

## EnzyChrom™ Ascorbic Acid Assay Kit (EASC-100)

### Quantitative Colorimetric/Fluorometric Ascorbic Acid Determination

#### DESCRIPTION

Ascorbic acid (the *L*-enantiomer commonly known as vitamin C) is an important antioxidant found in living organisms and applied as additives in food and other industrial processes. By reacting with reactive oxygen species, it protects the cell from oxidative damages. BioAssay Systems' method provides a simple, direct and high-throughput assay for measuring ascorbic acid. In this assay, ascorbic acid is oxidized by ascorbate oxidase resulting in the production of H<sub>2</sub>O<sub>2</sub> which reacts with a specific dye to form a pink colored product. The color intensity at 570nm or fluorescence intensity (530/585 nm) is directly proportional to the ascorbic acid concentration in the sample.

#### KEY FEATURES

Use 20 μL samples. Linear detection range: colorimetric assay 6 to 1000 μM, fluorimetric assay 1 to 100 μM ascorbic acid.

#### APPLICATIONS

**Assays:** ascorbic acid in biological samples such as serum, plasma, urine, saliva, milk, tissue, and cell culture.

**Drug Discovery/Pharmacology:** effects of drugs on ascorbic acid metabolism.

#### KIT CONTENTS

**Assay Buffer:** 10 mL                      **Enzyme Mix:** 120 μL  
**Dye Reagent:** 120 μL                    **Standard:** 400 μL 10 mM ascorbic acid

**Storage conditions.** The kit is shipped on ice. Store all components at -20°C. Shelf life of 12 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.

#### COLORIMETRIC ASSAY

*Note: SH-containing reagents (e.g. β-mercaptoethanol, dithiothreitol, > 5 μM) are known to interfere in this assay and should be avoided in sample preparation.*

**Sample treatment:** liquid samples such as serum and plasma can be assayed directly. Tissue and cell (10<sup>6</sup>-10<sup>7</sup>) lysates can be prepared by homogenization in cold 1 x PBS and centrifugation (5 min at 14,000 rpm). Use clear supernatants for assay. Milk samples should be cleared by mixing 600 μL milk with 100 μL 6 N HCl. Centrifuge 5 min at 14,000 rpm. Transfer 300 μL supernatant into a clean tube and neutralize with 50 μL 6 N NaOH. The neutralized supernatant is ready for assay (dilution factor *n* = 1.36).

- Equilibrate all components to room temperature. Briefly centrifuge the tubes before opening. Keep thawed tubes on ice during assay.
- Standards:** mix 22 μL 10 mM Standard with 198 μL dH<sub>2</sub>O (final 1000 μM). Dilute standard in dH<sub>2</sub>O as follows.

No	1000 μM STD + H <sub>2</sub> O	Vol (μL)	Ascorbic acid (μM)
1	100 μL + 0 μL	100	1000
2	60 μL + 40 μL	100	600
3	30 μL + 70 μL	100	300
4	0 μL + 100 μL	100	0

Transfer 20 μL diluted standards into separate wells of a clear flat-bottom 96-well plate.

**Samples:** transfer 20 μL of each sample into separate wells of the plate.

- Color reaction.** Prepare enough Working Reagent by mixing, for each reaction well, 85 μL Assay Buffer, 1 μL Enzyme Mix and 1 μL Dye Reagent. Add 80 μL Working Reagent to each well. Tap plate to mix. Incubate 10 min at room temperature.

- Read optical density at 570nm (550-585nm).

#### FLUORIMETRIC ASSAY

The fluorimetric assay procedure is similar to the Colorimetric Assay except that (1) 0, 30, 60 and 100 μM ascorbic acid standards and (2) a black 96-well plate are used. Read fluorescence intensity at λ<sub>ex</sub> = 530 nm and λ<sub>em</sub> = 585 nm.

*Note:* if the calculated Ascorbic acid concentration of a sample is higher than 1000 μM in the Colorimetric Assay or 100 μM in the Fluorimetric Assay, dilute sample in water and repeat the assay. Multiply result by the dilution factor *n*.

#### CALCULATION

Subtract blank value (#4) from the standard values and plot the ΔOD or ΔF against standard concentrations. Determine the slope and calculate the ascorbic acid concentration of Sample,

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

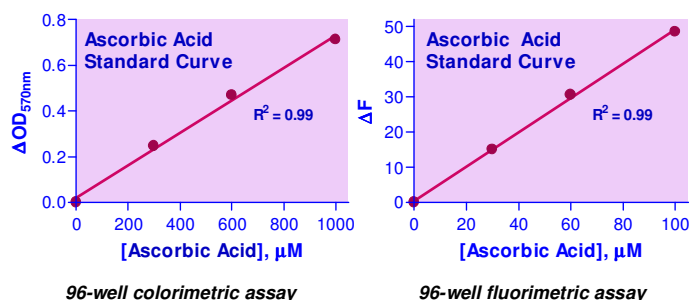
R<sub>SAMPLE</sub> and R<sub>BLANK</sub> are optical density or fluorescence intensity readings of the Sample and H<sub>2</sub>O Blank, respectively. *n* is the sample dilution factor.

**Conversions:** 1 mM ascorbic acid equals 17.6 mg/dL, 0.0176% or 176 ppm.

#### MATERIALS REQUIRED, BUT NOT PROVIDED

Pipetting devices, centrifuge tubes, clear flat-bottom uncoated 96-well plates, optical density plate reader; black flat-bottom uncoated 96-well plates, fluorescence plate reader.

#### Ascorbic acid Standard Curves



#### PUBLICATIONS

- Susanto, L., Y. Siregar, and L. Kusumawati (2018). Vitamin C supplementation improve the sputum conversion culture rate in pulmonary tuberculosis treatment while rifampicin susceptible. IOP Conference Series: Earth and Environmental Science 125(1).
- Kantar, Michael B., et al (2016). Vitamin variation in Capsicum Spp. provides opportunities to improve nutritional value of human diets. PloS one 11.8: e0161464.
- rezzi, Jean-Pierre, et al (2016). LacaScore: a novel plasma sample quality control tool based on ascorbic acid and lactic acid levels. Metabolomics 12.6: 96.

