

Version 3a, Last updated 6 June 2025

# **ab204713**

## **Phosphoenolpyruvic acid (PEP) Assay Kit**

For the measurement of PEP levels in various samples.

This product is for research use only and is not intended for diagnostic use.

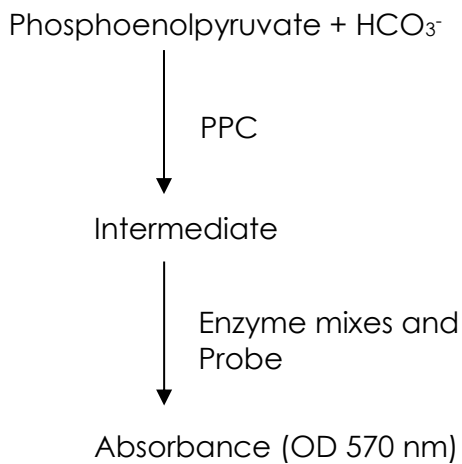
## Table of Contents

1. Overview	3
2. Protocol Summary	4
3. General guidelines, precautions, and troubleshooting	5
4. Materials Supplied, and Storage and Stability	6
5. Materials Required, Not Supplied	6
6. Reagent Preparation	7
7. Standard Preparation	8
8. Sample Preparation	10
9. Assay Procedure	11
10. Calculations	11
11. Typical Data	13
12. Notes	14

## 1. 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 ( $\lambda_{\text{max}} = 570 \text{ nm}$ ) or fluorometric methods (Ex/Em = 535/587 nm). The assay is simple, sensitive and reliable. The detection limit is approximately 1  $\mu\text{M}$  PEP in biological samples.

## 2. Protocol Summary



### **3. General guidelines, precautions, and troubleshooting**

- Please observe safe laboratory practice and consult the safety datasheet.
- For general guidelines, precautions, limitations on the use of our assay kits and general assay troubleshooting tips, particularly for first time users, please consult our guide:  
[www.abcam.com/assaykitguidelines](http://www.abcam.com/assaykitguidelines)
- For typical data produced using the assay, please see the assay kit datasheet on our website.

## 4. Materials Supplied, and Storage and Stability

- Store kit at -20°C in the dark immediately upon receipt and check below in Section 6 for storage for individual components. Kit can be stored for 1 year from receipt, if components have not been reconstituted.
- Aliquot components in working volumes before storing at the recommended temperature (see section 6).

Item	Quantity	Storage condition
Assay Buffer 4	25 mL	-20°C
OxiRed™ Probe	0.2 mL	-20°C
PEP	1 vial	-20°C
PEP Converter Mix	1 vial	-20°C
Developer Mix A	1 vial	-20°C

PLEASE NOTE: Assay Buffer 4 was previously labelled as Assay Buffer IV and PEP Assay Buffer, and OxiRed™ Probe as OxiRed Probe and PEP Probe (in DMSO). Developer Mix A was previously labelled as Development Enzyme Mix I and PEP Developer Mix. The composition has not changed.

## 5. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- 3M HClO<sub>4</sub>
- 3M KHCO<sub>3</sub>
- activated charcoal

## 6. Reagent Preparation

- Before using the kit, spin tubes and bring down all components to the bottom of tubes.
- Prepare only as much reagent as is needed on the day of the experiment.

### 6.1 Assay Buffer 4:

Bring to room temperature before use.

### 6.2 OxiRed™ Probe:

Ready to use as supplied. Warm for 1-2 min at 37°C to melt the frozen DMSO before use. Mix well. Store at –20°C, protect from light and moisture. Use within two months.

### 6.3 PEP:

Dissolve in 100 µL dH<sub>2</sub>O to generate 10 mM stock solution. Keep cold while in use. Store at -20°C.

### 6.4 PEP Converter Mix:

Dissolve with 220 µL Assay Buffer 4. Pipette up and down to dissolve. Store at –20°C. Use within two months.

