

## PractiChrom™ L-Lactate Assay Kit (PLLC-25)

### Quantitative L-Lactate Determination Using PICOEXPLORER™

#### DESCRIPTION

L-LACTATE, or L-lactic acid, is generated by lactate dehydrogenase (LDH) under hypoxic or anaerobic conditions. L-lactate is added to many foods and beverages to give it a tart flavor. Increased levels of L-lactate in milk, egg, and fruit juices can be an indication of spoilage. In the wine industry, increasing levels of L-lactate and the decreasing levels of L-malic acid are monitored (Malolactic fermentation). In this process, the overall acidity of the wine is reduced and can lead to the improvement of the flavor of the wine.

BioAssay Systems' L-lactate assay is based on L-lactate dehydrogenase catalyzed oxidation of L-lactate in which the formed NADH reduces a chromogenic reagent. The intensity of product color is directly proportional to the L-lactate concentration in the sample.

#### KEY FEATURES

**Sensitive and accurate.** Detection limit of 0.05 mM (4.5 ppm, 0.45 mg/dL) and linearity to 2 mM (178 ppm, 17.8 mg/dL) L-lactate.

**Convenient.** Assay performed with portable PICO Explorer device.

**Cost efficient.** No need for expensive plate readers.

#### APPLICATIONS

**Direct Assays:** L-lactate in beverage samples (e.g. wine, beer, fruit juices, milk, etc) and biological samples (e.g. serum, plasma, etc).

#### KIT CONTENTS (25 TESTS)

<b>Assay Buffer:</b>	5 mL	<b>Standard:</b>	1.0 mL 20 mM L-Lactate
<b>Enzyme A:</b>	30 µL	<b>NAD/MTT:</b>	1.0 mL
<b>Enzyme B:</b>	120 µL	<b>ALT Enzyme:</b>	30 µL

**Storage conditions.** The kit is shipped on ice. Store all components at -20°C upon receiving. Shelf life: 6 months after receipt.

**Precautions:** reagents are for research use only. Equilibrate all components to room temperature prior assay. Normal precautions for laboratory reagents should be exercised while using the reagents. Please refer to Material Safety Data Sheet for detailed information.

#### PROCEDURES

##### Sample Preparation:

**White Wine, Serum, & Plasma.** Dilute 10× in dH<sub>2</sub>O by adding 10 µL sample to 90 µL dH<sub>2</sub>O.

**Beer & Juice.** Dilute 5× in dH<sub>2</sub>O by adding 20 µL sample to 80 µL dH<sub>2</sub>O.

**Red Wine.** Dilute 20× in dH<sub>2</sub>O by adding 10 µL sample to 190 µL dH<sub>2</sub>O.

**Milk samples** should be cleared by mixing 600 µL milk with 100 µL 6 M HCl. Centrifuge 5 min at 14,000 rpm. Transfer 300 µL supernatant into a clean tube and neutralize with 50 µL 6 M NaOH. The neutralized supernatant is ready for assay (dilution factor n = 1.36).

Other colored samples or samples with high levels of L-lactate will require a dilution. For preparation protocols for other samples, please contact our technical support at [info@bioassaysys.com](mailto:info@bioassaysys.com)

##### Procedure

1. Prepare 2 mM L-lactate Standard by mixing 10 µL of the provided 20 mM Standard and 90 µL dH<sub>2</sub>O in an Eppendorf tube.
2. In separate PCR tubes, add 10 µL dH<sub>2</sub>O and 10 µL 2 mM L-lactate Standard.

**Samples.** Add 10 µL sample to one PCR tube. For samples that are not colorless (e.g. beer, red wine, most fruit juices, etc), add 10 µL sample to two separate tubes, one serving as the Sample tube and one as the Sample Blank tube.

