

## EnzyChrom™ L-Carnitine Assay Kit (ELCR-100)

### Quantitative Colorimetric/Fluorimetric L-Carnitine Determination

#### DESCRIPTION

L-CARNITINE is found in most mammals, plants, and some bacteria. L-Carnitine's most crucial role is in the transport of long-chain fatty acids from the cytosol into the mitochondria for fatty acid oxidation. BioAssay Systems' method provides a simple and high-throughput assay for measuring L-Carnitine. In this assay, a multistep reaction produces H<sub>2</sub>O<sub>2</sub> which reacts with a specific dye to form a pink colored product. The optical density at 570nm or fluorescence intensity at  $\lambda_{ex/em} = 530/585$  nm is directly proportional to the L-Carnitine concentration in the sample.

#### KEY FEATURES

**Sensitive.** Uses 10  $\mu$ L samples. Detection range: colorimetric assay 12 - 1000  $\mu$ M, fluorimetric assay 1 - 100  $\mu$ M L-Carnitine.

**Convenient.** Room temperature "mix-incubate-read" procedure can be readily automated for high-throughput assay of thousands of samples per day.

#### APPLICATIONS

**Assays:** L-Carnitine in urine, serum, plasma, cell lysates, tissue lysates and food products (e.g. milk).

#### KIT CONTENTS

<b>Assay Buffer:</b>	10 mL	<b>Dye Reagent:</b>	120 $\mu$ L
<b>Enzyme A:</b>	Dried	<b>Enzyme B:</b>	100 $\mu$ L
<b>Enzyme C:</b>	100 $\mu$ L	<b>Standard:</b>	50 $\mu$ L
<b>Substrate Mix:</b>	400 $\mu$ L		

**Storage conditions.** The kit is shipped on ice. Store all 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 Safety Data Sheet for detailed information.

#### PROCEDURES

##### Reagent Preparation:

Reconstitute Enzyme A by adding 120  $\mu$ L dH<sub>2</sub>O to the Enzyme A tube. Make sure Enzyme A is fully dissolved by pipetting up and down and incubating at RT for 15 min. Store reconstituted Enzyme A at -20°C and use within 2 months.

##### Sample Preparation:

Serum, plasma, and milk should be clarified by perchloric acid:potassium hydroxide precipitation (see below). Serum and Plasma should be diluted no more than 1:5 with an internal standard. Cell lysates may be assayed directly. Milk can be measured directly or with an internal standard.

For perchloric acid:potassium hydroxide clarification, add 10  $\mu$ L of 60% perchloric acid to 100  $\mu$ L of sample, mix well, and centrifuge for 5 minutes at 14k rpm. Add 40  $\mu$ L of 2 M KOH to 100  $\mu$ L of supernatant, allow to precipitate on ice, and then centrifuge for 2 minutes at 14k rpm to collect the supernatant. The dilution factor  $n$  is 1.54.

*Note: SH-containing reagents (e.g.  $\beta$ -mercaptoethanol, or dithiothreitol, > 5  $\mu$ M), sodium azide, EDTA, and sodium dodecyl sulfate are known to interfere in this assay and should be avoided in sample preparation.*

##### Fluorimetric Assay:

1. Equilibrate all components to room temperature. Briefly centrifuge the tubes before opening. Keep thawed tubes on ice during assay. *Important: The Substrate Mix may be slightly turbid, but the turbidity does not interfere with the reaction.*

2. **Standards:** Prepare a 100  $\mu$ M stock of standard by diluting 1  $\mu$ L of the 100 mM Standard with 999  $\mu$ L of dH<sub>2</sub>O. Dilute the 100  $\mu$ M standard in Assay Buffer as follows:

No	100 $\mu$ M STD + dH <sub>2</sub> O	Vol ( $\mu$ L)	L-Carnitine ( $\mu$ M)
1	100 $\mu$ L + 0 $\mu$ L	100	100
2	60 $\mu$ L + 40 $\mu$ L	100	60
3	30 $\mu$ L + 70 $\mu$ L	100	30
4	0 $\mu$ L + 100 $\mu$ L	100	0

Transfer 10  $\mu$ L of each standard into separate wells of a black 96-well plate.

**Samples:** Transfer 10  $\mu$ L of each sample into separate wells of the plate.

Samples requiring an internal standard will need three separate reactions: 1) Sample plus Standard, 2) Sample alone and 3) Sample Blank. For the internal standard, prepare 500  $\mu$ L of 250  $\mu$ M L-Carnitine standard by mixing 2.5  $\mu$ L of 100 mM Standard and 997.5  $\mu$ L of dH<sub>2</sub>O. For the Sample and Sample Blank, add 10  $\mu$ L of sample to the Sample and Sample Blank wells, respectively. For the Sample plus Standard well, add 1  $\mu$ L of 1 mM Standard to 40  $\mu$ L of sample, mix well, and transfer 10  $\mu$ L to the well.