### 6.5 Developer Mix A:

Dissolve with 220 µL Assay Buffer 4. Pipette up and down to dissolve. Store at –20°C. Use within two months.

## 7. Standard Preparation

- Always prepare a fresh set of standards for every use.
- Discard working standard dilutions after use as they do not store well.

### 7.1 For Colorimetric Assay:

- 7.1.1 Prepare 1 mM (1 nmol/μL) PEP Assay Standard by adding 10 μL of the 10 mM Standard to 90μL of Assay Buffer 4, mix well.
- 7.1.2 Add 0, 2, 4, 6, 8, 10 μL into a series of wells.
- 7.1.3 Adjust volume to 50 μL/well with Assay Buffer 4 to generate 0, 2, 4, 6, 8, 10 nmol/well of PEP.

Standard #	1 mM PEP Assay Standard (μL)	Assay Buffer 4 (μL)	PEP nmoles/well
1	10	40	10
2	8	42	8
3	6	44	6
4	4	46	4
5	2	48	2
6	0	50	0

### 7.2 For Fluorometric Assay:



- 7.2.1 Prepare 0.1 mM (0.1 nmol/ $\mu$ L) PEP Assay Standard by adding 1  $\mu$ L of the 10 mM Standard to 99  $\mu$ L of Assay Buffer 4, mix well (the fluorometric assay is 10 to 100 fold more sensitive than the colorimetric assay).
- 7.2.2 Add 0, 2, 4, 6, 8, 10  $\mu$ L into a series of standards wells.
- 7.2.3 Adjust volume to 50  $\mu$ L/well with Assay Buffer 4 to generate 0, 0.2, 0.4, 0.6, 0.8, 1.0 nmol/well of PEP.

Standard #	0.1 mM PEP Assay Standard ( $\mu$ L)	Assay Buffer 4 ( $\mu$ L)	PEP nmoles/well
1	10	40	1.0
2	8	42	0.8
3	6	44	0.6
4	4	46	0.4
5	2	48	0.2
6	0	50	0

## 8. Sample Preparation

- 8.1** Tissue (20-50 mg) should be frozen in liquid N<sub>2</sub> or dry ice/MeOH then powdered thoroughly with mortar and pestle at -80°C and transferred to an Eppendorf tube.
- 8.2** Add 100 µL ice cold HClO<sub>4</sub> and vortex until the contents are thoroughly mixed.
- 8.3** Neutralize carefully by adding repeated small aliquots (~10 µL per aliquot) of KHCO<sub>3</sub> followed by vortexing. Final pH should be 6.5-7.5.
- 8.4** Centrifuge at 12,000 g for 3 min.

**Δ Note:** Some plant extracts need to be decolorized with activated charcoal (5 mg per tube) which can be added and vortexed prior to the centrifugation step.

- 8.5** Use samples immediately or store at -80°C.

- 8.6** Add up to 50 µL sample per well in a 96-well plate; bring the volume to 50 µL with Assay Buffer 4.

**Δ Note:** Intracellular PEP level is usually in the range of 0.05 - 0.3 mM. We suggest testing several doses of your sample to ensure the readings are within the standard curve range.

## 9. Assay Procedure

- 9.1** Mix enough reagents for the number of assays to be performed. For each well, prepare 50 µL mix containing:

	Reaction Mix	
	Colorimetric Assay	Fluorometric Assay
Assay Buffer 4	44 µL	45.8 µL
OxiRed™ Probe	2 µL	0.2 µL*
PEP Converter Mix **	2 µL	2 µL
Developer Mix A	2 µL	2 µL

- 9.2** Mix and add 50 µL of the Reaction Mix to each well containing the Standard Curve and test samples.
- 9.3** Incubate at room temperature for 1 h, protected from light.
- 9.4** Measure O.D. 570 nm for colorimetric assay or Ex/Em 535/587 nm for fluorometric assay.

