

## Product Information

### Citrate Assay Kit

Catalog Number **MAK333**  
Storage Temperature  $-20\text{ }^{\circ}\text{C}$

## TECHNICAL BULLETIN

### Product Description

Citrate is an intermediate in the citric acid cycle and is involved in fatty acid synthesis. The Citrate Assay Kit provides a simple and automation-ready procedure for measuring citrate concentration. Citrate is converted into pyruvate which is then oxidized with the conversion of the dye into a colored, fluorescent form. The color intensity at 570 nm or fluorescence intensity at  $\lambda_{\text{ex}} = 530\text{ nm}/\lambda_{\text{em}} = 585\text{ nm}$  is directly proportional to the citrate concentration in the sample.

The Citrate Assay Kit has a linear detection range of 4–400  $\mu\text{M}$  citrate for the colorimetric assay and 0.5–40  $\mu\text{M}$  for the fluorimetric assay.

Suitable for citrate determination in biological samples (e.g. plasma, serum, urine, tissue and culture media.)

### Components

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

Developer Catalog Number MAK333A	10 mL
CL Enzyme (Dried) Catalog Number MAK333B	1 Vial
Dye Reagent Catalog Number MAK333C	120 $\mu\text{L}$
ODC Enzyme Catalog Number MAK333D	120 $\mu\text{L}$
Citrate Standard (10 mM) Catalog Number MAK333E	500 $\mu\text{L}$

### 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 black plates with clear bottoms for fluorescence assays and clear plates for colorimetric assays.
- Fluorescence or spectrophotometric multiwell plate reader
- Reagents and equipment for sample preparation

### Precautions and Disclaimer

This product is for Research Use Only. Not for Use in Diagnostic Procedures. 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.

### Preparation Instructions

#### Reagent Preparation

Dissolve the CL Enzyme with 120  $\mu\text{L}$  of Developer. Pipette up and down to assure the enzyme is fully dissolved. Reconstituted CL Enzyme Solution is stable for 4 weeks stored at  $-20\text{ }^{\circ}\text{C}$ . Before each use of the CL Enzyme Solution, pipette up and down to assure the enzyme is resuspended.

### Sample Preparation

Tissue or cell samples ( $2 \times 10^6$ ) can be homogenized in 100  $\mu\text{L}$  of phosphate buffered saline (PBS). Centrifuge at 14,000 rpm for 5 minutes. Use clear supernatant for assay. If planning to measure citrate in culture media, avoid media with high pyruvate concentrations (e.g. DMEM, L-15, F12, etc.).

Serum and plasma samples should be deproteinated using a 10 kDa spin filter (e.g. Amicon<sup>®</sup> Ultra-0.5 filter). Alternatively, untreated serum and plasma can be measured directly if an internal standard is used.

Urine samples should be diluted at least 5-fold and an internal standard should be used.

### **Procedures**

#### Colorimetric Procedure

For the colorimetric assay, the linear detection range is 4–400  $\mu\text{M}$  citrate.

#### Standard Preparation

Prepare the 400  $\mu\text{M}$  Citrate Standard Premix by mixing 10  $\mu\text{L}$  of 10 mM Citrate Standard with 240  $\mu\text{L}$  of ultrapure water. Next, dilute standards in 1.5-mL centrifuge tubes according to Table 1.

Note: If assaying culture medium with phenol red, dilute the Citrate Standard in culture medium.

**Table 1.**

Preparation of Citrate Standards

Tube	Standard Premix	Ultrapure Water	Citrate ( $\mu\text{M}$ )
1	100 $\mu\text{L}$	0 $\mu\text{L}$	400
2	60 $\mu\text{L}$	40 $\mu\text{L}$	240
3	30 $\mu\text{L}$	70 $\mu\text{L}$	120
4	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:

- 85  $\mu\text{L}$  of Developer
- 1  $\mu\text{L}$  of CL Enzyme Solution
- 1  $\mu\text{L}$  of ODC Enzyme
- 1  $\mu\text{L}$  of Dye Reagent

#### Sample Blank Reaction Mix

For each sample blank, prepare Sample Blank Reaction Mix by mixing into a clean tube:

- 85  $\mu\text{L}$  of Developer
- 1  $\mu\text{L}$  of ODC Enzyme
- 1  $\mu\text{L}$  of Dye Reagent

### 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 of a clear flat bottom 96 well plate. Each sample being tested requires a sample blank.

Note: If using an internal standard, samples will need three separate reactions:

- Sample plus Internal Standard
- Sample (alone)
- Sample Blank

For the internal standard prepare 500  $\mu\text{L}$  of 1,000  $\mu\text{M}$  Citrate Standard by mixing 50  $\mu\text{L}$  of 10 mM Citrate Standard and 450  $\mu\text{L}$  of ultrapure water.

