

ab83426

Glucose 6 Phosphate Assay Kit (Colorimetric)

Instructions for Use

For the rapid, sensitive and accurate measurement of Glucose 6 Phosphate levels in various samples

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

PLEASE NOTE: With the acquisition of BioVision by Abcam, we have made some changes to component names and packaging to better align with our global standards as we work towards environmental-friendly and efficient growth. You are receiving the same high-quality products as always, with no changes to specifications or protocols.

Table of Contents

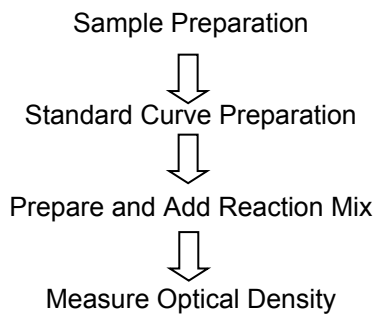
| | |
|---------------------------|----|
| 1. Overview | 3 |
| 2. Protocol Summary | 4 |
| 3. Components and Storage | 4 |
| 4. Assay Protocol | 6 |
| 5. Data Analysis | 10 |
| 6. Troubleshooting | 10 |

1. Overview

Glucose 6 Phosphate is a key sugar intermediate for glucose to get into cells, and then enter either metabolic pathways or storage. Glucose 6 Phosphate can enter the glycolytic pathway, the pentose phosphate shunt or be stored as glycogen or starch. Glucose 6 Phosphate is utilized by its dehydrogenase to generate reducing equivalents in the form of NADPH. This is particularly important in red blood cells where a G6PDH deficiency leads to hemolytic anemia.

Abcam's Glucose 6 Phosphate Assay Kit is a simple, sensitive and rapid means of quantifying Glucose 6 Phosphate in a variety of samples. In the assay, Glucose 6 Phosphate is oxidized with the generation of a product which is utilized to convert a nearly colorless probe to an intensely colored product with an absorbance at 450 nm. The Glucose 6 Phosphate Assay Kit can detect Glucose 6 Phosphate in the range of 1 to 30 nmol with detection sensitivity ~10 μ M of Glucose 6 Phosphate.

2. Protocol Summary



3. Components and Storage

A. Kit Components

| Item | Quantity |
|---|----------|
| Assay Buffer II/Glucose 6 Phosphate Assay Buffer | 25 mL |
| Development Enzyme Mix IX/Glucose 6 Phosphate Enzyme Mix (Lyophilized) | 1 vial |
| Developer Solution III/Glucose 6 Phosphate Substrate Mix (Lyophilized) | 1 vial |
| G6P Standard/Glucose 6 Phosphate Standard (10 μ mol; Lyophilized) | 1 vial |

Store kit at -20°C , protect from light. Warm Assay Buffer II/Glucose 6 Phosphate Assay Buffer to room temperature before use. Briefly centrifuge all small vials prior to opening. Keep enzyme mix on ice while in use.

G6P STANDARD/GLUCOSE 6 PHOSPHATE STANDARD: Dissolve in 100 μL dH_2O to generate 100 mM (100 nmol/ μL) G6P Standard/Glucose 6 Phosphate Standard solution. Keep cold while in use. Store at -20°C .

DEVELOPER SOLUTION III/GLUCOSE 6 PHOSPHATE
SUBSTRATE MIX: Dissolve with 220 μL of Assay Buffer II/Glucose 6 Phosphate Assay Buffer. Pipette up and down to dissolve. Stable for 2 months at $+4^{\circ}\text{C}$.

DEVELOPMENT ENZYME MIX IX/GLUCOSE 6 PHOSPHATE
ENZYME MIX: Dissolve with 220 μL dH_2O . Pipette up and down to dissolve. Aliquot into portions and store at -20°C . Avoid repeated freeze/thaw cycles. Use within two months.