### Δ Note:

- \*For the fluorometric assay, use 1/10 of OxiRed™ Probe to reduce fluorescence background.
- \*\*Pyruvate generates background.
  - If significant amount of pyruvate is suspected in your samples, a sample pyruvate background control needs to be performed by replacing the PEP Converter Mix with 2 µL of Assay Buffer 4.
  - Then follow the same protocol as the sample.
  - In the absence of PEP Converter Mix, the assay detects only pyruvate, not PEP. The pyruvate background reading can be subtracted from the PEP readings.

## 10. Calculations

- 10.1** Subtract 0 PEP control reading from all readings. Plot the PEP Assay Standard Curve.

**10.2** If a sample pyruvate background control was used, subtract the background control reading from the PEP readings.

**10.3** Apply the sample readings to the standard curve to get PEP amount in the sample wells.

**10.5** The PEP concentrations in the test samples:

$C = A_y/S_v$  (nmol/  $\mu$ L,  $\mu$ mol/mL or mM)

Where:

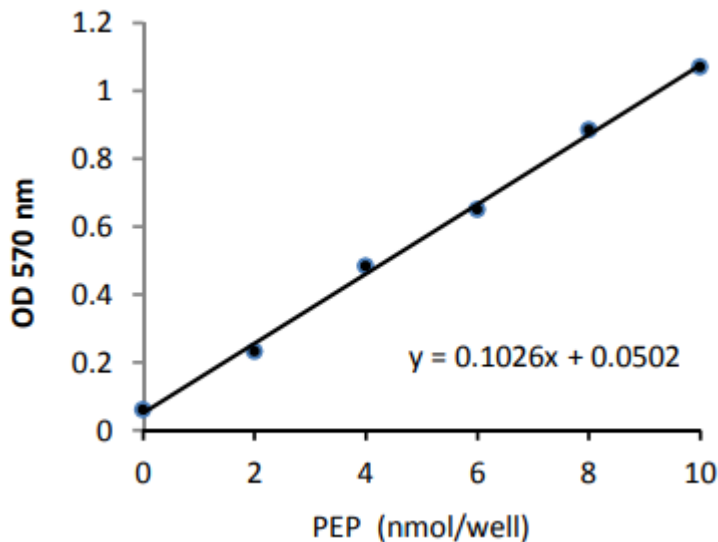
$A_y$  is the PEP amount (nmol) in your sample from the standard curve.

$S_v$  is the sample volume ( $\mu$ L) added to the assay well.

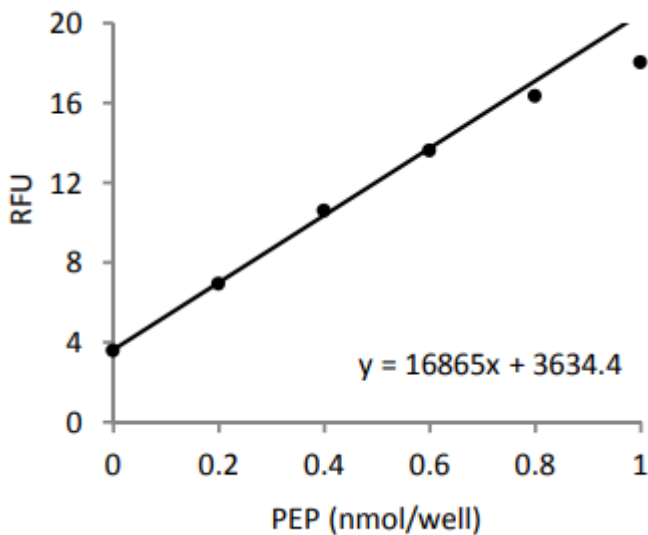
PEP molecular weight: 168.04.

# 11. Typical Data

Typical data provided for demonstration purposes only.



**Figure 1.** Colorimetric Assay standard curve.



**Figure 2.** Fluorometric Assay standard curve.

## 12. Notes



## **Technical Support**

Copyright © 2025 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)