- Fluorimetric Reaction.** Prepare enough Working Reagent by mixing, for each well, 95  $\mu$ L Assay Buffer, 1  $\mu$ L Enzyme A, 1  $\mu$ L Enzyme B, 1  $\mu$ L Enzyme C, 1  $\mu$ L Dye Reagent and 4  $\mu$ L Substrate Mix. Add 90  $\mu$ L of WR to each well, tap to mix and incubate at room temperature (RT) for 30 min protected from light. If using an Internal Standard, prepare a Sample Blank Working Reagent with all the components for the Working Reagent except Enzyme A. Add 90  $\mu$ L of Sample Blank Working Reagent to the Sample Blank wells.
- Read fluorescence ( $\lambda_{ex/em} = 530/585$  nm) at immediately ( $t_{0min}$ ) and at 30 min ( $t_{30min}$ ), or record kinetics for 30 min.

##### Colorimetric Assay:

The colorimetric assay procedure is similar to the fluorimetric procedure except that (1) 0, 300, 600 and 1000  $\mu$ M Standards and (2) a clear 96-well plate are used. Read optical density at 570 nm.

#### CALCULATION

For samples without background matrix effects, use values obtained at 30min. Subtract Blank value (Standard #4) from the standard values and plot the  $\Delta OD$  or  $\Delta F$  against standard concentrations. Determine the slope and calculate the L-Carnitine concentration of Sample,

$$[\text{L-Carnitine}] = \frac{R_s - R_{BL}}{\text{Slope } (\mu\text{M}^{-1})} \times n (\mu\text{M})$$

For samples with background matrix effects (e.g. milk and serum), an internal standard should be used, the sample L-Carnitine concentration is calculated as follows:

$$[\text{L-Carnitine}] = \frac{\Delta R_s - \Delta R_{SB}}{\Delta R_{IS} - \Delta R_s} \times n \times 25 (\mu\text{M})$$

Where  $\Delta R_s$ ,  $\Delta R_{SB}$ , and  $\Delta R_{IS}$  are the fluorescence or optical density readings of the Sample, Sample Blank and Sample plus Internal Standard, respectively, at  $t_{30min} - t_{0min}$ .  $n$  is the sample dilution factor. If the standard curve is non-linear, use polynomial fitting to solve for the concentration.

*Note:* If the calculated L-Carnitine concentration of a sample is higher than 1000  $\mu$ M in the Colorimetric Assay or 100  $\mu$ M in the Fluorimetric Assay, dilute sample in dH<sub>2</sub>O and repeat the assay. Multiply result by the dilution factor,  $n$ .

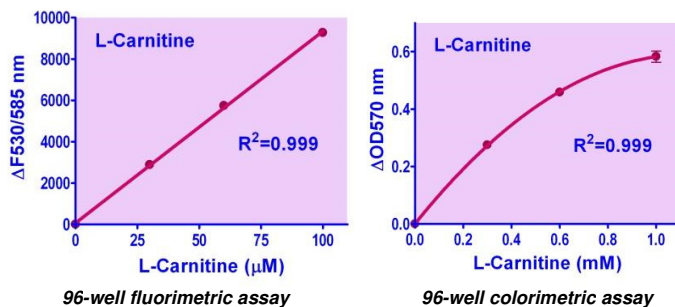
#### MATERIALS REQUIRED, BUT NOT PROVIDED

Pipetting devices, centrifuge tubes, clear flat-bottom uncoated 96-well plates (e.g. VWR cat# 82050-760), optical density plate reader, black flat-bottom uncoated 96-well plates (e.g. VWR cat# 82050-676), fluorescence plate reader. For milk, serum, and plasma samples: 60% perchloric acid and potassium hydroxide.



## EXAMPLES

A PANC-1 cell lysate sample was assayed directly at 0.99 mg/mL protein concentration. The calculated L-Carnitine concentration was 23.2 +/-5.2 nmoles/mg total protein. Human serum and whole cow's milk were assayed using the perchloric acid:potassium hydroxide precipitation and internal standard method. The L-Carnitine concentrations were found to be 46.3 ± 5.9 µM (human serum) and 313.7 ± 85.1 µM (whole cow's milk), respectively.



## LITERATURE

1. Pekala J, et al. (2011). L-carnitine-metabolic functions and meaning in human life. *Curr. Drug Metab.* 12:667-78
2. Almannai M, et al. (1980). Carnitine inborn errors of metabolism. *Molecules.* 24: 3251-3267.
3. Longo N, et al. (2016). Carnitine transport and fatty acid oxidation. *Biochim Biophys Acta.* 1863:2422-35.

