

Phosphoenolpyruvic acid (PEP) Assay Kit ab204713

[1 References](#) [2 Images](#)

Overview

Product name	Phosphoenolpyruvic acid (PEP) Assay Kit
Detection method	Colorimetric/Fluorometric
Sample type	Urine, Serum, Plasma, Other biological fluids, Cell Lysate, Tissue Lysate
Assay type	Enzyme activity
Sensitivity	1 μ M
Product overview	Phosphoenolpyruvic acid (PEP) Assay Kit (ab204713) provides a convenient colorimetric and fluorometric means to measure PEP levels in various samples. In the assay, PEP is converted to ATP and pyruvate. The generated pyruvate is quantified by colorimetric ($\lambda_{max} = 570$ nm) or fluorometric methods (Ex/Em = 535/587 nm). The assay is simple, sensitive and reliable. The detection limit is approximately 1 μ M PEP in biological samples.

Notes This product is manufactured by BioVision, an Abcam company and was previously called K365 PEP Colorimetric/Fluorometric Assay Kit. K365-100 is the same size as the 100 test size of ab204713.

Platform Microplate reader

Properties

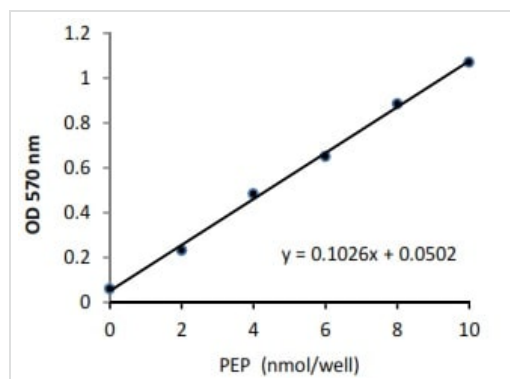
Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests
PEP Assay Buffer	1 x 25ml
PEP Converter	1 vial
PEP Developer Mix	1 vial
PEP Probe (in DMSO)	1 x 0.2ml
PEP Standard (1 μ mole)	1 vial

Relevance

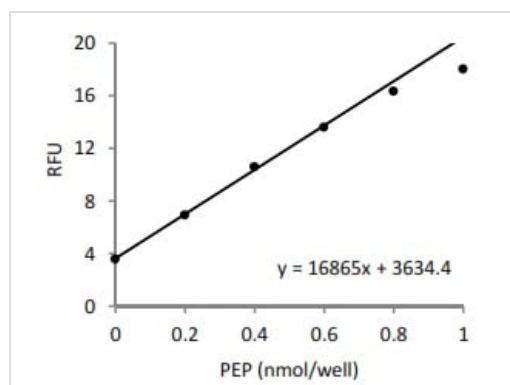
Phosphoenolpyruvate (PEP) is an important intermediate in carbohydrate metabolism. Containing a high-energy phosphate bond, PEP is involved in glycolysis and gluco-neogenesis and in the shikimate pathway as well as in carbon fixation in plants. Bacteria utilize PEP in the phosphor-transferase system to acquire sugars from the environment. In the glycolytic pathway, PEP is formed from 2-phosphoglycerate by enolase and generates ATP through the action of pyruvate kinase.

Images



Typical PEP Standard calibration curve using colorimetric reading.

Phosphoenolpyruvic acid (PEP) Assay Kit
(ab204713)



Typical PEP Standard calibration curve using fluorometric reading.

Phosphoenolpyruvic acid (PEP) Assay Kit
(ab204713)

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you

- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors