

Version 3b, Last updated 16 July 2026

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.

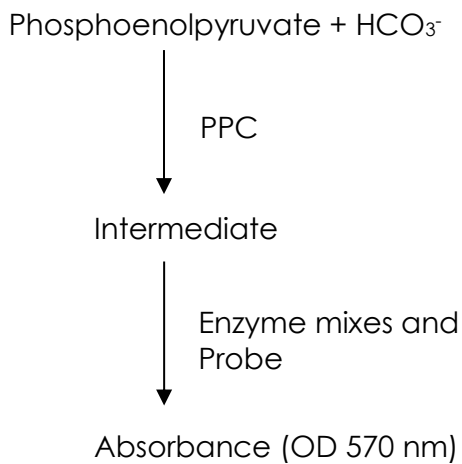
Table of Contents

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 μM 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
- For typical data produced using the assay, please see the assay kit datasheet on our website.
- If applicable, please refer to the current Safety Data Sheet (SDS) provided with this product for safety, handling, and disposal information. The most up to date and current versions are available on our website <https://www.abcam.com/en-us>.

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.
- 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 HClO₄
- 3M KHCO₃
- 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 the DMSO solution at 37 °C until completely thawed (~ 5 minutes), vortex briefly to create a homogenous solution, and spin down briefly to collect any contents trapped in the cap or seal. Keep at room temperature during use and reseal promptly to minimize moisture exposure.

For multiple use aliquot and store at –20 °C, protected from light and moisture. Stable for up to 2 months after thawing (≤3 freeze–thaw cycles).

6.3 PEP:

Dissolve in 100 µL dH₂O 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/ μ L) PEP Assay Standard by adding 1 μ L of the 10 mM Standard to 99 μ 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 μ L into a series of standards wells.
- 7.2.3 Adjust volume to 50 μ 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 (μ L)	Assay Buffer 4 (μ 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 N₂ 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 HClO₄ and vortex until the contents are thoroughly mixed.
- 8.3** Neutralize carefully by adding repeated small aliquots (~10 µL per aliquot) of KHCO₃ 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/ μ L, μ 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 (μ L) added to the assay well.

PEP molecular weight: 168.04.

11. Typical Data

Typical data provided for demonstration purposes only.

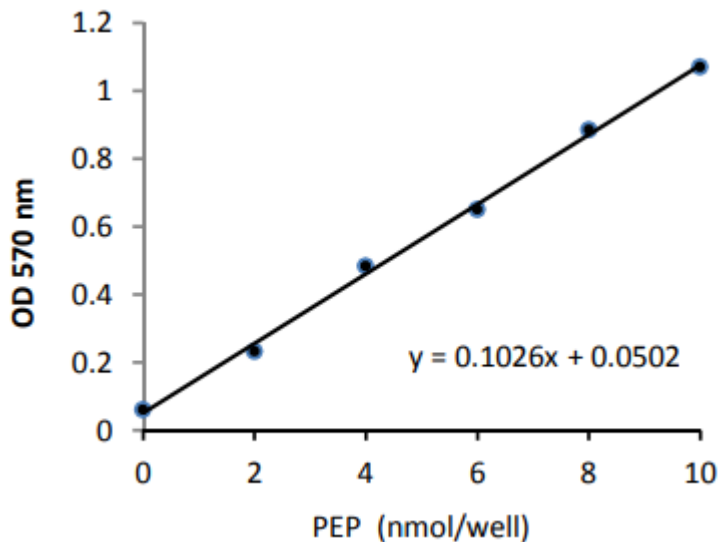


Figure 1. Colorimetric Assay standard curve.

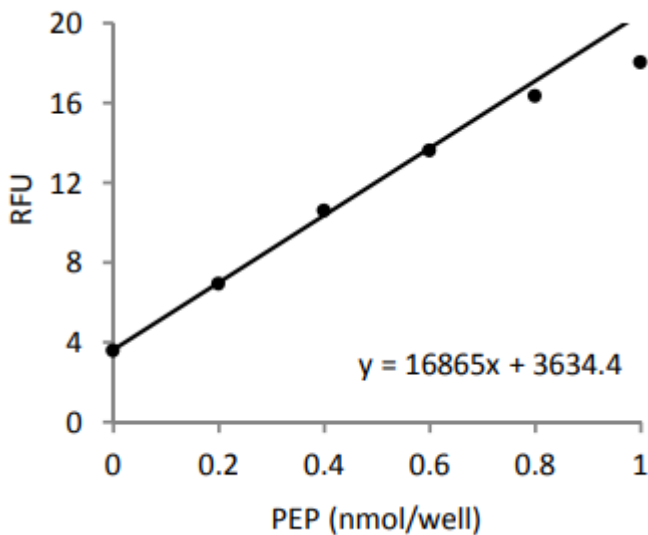


Figure 2. Fluorometric Assay standard curve.

12. Notes

Technical Support

Copyright © 2026 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:

<https://www.abcam.com/en-us/contact-us>

<https://www.abcam.cn/contact-us> (China)

<https://www.abcam.co.jp/contact-us> (Japan)