**Reagent Preparation.** Prepare sufficient Working Reagent for all dH<sub>2</sub>O, Standard, and Sample tubes by mixing, for each tube: 40 µL Assay Buffer, 4 µL NAD/MTT, 0.5 µL Enzyme A, 0.5 µL Enzyme B, and 0.5 µL ALT Enzyme. For each Sample Blank tube, prepare Blank Working Reagent (without Enzyme A): 40 µL Assay Buffer, 4 µL NAD/MTT, 0.5 µL Enzyme B, and 0.5 µL ALT Enzyme.

Quickly add 40 µL Working Reagent to all dH<sub>2</sub>O, Standard, and Sample tubes. To each Sample Blank tube, add 40 µL Blank Working Reagent (No Enzyme A). Close the tubes, briefly vortex or tap to mix. Tap tube on bench to settle liquid to the bottom of the tube if needed. Incubate for 10 min at room temperature in the dark.

3. Please refer to the PICOEXPLORER™ User's Manual for detailed instructions for operating the device.

Download the PAS-110 application. Turn on Bluetooth.

Push the Power button on the device. Then, open the app and tap the Connection Setting button and connect the device.

##### Measuring a Standard Curve (See pg 17-19 in User's Manual)

Return to the main menu and tap the Standard Curve button. Set the following:

LED Output: 10%

Unit: mM

RBG Selection: G

Tap the first Known Concentration Data Input Area box and input 0.0. Then, tap on the second box and input 2.0 (this represents the 0 and 2 mM L-lactate Standards). Then, place the dH<sub>2</sub>O tube into the measurement chamber of the photo absorbance sensor. Tap the Known Concentration Measurement Input Area (the box below 0.0), and click Measure. Remove the tube, then place the 2 mM L-lactate Standard into the measurement chamber. Tap the box below 2.0 and click Measure. Click Graph to view the standard curve.

##### Measuring Sample Concentrations

Return to the main menu and tap the Measure button. Edit the LED output, Units, and RBG selection as done above for the standard curve.

Place each Sample and Sample Blank tube into the measurement chamber of the photo absorbance sensor and tap Measure.

#### CALCULATION

The "concentration" will be displayed on the PICOEXPLORER™ for each Sample and Sample Blank. To calculate the L-lactate concentration in the sample, subtract the Sample Blank concentration from the Sample concentration and multiply by the dilution factor used (e.g. 5, 10, 20, etc). If no sample blank was used, simply multiply the Sample concentration by the dilution factor.

$$[\text{L-Lactate}] = ([\text{Sample}] - [\text{Sample Blank}]) \times n \text{ (mM)}$$

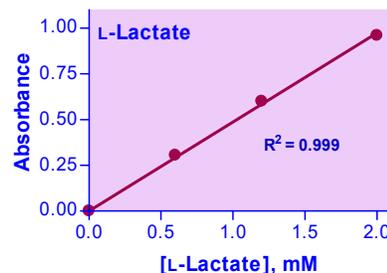
where [Sample] is the concentration of sample plus Enzyme A, [Sample Blank] is the concentration of sample *without* Enzyme A and *n* is the dilution factor.

*Note: if the sample concentration says "Out of range" the sample is not within the linear range of the assay. If the color of the tube is yellow like the dH<sub>2</sub>O tube, then the sample has low levels of L-lactate that cannot be detected by the assay. If the sample is very dark, dilute further in water and repeat the assay. Multiply the results by the dilution factor.*

**Conversions:** 1 mM L-lactate equals 8.9 mg/dL, or 89 ppm.

#### MATERIALS REQUIRED, BUT NOT PROVIDED

Pipetting devices, PCR tubes (e.g. Watson 137-211c 0.2 mL; or Cat# PCR-50 from BioAssay Systems), Eppendorf tubes (e.g. Phenix Cat# MAX-715, or Cat # EPP-50 from BioAssay Systems), and PICOEXPLORER™ (Cat # PICO001).



Standard Curve in water measured with PICOEXPLORER™