

# L-lactate Colorimetric Assay Kit (Lactate Oxidase Method)

ISO 13485:2016

Catalog #: MA-LAC

## Introduction

Lactate is an intermediary product of carbohydrate metabolism and exists in the human body as two optical isomers: L-lactate and D-lactate. L-lactate is produced from pyruvate during anaerobic glycolysis and is present in blood at concentrations 100 times greater than D-lactate. Its concentration in the blood is dependent on the liver's metabolic rate, and its production by the erythrocytes and muscle cells. L-lactate has often been used as a physiological indicator for stress. Its determination in serum is very important in the diagnosis and medical management of various diseases, such as tissue hypoxia, diabetes, circulatory failure, and hematological disorders.

L-lactate Colorimetric Assay Kit provides a simple, reproducible, and sensitive tool of measuring L-lactate concentration in plasma, serum, cell lysates, urine and other biological liquid samples. This assay employs a coupled enzymatic reaction system that combines multiple steps. Initially, lactate oxidase catalyzes the oxidation of L-lactate to pyruvate and hydrogen peroxide (H<sub>2</sub>O<sub>2</sub>). Peroxidase then catalyzes the reaction of hydrogen peroxide with the Colorimetric Probe to form a pink colored product. The optical density measured at 510nm is directly proportional to the concentration of L-lactate present in the sample.

## Storage

The entire kit should be stored at < -20 °C for up to 6 months from the date of shipment. For prepared reagent storage, see table below.

Component	Size / Description	Storage After Preparation
Microplate (Item A)	A 96-well (12 strips x 8 wells) plate	RT*
Assay Dilution Buffer	10 ml	2–8 °C, 6 months
L-lactate Standard	1 vial (100 µl of 4.9mM)	2–8 °C, 1 week
Enzyme Mix	2 amber vials (Lyophilized)	Do not store and reuse
Enzyme Buffer	25 ml	2–8 °C, 6 months
Colorimetric Probe	2 amber vials (Lyophilized)	Do not store and reuse

RT = room temperature

\*Return unused wells to the pouch containing desiccant pack, reseal along entire edge.

## Additional Materials Required

1. Microplate reader capable of measuring absorbance at 510 nm
2. Precision pipettes to deliver 2 µl to 1 ml volumes
3. Tubes to prepare sample dilutions
4. Incubator at 37°C
5. 50 ml conical tubes

## Sample Tips and General Considerations

NOTE: Optimal methods of sample preparation will need to be determined by each researcher empirically based on researched literature and knowledge of the samples.

- If not using fresh samples, freeze samples as soon as possible after collection.
- Avoid multiple freeze-thaw cycles. If possible, sub-aliquot samples prior to initial storage.
- It is strongly recommended to add a protease inhibitor cocktail to cell and tissue lysate samples.
- Avoid sonication of 1 ml or less as this can quickly heat and denature proteins.
- Most samples will not need to be concentrated. If concentration is required, a spin column concentrator with a chilled centrifuge is recommended.

### 1. Cell lysates can be prepared as follows:

For attached cells, remove supernatant from cell culture, wash cells twice with cold 1X PBS (for suspension cells, pellet the cells by spinning down the cells at 1,000 x g for 10 min) making sure to remove any remaining PBS before adding lysis buffer. Solubilize the cells at  $2 \times 10^7$  cells/ml in lysis buffer containing protease inhibitors. Pipette up and down to resuspend cells and rock the lysates gently at 2–8 °C for 30 minutes. Transfer extracts to microfuge tubes and centrifuge at 14,000 x g for 10 minutes.

It is recommended that sample protein concentrations should be determined using a total protein assay. Lysates should be used immediately or aliquot and stored at -70 °C. Thawed lysates should be kept on ice prior to use.

General tips for preparing lysate samples can be viewed on the online Resources page of the website:

<https://www.raybiotech.com/tips-on-sample-preparation/>

### 2. Plasma samples:

Collect blood with an anticoagulant such as citrate, EDTA or oxalate and mix by inversion. Centrifuge the blood at 1000 x g at 4°C for 10 minutes. Collect plasma supernatant without disturbing the white buffy layer. Sample should be tested immediately or frozen at -80°C for storage. Typically, normal human plasma has L-lactate concentrations in the range of 750-2000 µM. It is recommended to dilute the sample at least 1:2 with Assay Dilution Buffer before assaying.

