

## Product Information

### Pyruvate Assay Kit

Catalog Number **MAK332**  
Storage Temperature  $-20\text{ }^{\circ}\text{C}$

## TECHNICAL BULLETIN

### Product Description

Pyruvate is a key intermediate in cellular metabolic pathways. Pyruvate can be converted to carbohydrates via gluconeogenesis, to fatty acids or energy through acetyl-CoA, to the amino acid alanine, and to ethanol. Abnormal levels of pyruvate have been linked to liver diseases and metabolic disorders.

Simple, direct, and automation-ready procedures for measuring pyruvate concentrations find wide applications in research and drug discovery. The Pyruvate Assay Kit uses a single Reaction Mix that combines pyruvate oxidase and hydrogen peroxide determination in one step. The color intensity of the reaction product at 570 nm or fluorescence intensity at  $\lambda_{\text{ex}} = 530\text{ nm}/\lambda_{\text{em}} = 585\text{ nm}$  is directly proportional to the pyruvate concentration in the sample.

The Pyruvate Assay Kit has a linear detection range for pyruvate in a 96 well plate assay of 2–500  $\mu\text{M}$  (17  $\mu\text{g}/\text{dL}$  to 4.4  $\text{mg}/\text{dL}$ ) for colorimetric assays and 0.2–50  $\mu\text{M}$  for fluorometric assays.

Suitable for pyruvate detection in biological samples and for studying the effects of drugs on pyruvate metabolism.

### Components

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

Enzyme Mix Catalog Number MAK332A	10 mL
Dye Reagent Catalog Number MAK332B	120 $\mu\text{L}$
Standard (25 mM Pyruvate) Catalog Number MAK332C	400 $\mu\text{L}$

### Reagents and Equipment Required but Not Provided.

- Pipetting devices and accessories (e.g., multichannel pipettor)
- Centrifuge tubes
- 96 well flat bottom plates— It is recommended to use black plates with clear bottoms for fluorescence assays and clear plates for colorimetric assays.
- Fluorescence or 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.

### Procedures

Equilibrate all components to room temperature prior to assay.

#### Sample Preparation

SH-group containing reagents (e.g. mercaptoethanol, DTT) may interfere with this assay and should be avoided in sample preparation.

#### Standard Preparation for Colorimetric Procedure

Prepare a 500  $\mu\text{M}$  Standard Premix by mixing 10  $\mu\text{L}$  of the 25 mM Standard and 490  $\mu\text{L}$  of ultrapure water. Dilute 500  $\mu\text{M}$  Standard Premix according to Table 1.

**Table 1.**

Preparation of Pyruvate Standards

Tube	Standard Premix	Ultrapure Water	Pyruvate ( $\mu\text{M}$ )
1	100 $\mu\text{L}$	0 $\mu\text{L}$	500
2	80 $\mu\text{L}$	20 $\mu\text{L}$	400
3	60 $\mu\text{L}$	40 $\mu\text{L}$	300
4	40 $\mu\text{L}$	60 $\mu\text{L}$	200
5	30 $\mu\text{L}$	70 $\mu\text{L}$	150
6	20 $\mu\text{L}$	80 $\mu\text{L}$	100
7	10 $\mu\text{L}$	90 $\mu\text{L}$	50
8	0 $\mu\text{L}$	100 $\mu\text{L}$	0

#### Reaction Mix

For each well of reaction, prepare Reaction Mix by mixing into a clean tube:

94  $\mu\text{L}$  of Enzyme Mix  
1  $\mu\text{L}$  of Dye Reagent

#### Colorimetric Assay Reaction

1. Transfer 10  $\mu\text{L}$  of standards and 10  $\mu\text{L}$  of samples into separate wells of a clear flat bottom 96 well plate.
2. Add 90  $\mu\text{L}$  of the Reaction Mix to each standard and sample well. Tap plate to mix well. Freeze unused reagents for future use.
3. Incubate 30 minutes at room temperature.
4. Measure the absorbance at 570 nm ( $A_{570}$ ).  
Note: If the Sample ( $A_{570}$ ) is higher than the 500  $\mu\text{M}$  Standard ( $A_{570}$ ), dilute sample in ultrapure water and repeat the assay. Multiply result by the dilution factor.

#### Fluorometric Procedure

For the fluorometric assay, the linear detection range is 0.2–50  $\mu\text{M}$  pyruvate.

1. Dilute the Standards prepared in Colorimetric Procedure 10-fold in ultrapure water.
2. Transfer 10  $\mu\text{L}$  of standards and 10  $\mu\text{L}$  of samples into separate wells of a black 96 well plate.
3. Add 90  $\mu\text{L}$  of the Reaction Mix to each standard and sample well. Tap plate to mix well. Freeze unused reagents for future use.
4. Tap plate to mix. Incubate 30 minutes at room temperature.
5. Read fluorescence at  $\lambda_{\text{ex}} = 530 \text{ nm} / \lambda_{\text{em}} = 585 \text{ nm}$ .  
Note: If assays in 384 well plate are desired, use 5  $\mu\text{L}$  of Standards or 5  $\mu\text{L}$  of samples, and 45  $\mu\text{L}$  of the Reaction Mix to each standard or sample well.

## Results

Subtract the blank  $A_{570}$  (Tube 8) from the standard  $A_{570}$  values and plot the  $A_{570}$  against standard concentrations. Determine the slope using linear regression fitting. Calculation of the pyruvate concentration for the Sample in the colorimetric assay:

$$[\text{Pyruvate } \mu\text{M}] = \frac{A_{570 \text{ Sample}} - A_{570 \text{ Blank}}}{\text{Slope}}$$

Calculation of the pyruvate concentration for the Sample in the fluorometric assay:

$$[\text{Pyruvate } \mu\text{M}] = \frac{F_{\text{Sample}} - F_{\text{Blank}}}{\text{Slope}}$$

## References

1. Hansen, J.L., and Freier, E.F., Direct assays of lactate, pyruvate,  $\beta$ -hydroxybutyrate, and acetoacetate with a centrifugal analyzer. *Clin. Chem.*, **24(3)**, 475-9 (1978).
2. Sutherland, D.V. et al., *Trypanosoma evansi*: measurement of pyruvate production as an indicator of the drug sensitivity of isolates *in vitro*. *Trop. Med. Parasitol.*, **46(2)**, 93-8 (1995).
3. Chariot, P. et al., Optimal handling of blood samples for routine measurement of lactate and pyruvate. *Arch. Pathol. Lab. Med.*, 118(7), 695-7 (1994).

HM,MAM 11/18-1

## Figure 1.

Serum samples were run in duplicate according to the standard procedure.

