, 1 second integration and 5 minute delay. Select the wells you wish to read and click on "Finish".

3.2.4 Enter your information into the "Experiment", "Operator", "Plate No.", and "Notes" fields in the "Main Dialog Box".

### 3.3 Sample Analysis

3.3.1 Equilibrate alkaline phosphatase, samples and APS-5 to room temperature.

3.3.2 Prepare a 10-fold serial dilution of alkaline phosphatase in your assay buffer.

3.3.3 Add 10  $\mu$ L of sample or standard to the 96-well plate.

3.3.4 Add 90  $\mu$ L of APS-5 reagent to the sample or standard.

**Note:** The recommended ratio of sample volume to APS-5 volume is 1:10 and should not exceed 1:1

3.3.5 Insert the plate into the Veritas and click on "Start" to begin assay. After the programmed 5-minute delay, the Veritas will begin measurement. RLU values measured by the Veritas will appear in the Excel spreadsheet after all the selected wells in each row have been read. If you encounter an error message, refer to the troubleshooting guide in the Veritas Operating Manual for more information.

**Note:** It is not recommended to open another Excel spreadsheet while the Veritas reads your sample plate.

3.3.6 Once the measurements are complete, you can access Excel to analyze your data.

3.3.7 Please make sure to remove your plate after measurement.

### 4. ABOUT LUMIGEN

Orders for Lumigen's products may be placed by:

Phone: (248) 351-5600  
Fax: (248) 351-0518  
E-mail: [marketing@lumigen.com](mailto:marketing@lumigen.com)  
Internet: [www.lumigen.com](http://www.lumigen.com)

Mailing address:

Lumigen, Inc.  
22900 West Eight Mile  
Southfield, MI 48034  
USA

### 5. ABOUT BIOZYME

Orders for Biozyme's products may be placed by:

Phone: (800) 423-8199  
Fax: (858) 549-0138

E-mail: [bioinfo@biozyme.com](mailto:bioinfo@biozyme.com)

Web Site: [www.biozyme.com](http://www.biozyme.com)

Mailing address:

Biozyme Laboratories International Ltd.  
9939 Hibert Street, Suite 101  
San Diego, CA 92131  
USA

### 6. ABOUT TURNER BIOSYSTEMS

Veritas is a trademark of Turner BioSystems.

Orders for Turner BioSystems' products may be placed by:

Phone: (408) 636-2400 or  
Toll Free: (888) 636-2401 (US and Canada)  
Fax: (408) 737-7919

Web Site: [www.turnerbiosystems.com](http://www.turnerbiosystems.com)

E-Mail: [sales@turnerbiosystems.com](mailto:sales@turnerbiosystems.com)

Mailing Address:

Turner BioSystems, Inc.  
645 N. Mary Avenue  
Sunnyvale, CA 94085 USA