

Amplite™ Fluorimetric L-Lactate Assay Kit

Catalog number: 13814

Unit size: 200 Tests

Component	Storage	Amount
Component A: Enzyme Mix	Freeze (<-15 °C), Dessicated, Avoid Light	2 bottles (lyophilized powder)
Component B: Assay Buffer	Freeze (<-15 °C), Avoid Light	1 bottle (10 mL)
Component C: NAD	Freeze (<-15 °C), Dessicated, Avoid Light	1 vial
Component D: L-Lactate Standard	Freeze (<-15 °C), Dessicated, Avoid Light	2.25 mg/vial

OVERVIEW

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 (generated by bacteria) 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™ L-Lactate Assay Kit, we were able to detect as little as 1.4 μM L-lactate in a 100 μL reaction volume.

AT A GLANCE

Protocol summary

1. Prepare L-lactate working solution (50 μL)
2. Add L-lactate standards or test samples (50 μL)
3. Incubate at room temperature for 30 min - 2 hours
4. Monitor fluorescence increase at Ex/Em = 540/590 nm

Important Thaw one vial of each kit component at room temperature before starting the experiment.

KEY PARAMETERS

Instrument:	Fluorescence microplate reader
Excitation:	540 nm
Emission:	590 nm
Cutoff:	570 nm
Recommended plate:	Solid black

PREPARATION OF STOCK SOLUTIONS

Unless otherwise noted, all unused stock solutions should be divided into single-use aliquots and stored at -20 °C after preparation. Avoid repeated freeze-thaw cycles.

1. NAD stock solution (100X):

Add 100 μL of H₂O into the vial of NAD (Component C) to make 100X NAD stock solution.

2. L-Lactate standard solution (100 mM):

Add 200 μL of H₂O or 1x PBS buffer into the vial of L-Lactate Standard (Component D) to make 100 mM L-Lactate standard solution.

PREPARATION OF STANDARD SOLUTION

L-Lactate standard

For convenience, use the Serial Dilution Planner:

<https://www.aatbio.com/tools/serial-dilution/13814>

Add 10 μL of L-Lactate stock solution into 990 μL PBS buffer to generate 1 mM L-Lactate standard solution (Lac7). Take the 1 mM L-Lactate standard solution and perform 1:3 serial dilutions to get serial dilutions of L-Lactate standard (Lac6 - Lac1).

Note Diluted L-Lactate standard solution is unstable and should be used within 4 hours.

PREPARATION OF WORKING SOLUTION

Add 5 mL of Assay Buffer (Component B) into one bottle of Enzyme Probe (Component A). Add 50 μL NAD stock solution (100X) into the bottle of Component A, and mix well.

Note This L-lactate working solution is not stable, and should be used promptly.

SAMPLE EXPERIMENTAL PROTOCOL

Table 1. Layout of L-Lactate standards and test samples in a solid black 96-well microplate. Lac= L-Lactate Standards (Lac1 - Lac7, 1 μM to 1 mM), BL=Blank Control, TS=Test Samples.

BL	BL	TS	TS
Lac1	Lac1
Lac2	Lac2
Lac3	Lac3		
Lac4	Lac4		
Lac5	Lac5		
Lac6	Lac6		
Lac7	Lac7		

Table 2. Reagent composition for each well.

Well	Volume	Reagent
Lac1 - Lac7	50 μL	Serial Dilutions (1 μM to 1 mM)
BL	50 μL	Dilution Buffer
TS	50 μL	Test Sample

1. Prepare L-Lactate standards (Lac), blank controls (BL), and test samples (TS)

according to the layout provided in Tables 1 and 2. For a 384-well plate, use 25 μL of reagent per well instead of 50 μL .

2. Add 50 μL of L-Lactate working solution to each well of L-Lactate standard, blank control, and test samples to make the total L-Lactate assay volume of 100 μL /well. For a 384-well plate, add 25 μL of L-Lactate working solution into each well instead, for a total volume of 50 μL /well.
3. Incubate the reaction at room temperature for 30 minutes to 2 hours, protected from light.
4. Monitor the fluorescence increase with a fluorescence plate reader at Excitation = 530 - 570 nm, Emission = 590 - 600 nm (optimal Ex/Em = 540/590 nm, cut off at 570 nm).

EXAMPLE DATA ANALYSIS AND FIGURES

The reading (RFU) obtained from the blank standard well is used as a negative control. Subtract this value from the other standards' readings to obtain the baseline corrected values. Then, plot the standards' readings to obtain a standard curve and equation. This equation can be used to calculate L-Lactate samples. We recommend using the Online Linear Regression Calculator which can be found at:

<https://www.aatbio.com/tools/linear-logarithmic-semi-log-regression-online-calculator/>

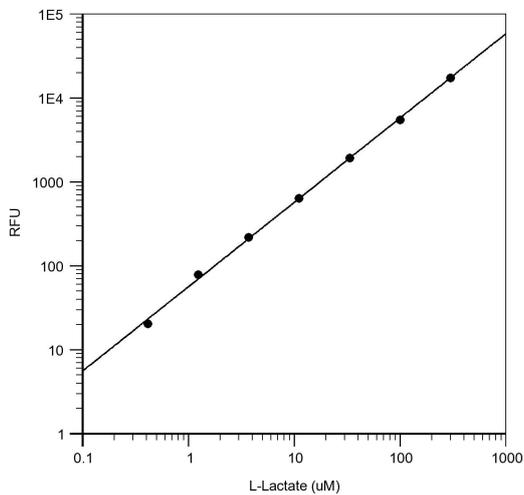


Figure 1. L-lactate dose response was measured with Amplite™ Fluorimetric L-Lactate Assay Kit in a 96-well solid black plate using a Gemini (Molecular Devices) microplate reader.

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.