### 3. Serum samples:

Collect blood in a tube with no anticoagulant. Allow the blood to clot at room temperature for 30 minutes. Centrifuge at 2500 x g for 20 minutes. Remove the yellow serum supernatant without disturbing the white buffy layer. Samples should be tested immediately or frozen at -80°C for storage. Typically, normal human serum has L-lactate concentrations in the range of 750-2000 µM. It is recommended to dilute the sample at least 1:2 with Assay Dilution Buffer before assaying.

### 4. Urine samples:

To remove insoluble particles, spin at 10,000 x g for 5 min. The supernatant can be assayed directly.

## Reagent Preparation

REAGENT	PREPARATION
Microplate (Item A)	No preparation
Assay Dilution Buffer	
Enzyme Buffer	
L-lactate Standard	See "Standard Preparation" section
Enzyme Mix	For each vial, reconstitute with 1 ml Enzyme Buffer.
Colorimetric Probe	For each vial, reconstitute with 1 ml Enzyme Buffer.
Working Solution	Add 1 mL reconstituted Enzyme Mix and 1 mL reconstituted Colorimetric Probe into 9 mL Enzyme Buffer to make the <b>Working Solution</b> . This is enough to assay 55 wells. Prepare the additional vial as needed. The Working Solution is stable for 1 day at 2-8°C in the dark.

## Standard Preparation

To prepare a dilution series of L-lactate Standard in the concentration range of 0  $\mu\text{M}$  – 2000  $\mu\text{M}$  (see Table below),

1. Label 8 microtubes #1 through 8 which with the following concentrations: 2000, 1000, 500, 250, 125, 62.5, 31.25, 0  $\mu\text{M}$ .
2. Pipette 59.5  $\mu\text{L}$  Assay Dilution Buffer into labeled tube #1, and 40  $\mu\text{L}$  Assay Dilution Buffer into labeled tube #2 – tube #8.
3. Pipette 40.5  $\mu\text{L}$  of L-lactate Standard into tube#1, mix thoroughly and quick spin, to make a 2000  $\mu\text{M}$  L-lactate Standard stock.
4. To make the 1000  $\mu\text{M}$  standard, pipette 40  $\mu\text{L}$  of tube #1 into the tube labeled #2. Mix thoroughly and quick spin.
5. Repeat this step with each successive concentration, preparing a dilution series as shown in the Table below. Each time, use 40  $\mu\text{L}$  of the prior concentration until the 31.25  $\mu\text{M}$  is reached. Mix each tube thoroughly before the next transfer. Standards should be prepared fresh and used immediately.

Labeled Tubes	L-lactate Standard ( $\mu\text{L}$ )	Assay Dilution Buffer ( $\mu\text{L}$ )	Standard Conc. ( $\mu\text{M}$ )	Standard Conc. (mg/dL)
1	40.5 $\mu\text{L}$ of L-lactate Standard	59.5 $\mu\text{L}$	2000	40.49
2	40 $\mu\text{L}$ of Tube #1	40 $\mu\text{L}$	1000	20.25
3	40 $\mu\text{L}$ of Tube #2	40 $\mu\text{L}$	500	10.12
4	40 $\mu\text{L}$ of Tube #3	40 $\mu\text{L}$	250	5.06
5	40 $\mu\text{L}$ of Tube #4	40 $\mu\text{L}$	125	2.53
6	40 $\mu\text{L}$ of Tube #5	40 $\mu\text{L}$	62.5	1.27
7	40 $\mu\text{L}$ of Tube #6	40 $\mu\text{L}$	31.25	0.63
8	0 $\mu\text{L}$	40 $\mu\text{L}$	0	0

## Assay Procedure

Each L-lactate standard and sample should be assayed in duplicate or triplicate. A freshly prepared standard curve should be used each time the assay is performed.