B. Additional Materials Required

- Microcentrifuge
- Pipettes and pipette tips
- colorimetric microplate reader
- 96 well plate
- Orbital shaker

4. Assay Protocol

1. Sample Preparation:

- a. **Liquid or solution samples** can be assayed directly.
- b. **For tissue or cell samples:** 10-100 mg tissue or 5 million cells should be rapidly homogenized with 2-3 volume of ice cold PBS or other buffer (pH 6.5-8). Centrifuge at top speed for 10 min to remove insoluble materials.

Add 1-50 μL samples into duplicate wells of a 96-well plate and bring volume to 50 μL with Assay Buffer.

For unknown samples, we suggest testing several doses of your samples to ensure readings are within the standard curve range.

Notes:

A. Enzymes in samples may interfere with the assay. We suggest deproteinizing samples using 10 kDa molecular weight cut off spin columns (**ab93349**) or using a perchloric acid/KOH protocol as follows:

- a) Tissue samples (20-1000 mg) should be frozen rapidly (liquid N_2 or methanol/dry ice), weighed and pulverized.
- b) Add 2 μL 1N perchloric acid/mg per sample. KEEP COLD!
- c) Homogenize or sonicate thoroughly. Spin homogenate at 10,000 x g for 5-10 minutes.

- d) Neutralize supernatant with 10N KOH to minimize Glucose 6 Phosphate conversion, adding repeated 1 μL aliquots/10 μL supernate while vortexing. Add until bubble evolution ceases (2-5 aliquots). Put on ice for 5 minutes
- e) Check pH (using 1 μL) is \sim 6-8. Spin 2 minutes at 10,000 x g to pellet KClO_4 .

For tissues or cells containing low level of free Glucose 6 Phosphate (5-60 μM), try to minimize sample dilutions.

B. NADH or NADPH in samples will generate background readings. If NADH or NADPH is in your sample, you may do a background control (omit Development Enzyme Mix IX/Glucose 6 Phosphate Enzyme Mix from the reaction mix) to read the background, then subtract the background from Glucose 6 Phosphate readings.

2. Standard Curve Preparation:

Dilute the G6P Standard/Glucose 6 Phosphate Standard to 1 nmol/ μL by adding 10 μL of the 100 nmol/ μL Standard to 990 μL of dH_2O , mix well. Add 0, 2, 4, 6, 8, 10 μL into a series of standards wells on a 96-well plate.

Adjust volume to 50 μL /well with Assay Buffer to generate 0, 2, 4, 6, 8, 10 nmol/well of G6P Standard/Glucose 6 Phosphate Standard.

3. Reaction Mix: Mix enough reaction mix for the number of samples and standards to be performed: For each well, prepare a total 50 μL Reaction Mix containing:

| | Reaction Mix | Background |
|--------------------|---------------------|-------------------|
| Assay Buffer | 46 μL | 48 μL |
| Enzyme Mix | 2 μL | --- |
| Developer Solution | | |
| III/Substrate Mix | 2 μL | 2 μL |

Add 50 μL of the Reaction Mix to each well containing the G6P Standard/Glucose 6 Phosphate Standard and samples. Add 50 μL of the background mix into background control wells.

4. Incubate for 30 min at room temperature, protect from light.

5. Measure OD at 450 nm

5. Data Analysis

Correct background by subtracting the value of the zero Glucose 6 Phosphate blank from all sample and standard readings. If background control reading is significant, subtract the background reading from sample reading.

