

## Product Information

### L-Lactate Assay Kit

Catalog Number **MAK329**  
 Storage Temperature  $-20\text{ }^{\circ}\text{C}$

## TECHNICAL BULLETIN

### Product Description

Lactate is generated by lactate dehydrogenase (LDH) under hypoxic or anaerobic conditions. Monitoring lactate levels is, therefore, a good indicator of the balance between tissue oxygen demand and utilization, and is useful when studying cellular and animal physiology.

Simple, direct, and automation-ready procedures for measuring lactate concentration are very desirable. The Lactate Assay Kit is based on lactate dehydrogenase catalyzed oxidation of lactate, in which the formed NADH reduces a formazan (MTT) reagent. The intensity of the product color, measured at 565 nm, is proportionate to the lactate concentration in the sample.

Detection limit of 0.05 mM and linearity up to 2 mM L-Lactate in the 96 well plate assay. For cell culture samples containing phenol red: detection limit of 0.1 mM and linearity up to 1 mM L-Lactate using the 96 well plate assay.

This kit is suitable for lactate determination in serum, plasma, and cell media samples.

### Components

The kit is sufficient for 100 colorimetric assays in 96 well plates.

Assay Buffer Catalog Number MAK329A	10 mL
Enzyme A Catalog Number MAK329B	120 $\mu\text{L}$
Enzyme B Catalog Number MAK329C	120 $\mu\text{L}$
NAD Solution Catalog Number MAK329D	1 mL

MTT Solution Catalog Number MAK329E	1.5 mL
--	--------

Standard (20 mM L-Lactate) Catalog Number MAK329F	1 mL
--	------

### Reagents and Equipment Required but Not Provided.

- Pipetting devices and accessories (e.g., multichannel pipettor)
- Centrifuge tubes
- 96 well flat bottom plate. It is recommended to use clear plates for colorimetric assays
- Spectrophotometric multiwell plate reader

### Precautions and Disclaimer

This product is for R&D use only, not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

### Storage/Stability

The kit is shipped on dry ice. Store all components at  $-20\text{ }^{\circ}\text{C}$  upon receiving.

### Procedure

#### Sample Preparation

For samples with potential endogenous enzyme activity (i.e. serum, plasma, tissue extracts), two reactions should be run: one with added Enzyme A and a control reaction with No Enzyme A.

Serum and Plasma should be diluted at least 2 $\times$  with ultrapure water prior to the assay.

The following substances interfere with the assay and should be avoided in sample preparation: ascorbic acid, sodium dodecyl sulfate (>0.2%), sodium azide, NP-40 (>1%) and TWEEN® 20 (>1%).

### Standard Curve

Prepare 1,000  $\mu\text{L}$  of 2.0 mM L-Lactate Premix by mixing 100  $\mu\text{L}$  of 20 mM Standard and 900  $\mu\text{L}$  of ultrapure water. For cell culture samples containing phenol red, prepare 1,000  $\mu\text{L}$  of 1.0 mM lactate Premix by mixing 50  $\mu\text{L}$  of 20 mM Standard and 950  $\mu\text{L}$  of culture medium without serum. Dilute standard according to Table 1 (samples excluding cell culture containing phenol red) or Table 2 (cell culture samples containing phenol red).

**Table 1.**

Preparation of L-Lactate Standards for all samples excluding cell culture samples containing phenol red

Tube	Premix	Ultrapure Water	L-Lactate (mM)
1	100 $\mu\text{L}$	0 $\mu\text{L}$	2.0
2	80 $\mu\text{L}$	20 $\mu\text{L}$	1.6
3	60 $\mu\text{L}$	40 $\mu\text{L}$	1.2
4	40 $\mu\text{L}$	60 $\mu\text{L}$	0.8
5	30 $\mu\text{L}$	70 $\mu\text{L}$	0.6
6	20 $\mu\text{L}$	80 $\mu\text{L}$	0.4
7	10 $\mu\text{L}$	90 $\mu\text{L}$	0.2
8	0 $\mu\text{L}$	100 $\mu\text{L}$	0

**Table 2.**

Preparation of L-Lactate Standards for cell culture samples containing phenol red

Tube	Premix	Culture Medium	L-Lactate (mM)
1	100 $\mu\text{L}$	0 $\mu\text{L}$	1.0
2	80 $\mu\text{L}$	20 $\mu\text{L}$	0.8
3	60 $\mu\text{L}$	40 $\mu\text{L}$	0.6
4	40 $\mu\text{L}$	60 $\mu\text{L}$	0.4
5	30 $\mu\text{L}$	70 $\mu\text{L}$	0.3
6	20 $\mu\text{L}$	80 $\mu\text{L}$	0.2
7	10 $\mu\text{L}$	90 $\mu\text{L}$	0.1
8	0 $\mu\text{L}$	100 $\mu\text{L}$	0

### Reaction Mix

**Note:** Briefly centrifuge enzyme tubes before opening. For each Sample and Standard well, prepare Reaction Mix by mixing:

60  $\mu\text{L}$  of Assay Buffer  
 1  $\mu\text{L}$  of Enzyme A  
 1  $\mu\text{L}$  of Enzyme B  
 10  $\mu\text{L}$  of NAD  
 14  $\mu\text{L}$  of MTT

Fresh preparation just prior to use is recommended.

For the samples which are control reactions with No Enzyme A (see Sample Preparation, the Reaction Mix includes:

60  $\mu\text{L}$  of Assay Buffer  
 1  $\mu\text{L}$  of Enzyme B  
 10  $\mu\text{L}$  of NAD  
 14  $\mu\text{L}$  of MTT

### Assay Reaction

1. Transfer 20  $\mu\text{L}$  of standards into separate wells of a clear, flat bottom 96 well plate.
2. Transfer 20  $\mu\text{L}$  of each sample into separate wells.
3. Add 80  $\mu\text{L}$  of Reaction Mix per reaction well quickly. Tap plate to mix briefly and thoroughly. **Note:** This assay is based on an enzyme-catalyzed kinetic reaction. Addition of Reaction Mix should be quick and mixing should be brief but thorough. Use of a multichannel pipettor is recommended.
4. Immediately measure the initial absorbance at 565 nm ( $A_{565}$ ).
5. Incubate plate for 20 minutes at room temperature.
6. Measure the final absorbance at 565 nm ( $A_{565}$ ).

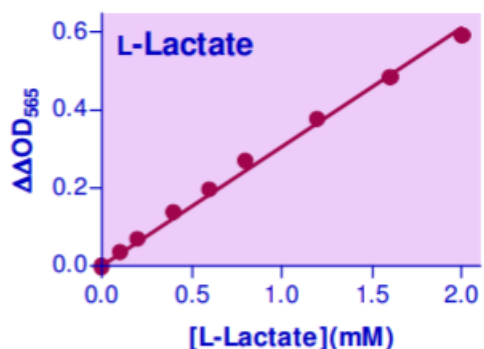
## Results

Subtract initial  $A_{565}$  from the final  $A_{565}$  for the standard and sample wells. Use the  $\Delta A_{565}$  values to determine the sample L-lactate concentration from the standard curve.

For samples requiring a No Enzyme A control reaction, subtract the  $\Delta A_{565}$  No Enzyme value from the  $\Delta A_{565}$  Sample and use this value to determine the sample L-lactate concentration from the standard curve.

**Note:** If the sample  $A_{565}$  value is higher than  $A_{565}$  for 2 mM L-lactate standard, dilute sample in ultrapure water and repeat the assay. Multiply the results by the dilution factor.

**Figure 1.**  
Typical Standard Curve



## References

1. Senadheera, D. et al., Inactivation of VicK affects acid production and acid survival of *Streptococcus mutans*. J. Bacteriol., **191(20)**, 6415-24 (2009).
2. Le Nihouannen, D. et al., Ascorbic acid accelerates osteoclast formation and death. Bone, **46(5)**, 1336-43 (2009).
3. Milovanova, T.N. et al., Lactate stimulates vasculogenic stem cells via the thioredoxin system and engages an autocrine activation loop involving hypoxia-inducible factor 1. Mol. Cell Biol., **28(20)**, 6248-61 (2008).

TWEEN is a registered trademark of Croda International PLC

HM,MAM 11/18-1