

EnzyChrom™ Creatine Kinase Assay Kit (ECPK-100)

Colorimetric Determination of Creatine Kinase Activity at 340 nm

DESCRIPTION

CREATINE KINASE (CK), also known as creatine phosphokinase (CPK), is an enzyme (EC 2.7.3.2) expressed predominantly in skeletal muscle, smooth muscle and the brain. The CK enzyme consists of two subunits, which can be either B (brain type) or M (muscle type), and hence three different isoenzymes: CK-MM, CK-BB and CK-MB. CK catalyzes the conversion of creatine to phosphocreatine, consuming adenosine triphosphate (ATP) and generating adenosine diphosphate (ADP) and the reverse reaction. CK is often determined routinely in emergency patients with chest pain and acute renal failure. Elevation of CK is an indication of damage to muscle and has been associated with injury, rhabdomyolysis, myocardial infarction, myositis, myocarditis, malignant hyperthermia and neuroleptic malignant syndrome, etc. Lower levels can be an indication of alcoholic liver disease and rheumatoid arthritis.

Simple, direct and automation-ready procedures for measuring CK activity are very desirable. BioAssay Systems' QuantiChrom™ Creatine Kinase Assay Kit is based on enzyme coupled reactions in which creatine phosphate and ADP is converted to creatine and ATP by CK, the generated ATP is used to phosphorylate glucose by hexokinase to generate glucose-6-phosphate, which is then oxidized by NADP in the presence of glucose-6-phosphate dehydrogenase. The produced NADPH, measured at 340 nm, is proportionate to the CK activity in the sample.

APPLICATIONS

Direct Assays: CK in serum, plasma and other biological samples.

Pharmacology: effects of drugs on CK activity.

KEY FEATURES

Sensitive and accurate. Detection range: 5 to 300 U/L creatine kinase in 96-well plate assay.

Convenient. The procedure involves adding a single working reagent, and reading the optical density at 10 min and 40 min at room temperature or 37°C.

High-throughput. Can be readily automated as a high-throughput 96-well plate assay for thousands of samples per day.

KIT CONTENTS (100 tests in 96-well plates)

Assay Buffer: 12 mL **Substrate Solution:** 1.0 mL
Enzyme Mix: 120 µL **Calibrator:** 150 µL

Storage conditions. This kit is shipped on dry ice. Store all reagents at -20°C. Shelf life of at least 6 months (see expiry dates on labels).

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

Sample Preparation. Use non-hemolyzed samples. Samples should be assayed within 4 hours of blood collection if they remain at room temperature, or within 12 hours if stored at 4°C. Samples can be stored at -20 to -80°C for up to one month. If turbidity is observed, centrifuge sample and use clear supernatant for assay.

Reagent Reconstitution. Bring all components to room temperature. For each reaction well, mix 10 µL Substrate Solution, 100 µL Assay Buffer and 1 µL Enzyme Mix. Fresh reconstitution is recommended. If the assay is to be carried out at 37°C, warm up the Reconstituted Reagent at 37°C.

1. **Calibrator:** transfer 110 µL water and (10 µL Calibrator + 100 µL water) into separate wells of a clear bottom 96-well plate.

Samples: transfer 10 µL samples into separate wells. Add 100 µL Reconstituted Reagent and tap plate to mix.

2. **Reaction.** Incubate at room temperature or 37°C. CK is fully activated within 10 min by glutathione provided in the Substrate Solution. Read OD_{340nm} at 10 min and again at 40 min.

3. **Calculation.** Calculate sample CK activity using the equation,

$$\text{CK (U/L)} = \frac{\text{OD}_{40\text{min}} - \text{OD}_{10\text{min}}}{\text{OD}_{\text{CALIBRATOR}} - \text{OD}_{\text{H}_2\text{O}}} \times 100$$

OD_{40min} and OD_{10min} are OD_{340nm} values at 40 min and 10 min for the sample. OD_{CALIBRATOR} and OD_{H₂O} are OD_{340nm} values of the Calibrator and water blank at 40 min. The value 100 is the equivalent activity (U/L) of the Calibrator under the assay conditions.

Unit definition: one unit of CK will transfer 1 µmole of phosphate from phosphocreatine to ADP per min at pH 6.0.

Note: If the calculated CK activity is higher than 300 U/L, dilute sample in 0.9% saline and repeat this assay. Multiply the results by the dilution factor.

MATERIALS REQUIRED, BUT NOT PROVIDED

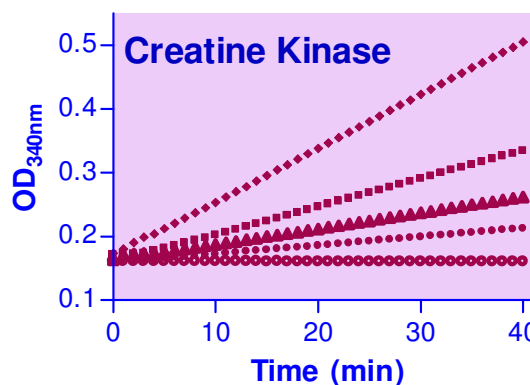
Pipeting (multi-channel) devices. Clear-bottom 96-well plates (e.g. Corning Costar) and plate reader.

GENERAL CONSIDERATIONS

This assay is based on an enzyme-catalyzed kinetic reaction. Addition of Reconstituted Reagent should be quick and mixing should be brief but thorough. Use of multi-channel pipettor is recommended.

EXAMPLES

Samples were assayed in duplicate using the 96-well protocol. The CK activity (U/L) was 12.0 ± 0.9 for a rat serum sample, 11.0 ± 0.5 for human serum, 28 ± 1 for human plasma, 9.0 ± 0.8 for mouse serum and 49 ± 2 for bovine serum.



Kinetics of CK Reaction at 25 (solid circle), 50 (triangle), 100 (square) and 200 (diamond) U/L. Control: open circle

LITERATURE

- [1]. C. Bishop, T. M. Chu, and Z. K. Shihabi (1971). Single Stable Reagent for Creatine Kinase Assay. *Clin. Chem.* 17 (6): 548-550.
- [2]. G Szasz, W Gerhardt, W Gruber, and E Bernt (1976). Creatine kinase in serum: 2. Interference of adenylate kinase with the assay. *Clin. Chem.* 22: 1806 - 1811.
- [3]. G Szasz, W Gruber, and E Bernt (1976). Creatine kinase in serum: 1. Determination of optimum reaction conditions. *Clin. Chem.*, 22: 650 - 656.