Plot the standard curve. Apply the corrected sample readings to the standard curve to get Glucose 6 Phosphate amount in the sample wells. The Glucose 6 Phosphate concentrations in the test samples:

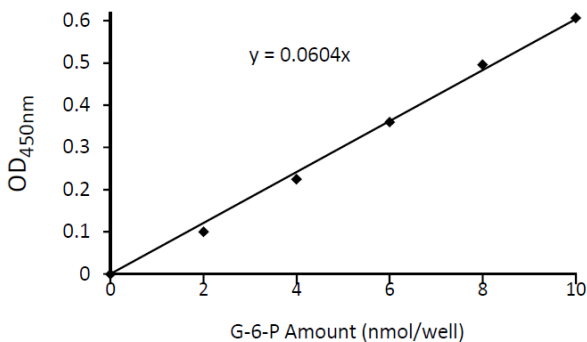
$$\text{Concentration} = \text{Ay} / \text{Sv} \text{ (nmol/}\mu\text{L; or } \mu\text{mol/mL; or mM)}$$

Where:

Ay is the amount of Glucose 6 Phosphate (nmol) in your sample from the standard curve.

Sv is the sample volume (μL) added to the sample well.

Glucose 6 Phosphate molecular weight: 260.14.



Glucose 6 Phosphate standard curve generated using this kit protocol

6. Troubleshooting

| Problem | Reason | Solution |
|--------------------|--|---|
| Assay not working | Assay buffer at wrong temperature | Assay buffer must not be chilled - needs to be at RT |
| | Protocol step missed | Re-read and follow the protocol exactly |
| | Plate read at incorrect wavelength | Ensure you are using appropriate reader and filter settings (refer to datasheet) |
| | Unsuitable microtiter plate for assay | Fluorescence: Black plates (clear bottoms); Luminescence: White plates; Colorimetry: Clear plates. If critical, datasheet will indicate whether to use flat- or U-shaped wells |
| Unexpected results | Measured at wrong wavelength | Use appropriate reader and filter settings described in datasheet |
| | Samples contain impeding substances | Troubleshoot and also consider deproteinizing samples |
| | Unsuitable sample type | Use recommended samples types as listed on the datasheet |
| | Sample readings are outside linear range | Concentrate/ dilute samples to be in linear range |

| | | |
|---|---|---|
| Samples with inconsistent readings | Unsuitable sample type | Refer to datasheet for details about incompatible samples |
| | Samples prepared in the wrong buffer | Use the assay buffer provided (or refer to datasheet for instructions) |
| | Samples not deproteinized (if indicated on datasheet) | Use the 10kDa spin column (ab93349) |
| | Cell/ tissue samples not sufficiently homogenized | Increase sonication time/ number of strokes with the Dounce homogenizer |
| | Too many freeze-thaw cycles | Aliquot samples to reduce the number of freeze-thaw cycles |
| | Samples contain impeding substances | Troubleshoot and also consider deproteinizing samples |
| | Samples are too old or incorrectly stored | Use freshly made samples and store at recommended temperature until use |
| Lower/ Higher readings in samples and standards | Not fully thawed kit components | Wait for components to thaw completely and gently mix prior use |
| | Out-of-date kit or incorrectly stored reagents | Always check expiry date and store kit components as recommended on the datasheet |
| | Reagents sitting for extended periods on ice | Try to prepare a fresh reaction mix prior to each use |
| | Incorrect incubation time/ temperature | Refer to datasheet for recommended incubation time and/ or temperature |
| | Incorrect amounts used | Check pipette is calibrated correctly (always use smallest volume pipette that can pipette entire volume) |

| Problem | Reason | Solution |
|------------------------------|--|--|
| Standard curve is not linear | Not fully thawed kit components | Wait for components to thaw completely and gently mix prior use |
| | Pipetting errors when setting up the standard curve | Try not to pipette too small volumes |
| | Incorrect pipetting when preparing the reaction mix | Always prepare a master mix |
| | Air bubbles in wells | Air bubbles will interfere with readings; try to avoid producing air bubbles and always remove bubbles prior to reading plates |
| | Concentration of standard stock incorrect | Recheck datasheet for recommended concentrations of standard stocks |
| | Errors in standard curve calculations | Refer to datasheet and re-check the calculations |
| | Use of other reagents than those provided with the kit | Use fresh components from the same kit |

Technical Support

Copyright © 2023 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

www.abcam.cn/contactus (China)

www.abcam.co.jp/contactus (Japan)