

## EnzyChrom™ Citrate Assay Kit (ECIT-100)

### Quantitative Colorimetric/Fluorimetric Citrate Determination

#### DESCRIPTION

**CITRATE** is an intermediate in the citric acid cycle and is involved in fatty acid synthesis. BioAssay Systems' 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 and fluorescent form. The color intensity at 570 nm or fluorescence intensity at  $\lambda_{ex/em} = 530/585$  nm is directly proportional to the citrate concentration in the sample.

#### KEY FEATURES

**Fast and sensitive.** Linear detection range: 4 to 400  $\mu\text{M}$  citrate for colorimetric assays and 0.5 to 40  $\mu\text{M}$  for fluorimetric assays.

**Convenient and high-throughput.** Homogeneous "mix-incubate-measure" type assay. Can be readily automated on HTS liquid handling systems for processing thousands of samples per day.

#### APPLICATIONS

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

#### KIT CONTENTS (100 TESTS IN 96-WELL PLATES)

**Developer:** 10 mL

**CL Enzyme:** Dried

**Dye Reagent:** 120  $\mu\text{L}$

**ODC Enzyme:** 120  $\mu\text{L}$

**Citrate Standard:** 500  $\mu\text{L}$

**Storage conditions.** The kit is shipped on ice. Store all kit components at -20 °C. Shelf life of six months after receipt.

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

#### PROCEDURES

##### Reagent Preparation

Dissolve the CL Enzyme with 120  $\mu\text{L}$  Developer. Pipette up and down to assure the enzyme is fully dissolved. Reconstituted CL enzyme is stable for 4 weeks stored at -20°C. Before each use of the CL Enzyme, 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  $\mu\text{L}$  PBS. Centrifuge at 14,000 rpm for 5 min. 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 Ultra-0.5). 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.

##### Colorimetric Procedure

**1. Standards.** Dilute the Citrate Standard to 400  $\mu\text{M}$  by mixing 10  $\mu\text{L}$  10 mM Standard with 240  $\mu\text{L}$  dH<sub>2</sub>O. Next, dilute standards in 1.5-mL centrifuge tubes as described in the table. *If assaying culture media with phenol red, dilute the Citrate Standard in culture media.*

No	Premix + dH <sub>2</sub> O	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

Transfer 20  $\mu\text{L}$  of each standard to separate wells in a 96 well plate.

**2. Samples.** Add 20  $\mu\text{L}$  of each sample to two separate wells in a 96 well plate (each sample requires a sample blank).

If using an internal standard, samples will need three separate reactions: 1) sample plus standard, 2) sample alone and 3) sample blank. For the internal standard prepare 500  $\mu\text{L}$  1000  $\mu\text{M}$  citrate standard by mixing 50  $\mu\text{L}$  10 mM Standard and 450  $\mu\text{L}$  dH<sub>2</sub>O. For the sample plus standard well, add 5  $\mu\text{L}$  1000  $\mu\text{M}$  citrate and 20  $\mu\text{L}$  sample. For the sample and sample blank wells, add 5  $\mu\text{L}$  dH<sub>2</sub>O and 20  $\mu\text{L}$  sample.

**3. Citrate Detection.** Prepare enough working reagent (WR) for all standards and samples. For each reaction combine the following: 85  $\mu\text{L}$  Developer, 1  $\mu\text{L}$  CL Enzyme, 1  $\mu\text{L}$  ODC Enzyme, and 1  $\mu\text{L}$  Dye Reagent. For the Sample Blanks, prepare a WR **without** the CL Enzyme. Add 80  $\mu\text{L}$  of the appropriate WR to each Standard and Sample well. Mix well and incubate protected from light for 15 min at RT.

**4. Read OD<sub>570nm</sub>.**

##### Fluorimetric Procedure

For fluorimetric assays, the linear detection range is 1 to 40  $\mu\text{M}$  citrate. Dilute the standards prepared in *Colorimetric Procedure* 1:10 in dH<sub>2</sub>O. If an internal standard is used, use 5  $\mu\text{L}$  of 100  $\mu\text{M}$  citrate.

Transfer 20  $\mu\text{L}$  standards and 20  $\mu\text{L}$  samples (2 wells per sample if a standard curve is used; 3 wells per sample if an internal standard is used, see *Colorimetric Procedure*) into separate wells of a **black** 96-well plate. Add 80  $\mu\text{L}$  of appropriate Working Reagent (see *Colorimetric Procedure*) to each well. Tap plate to mix.

Incubate protected from light for 15 min at RT and read fluorescence at  $\lambda_{ex/em} = 530/585$  nm.

#### CALCULATION

Subtract the blank value (#4) from the standard values and plot the  $\Delta\text{OD}$  or  $\Delta\text{F}$  against standard concentrations. Determine the slope and calculate the citrate concentration of the Samples as follows:

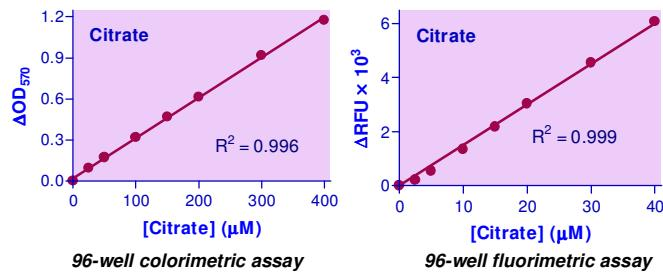
$$[\text{Citrate}] = \frac{R_{\text{SAMPLE}} - R_{\text{BLANK}}}{\text{Slope } (\mu\text{M}^{-1})} \times n \quad (\mu\text{M})$$

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

$$[\text{Citrate}] = \frac{R_{\text{SAMPLE}} - R_{\text{BLANK}}}{R_{\text{STANDARD}} - R_{\text{SAMPLE}}} \times \frac{[\text{Standard}]}{4} \times n \quad (\mu\text{M})$$

where  $R_{\text{SAMPLE}}$ ,  $R_{\text{BLANK}}$  and  $R_{\text{STANDARD}}$  are OD or fluorescence readings of the Sample, Sample Blank and the Sample plus Standard respectively.  $n$  is the sample dilution factor. Notes: The volume of the internal standard is 4x 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 dH<sub>2</sub>O and repeat assay. Multiply result by the dilution factor  $n$ .

**Conversions:** 100  $\mu\text{M}$  citrate equals 19.1 mg/L, 0.0019% or 19.1 ppm.



#### MATERIALS REQUIRED, BUT NOT PROVIDED

Pipetting devices, clear or black flat-bottom 96-well plates, plate reader or centrifuge tubes.

#### PUBLICATIONS

- Pant, A. et al (2021). Viral growth factor-and STAT3 signaling-dependent elevation of the TCA cycle intermediate levels during vaccinia virus infection. PLoS Pathogens, 17(2).
- Fu, X., et al (2018). Runx2/Osterix and zinc uptake synergize to orchestrate osteogenic differentiation and citrate containing bone apatite formation. Advanced Science 5:4: 1700755.
- Trivedi, A. K., et al. (2015). Adaptation of oxidative phosphorylation to photoperiod-induced seasonal metabolic states in migratory songbirds. Comparative Biochemistry and Physiology Part A: Molecular & Integrative Physiology 184): 34-40.

