

## Amplite™ Fluorimetric D-Lactate Assay Kit

### \*Red Fluorescence\*

Ordering Information	Storage Conditions	Instrument Platform
Product Number: 13810 (200 assays)	Keep in freezer Avoid exposure to light	Fluorescence microplate readers

### Introduction

Lactic acid is chiral and has two optical isomers: L-lactic acid and D-lactic acid. Lactate is constantly produced from pyruvate via the enzyme lactate dehydrogenase (LDH) in the process of metabolism and exercise. Monitoring lactate levels is a good way to evaluate the balance between tissue oxygen demand and utilization and is useful when studying cellular and animal physiology. D-lactate is not metabolized by mammals and its elimination from the body depends mainly on renal excretion. D- and L-lactic acid are found in many fermented milk products such as yoghurt and cheese, and also in pickled vegetables, and cured meats and fish. The D- and L-lactic acid content is a quality indicator of foods, such as egg, milk, fruit juice and wine. Abnormal high concentration of D-lactate in the blood is usually a reflection of bacterial overgrowth in the gastrointestinal tract.

AAT Bioquest's Amplite™ Lactate Assay Kits (Cat# 13814 and 13815 for L-lactate assay, and Cat# 13810 and 13811 for D-lactate assay) provide both fluorescence and absorbance-based method for detecting either L-lactate or D-lactate in biological samples such as serum, plasma, urine, as well as in cell culture samples. In the enzyme coupled assay, lactate is proportionally related to NADH, which is specifically monitored by a fluorogenic NADH sensor. The signal can be read by a fluorescence microplate reader at Ex/Em = 540 nm/590 nm. With this Fluorimetric Amplite™ D-Lactate Assay Kit, we were able to detect as little as 1.4 μM D-lactate in a 100 μL reaction volume. It is robust, and can be readily adapted for a wide variety of applications that require the measurement of D-lactate.

### Kit Components

Components	Amount
Component A: Enzyme mix	2 bottles (lyophilized powder)
Component B: Assay Buffer	1 bottle (10 mL)
Component C: NAD	1 vial
Component D: D-Lactate Standard	2.25 mg/vial

### Assay Protocol for One 96-Well Plate

#### Brief Summary

**Prepare D-lactate assay mixture (50 μL) → Add D-lactate standards or test samples (50 μL) → Incubate at room temperature for 30 min ~ 2 hours → Monitor fluorescence increase at Ex/Em = 540/590 nm**

*Note: Thaw one vial of each kit component at room temperature before starting the experiment.*

#### 1. Prepare NAD stock solution (100X):

Add 100 μL of H<sub>2</sub>O into the vial of NAD (Component C) to make 100X NAD stock solution.

#### 2. Prepare D-lactate stock solution:

Add 200 μL of H<sub>2</sub>O or 1xPBS buffer into the vial of D-lactate Standard (Component D) to make 100 mM D-lactate standard solution.

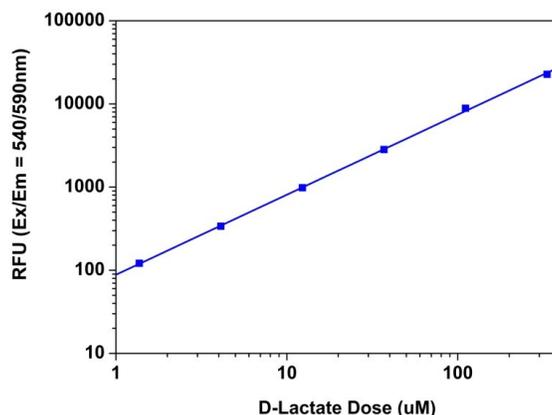
*Note: The unused D-lactate standard stock solution should be divided into single use aliquots and stored at -20°C.*



## Data Analysis

The fluorescence in blank wells (with the dilution buffer only) is used as a control, and is subtracted from the values for those wells with the D-lactate reactions. A typical D-lactate standard curve is shown in Figure 1.

*Note: The fluorescence background increases with time, thus it is important to subtract the fluorescence intensity value of the blank wells for each data point.*



**Figure 1.** D-lactate dose response was measured with Amplite™ Fluorimetric D-Lactate Assay Kit in a 96-well black plate using a Gemini (Molecular Devices) microplate reader. As low as 1.4  $\mu\text{M}$  D-lactate in 100  $\mu\text{L}$  volume can be detected with 1 hour incubation.

## References

1. McLellan A.C., et. al, *Analytical Biochemistry*, 1992, 206(1), 12-16.
2. Beaver W. L., et.al, Improved detection of lactate threshold during exercise using a log-log transformation, *Physiology*, 1985, 59 (6),1936-1940 .
3. Gérson F de Souza, et.al, Lactic acid levels in patients with chronic obstructive pulmonary disease accomplishing unsupported arm exercises, *Chronic Respiratory Disease*, 2010 7:(2) 75-82.
4. Garner H. E., et.al, Lactic acidosis: a factor associated with equine laminitis, *Journal of Animal Science*, 1977, 45:1037-1041.
5. Gladden, L.B. Lactate metabolism: A new paradigm for the third millenium. 2004, *J Physiol* **558(1)** 5-30.
6. Aguirre M., et.al, Lactic acid bacteria and human clinical infection, *Journal of Applied Microbiology*, 1993, 75 (2), 95-107.

**Warning: This kit is only sold to end users. Neither resale nor transfer to a third party is allowed without written permission from AAT Bioquest®, Inc. Chemical analysis of the kit components is strictly prohibited. Please call us at 408-733-1055 or e-mail us at [info@aatbio.com](mailto:info@aatbio.com) if you have any questions.**