For the Sample plus Internal Standard well, add 5  $\mu\text{L}$  of 1,000  $\mu\text{M}$  citrate and 20  $\mu\text{L}$  of sample.

For the Sample and Sample Blank wells, add 5  $\mu\text{L}$  of ultrapure water and 20  $\mu\text{L}$  of sample.

3. Add 80  $\mu\text{L}$  of the appropriate Reaction Mix to each Standard, Sample, and Sample Blank well. Mix well.
4. Incubate protected from light 15 minutes at room temperature.
5. Measure the absorbance at 570 nm ( $A_{570}$ ).

#### Fluorometric Procedure

For the fluorometric assay, the linear detection range is 1–40  $\mu\text{M}$  citrate.

1. Dilute the Standards prepared in Colorimetric Procedure 10-fold in ultrapure water. If an internal standard is used, use 5  $\mu\text{L}$  of 100  $\mu\text{M}$  citrate.
2. Transfer 20  $\mu\text{L}$  of standards and 20  $\mu\text{L}$  of samples (prepare 2 wells per sample if a standard curve is used; prepare 3 wells per sample if an internal standard is used, see Colorimetric Procedure) into separate wells of a black 96 well plate.
3. Add 80  $\mu\text{L}$  of appropriate Reaction Mix (see Colorimetric Procedure) to each well. Tap plate to mix.
4. Incubate protected from light for 15 minutes at room temperature.
5. Measure the fluorescence at  $\lambda_{\text{ex}} = 530 \text{ nm}$ /  
 $\lambda_{\text{em}} = 585 \text{ nm}$ .

## Results

Subtract the blank value (Standard Tube 4) from the standard values and plot the  $\Delta A_{570}$  or  $\Delta F$  against standard concentrations. Determine the slope using linear regression fitting. The citrate concentration of Sample is calculated as:

$$[\text{Citrate } (\mu\text{M})] = \frac{R_{\text{Sample}} - R_{\text{Blank}}}{\text{Slope } (\mu\text{M}^{-1})} \times n$$

If an internal standard was used, the sample citrate concentration is calculated as follows:

$$[\text{Citrate } (\mu\text{M})] = \frac{R_{\text{Sample}} - R_{\text{Blank}}}{R_{\text{Standard}} - R_{\text{Sample}}} \times \frac{[\text{Standard}] \times n}{4}$$

where:

$R_{\text{Sample}}$ ,  $R_{\text{Blank}}$ , and  $R_{\text{Standard}} = A_{570}$  or fluorescence readings of the Sample, Sample Blank, and the Sample plus Internal Standard, respectively.  
 $n$  = the sample dilution factor.

**Note:** The volume of the internal standard is 4× lower than the sample volume; thus, the internal standard concentration should be divided by 4.

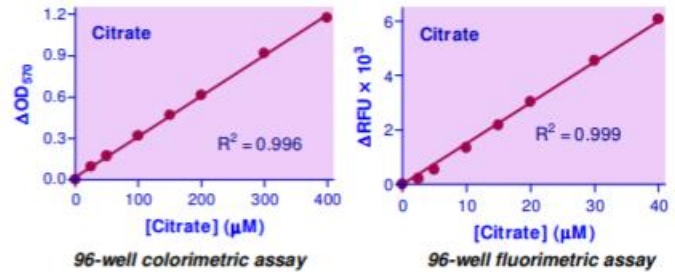
If the calculated citrate concentration is >400  $\mu\text{M}$  for the colorimetric assay or >40  $\mu\text{M}$  for the fluorimetric assay, dilute sample in ultrapure water and repeat assay. Multiply result by the dilution factor  $n$ .

## Conversions

100  $\mu\text{M}$  citrate = 19.1 mg/L, 0.0019%, or 19.1 ppm.

**Figure 1.**

Typical Standard Curves



## References

1. Tompkins, D., and Toffaletti, J., Enzymatic Determination of Citrate in Serum and Urine, with Use of the Worthington "Ultrafree" Device. *Clin. Chem.*, **28**, 192-5 (1982).
2. Costello, L.C., and O'Neill, J.J., A simplified and sensitive method for citrate determination in biological samples. *J. Appl. Physiol.*, **27**, 120-2 (1969).
3. Parvin, R. et al., Convenient rapid determination of picomole amounts of oxaloacetate and citrate. *Anal. Biochem.*, **104**, 296-9 (1980).

Amicon is a registered trademark of Merck KGaA, Darmstadt, Germany.

HM,VNC,MAM 11/18-1