## **Multilabel** Detection

# Photometric Assays with EnVision Multilabel Plate Reader



#### Introduction

Photometric assays are widely used for biochemical or cell based studies. Here we show on examples, NADH and Methyl-red, the possibility of performing those studies using absorbance mode on an EnVision<sup>®</sup> Multilabel Plate Reader.

NADH is a biomolecule, which participates in a number of biochemical reactions. Therefore it is widely studied among researchers.

Methyl-red is commonly used as a pH indicator. If the pH of the solution is less than 4.2, methyl-red appears as red. If the pH of the solution is above 6.3, it appears as yellow. This indicator can be used to follow, e.g., the bacteria medium growth.

#### **EnVision in photometric applications**

EnVision plate reader is an instrument with modular design, supporting all of the following technologies: AlphaScreen®/AlphaLISA®, luminescence flash and glow, fluorescence intensity, time-resolved fluorescence, fluorescence polarization and absorbance (UV/Visible). Absorbance measurements can be performed in the range of 240-850 nm with a speed of less than 3 min per 1536 well plate.

In absorbance measurements the light is directed through a light guide to the bottom-measuring head where the intensity of the light is measured using a reference photodiode. The light is then directed through the bottom of the plate and focused into the sample. The signal is detected with a photodiode, which is located above the plate.



The absorbance value is calculated by the equation

$$A = -\log(I/I_o)$$

where  $I_o$  is the light intensity without any sample and I is the intensity after an absorbing or reflecting medium.

### **Methods**

Dilution series of beta-NADH (Sigma-Aldrich<sup>®</sup>, N8129) was prepared in TSA, pH 7.4 buffer. The samples were pipetted into a black and clear 1536-well plate (Greiner<sup>®</sup>, 782096) (7.5  $\mu$ L/well), 6 replicates. The plate was measured with the EnVision instrument using a photometric filter for 340 nm (PerkinElmer, 2100-5190).

Dilution series of methyl-red was prepared in PBS-buffer. Samples were pipetted into a Greiner<sup>®</sup> black and clear 1536well plate (7.5  $\mu$ L/well), 6 replicates. The plate was measured with the EnVision instrument using a photometric filter for 492 nm (PerkinElmer, 2100-5220).

Dilution series of BSA was prepared into PBS buffer. Samples were pipetted into a 96-well plate (Corning<sup>®</sup>, 3615), 200  $\mu$ L/well. The plate was measured with the EnVision instrument using a photometric filter for 280 nm (PerkinElmer, 2100-5350).

### Label properties

Excitation filter	Absorbance 340/492/280
Number of flashes	3 flashes for 340 nm 100 flashes for 492 nm and 280

Number of flashes per A/D Conversion 1

#### Conclusion

In the two examples of NADH and Methyl-red, we show that the EnVision microplate reader is ideally suited to run photometric assays. Its optical range from 240-850 nm and the wide dynamic range of 4 OD combined with its short measurement times provide maximum flexibility, sensitivity and speed.







*Figure 2.* Methyl-red dilution series measured with EnVision, 100 flashes/well, abs. 492 nm, 1536-well plate.



*Figure 3.* BSA dilution series measured with EnVision, 100 flashes/ well, abs. 280 nm, 96-well plate.

#### PerkinElmer, Inc. 940 Winter Street Waltham, MA 02451 USA P: (800) 762-4000 or (+1) 203-925-4602 www.perkinelmer.com



For a complete listing of our global offices, visit www.perkinelmer.com/ContactUs

Copyright ©2002-2012, PerkinElmer, Inc. All rights reserved. PerkinElmer<sup>®</sup> is a registered trademark of PerkinElmer, Inc. All other trademarks are the property of their respective owners