1. Add 5  $\mu\text{L}$  of the diluted L-lactate Standards or samples to the 96-well microtiter plate.
2. Initiate the reaction by adding 200  $\mu\text{L}$  Working Solution (See “Reagent Preparation”) to each well.
3. Cover with the plate cover. Carefully shake the plate for a few seconds to mix.
4. Incubate the plate for 5 minutes at 37°C.
5. Measure the absorbance at 510nm using a plate reader.

## Calculation of Results

### Subtract the blanks

Average the absorbance value of the blank wells (Standard 0  $\mu\text{M}$ ) and subtract this from the absorbance values of all the other wells. These are the corrected absorbance.

### Plotting the standard curves

Make a plot of corrected absorbance at 510nm as a function of L-lactate concentration.

### Determination of sample L-lactate concentration

$$\text{L-lactate } (\mu\text{M}) = \frac{OD_{\text{Sample}} - OD_{\text{Blank}}}{\text{Slope}} \times DF$$

$OD_{\text{Sample}}$  = Optical density (OD) reading of the Sample

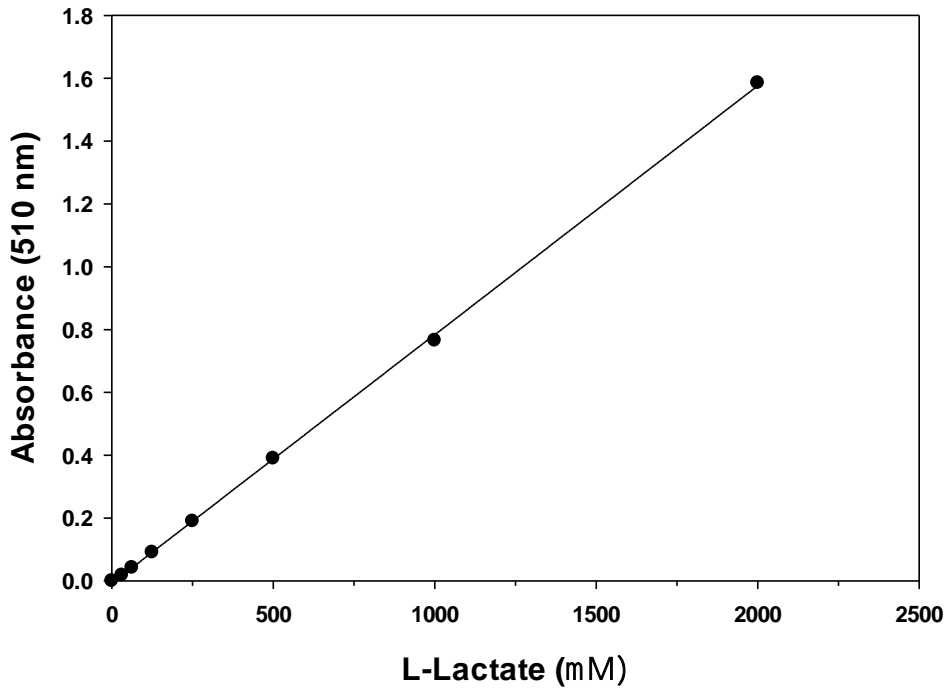
$OD_{\text{Blank}}$  = Optical density (OD) reading of the Blank (Standard 0  $\mu\text{M}$ )

Slope is from the plot of L-lactate concentration vs. Absorbance shown in Typical data below

DF = Sample Dilution factor (DF = 1 for undiluted Samples)

*Note:* If the calculated L-lactate concentration of the sample is higher than 2000  $\mu\text{M}$ , dilute the sample in assay diluent solution and repeat the assay.

## A. Typical Data



These standard curves are for demonstration only. A standard curve must be run with each assay.

## B. Reproducibility

**Intra-assay Precision (Precision within an assay):**

To assess intra-assay precision, 16 wells per sample (total of 4 samples) were tested on a single plate. The intra-assay coefficient of variation was found to be 1.4%.

**Inter-assay Precision (Precision between assays):**

To assess inter-assay precision, 4 samples were tested in separate assays (n=4). The inter-assay coefficient of variation was found to be 1.8%.

This product is for research use only.