

ab102517

Glucose Detection Kit

Instructions for Use

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

[View kit datasheet: www.abcam.com/ab102517](http://www.abcam.com/ab102517)

(use www.abcam.cn/ab102517 for China, or www.abcam.co.jp/ab102517 for Japan)

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

Table of Contents

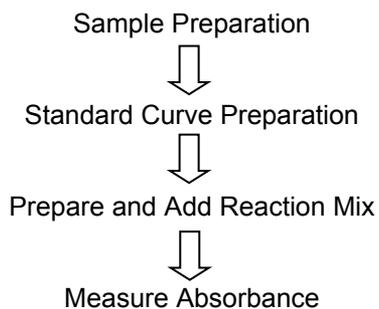
| | |
|---------------------------|----|
| 1. Overview | 3 |
| 2. Protocol Summary | 4 |
| 3. Components and Storage | 5 |
| 4. Assay Protocol | 7 |
| 5. Data Analysis | 9 |
| 6. Troubleshooting | 11 |

1. Overview

Glucose is an important fuel source to generate the universal energy molecule ATP. Serum glucose level is a key diagnostic parameter for many metabolic disorders.

Abcam's Glucose Detection Kit provides direct measurement of glucose in various biological samples (e.g., serum, plasma, other body fluids, food, growth media, etc.). In this assay, glucose is specifically oxidized to generate a product which reacts with a dye to generate color ($\lambda = 450 \text{ nm}$) whose intensity is proportional to glucose concentration. The method is rapid, simple, sensitive, and suitable for high throughput. This assay is particularly suitable for serum and urine samples since it is unaffected by reducing substances which can interfere with other suppliers offering oxidase-based kits. The assay is also suitable for monitoring glucose level during fermentation and glucose feeding in protein expression processes. The kit can detect glucose concentrations in the range of $20\mu\text{M}$ - 10mM .

2. Protocol Summary



3. Components and Storage

A. Kit Components

| Item | Quantity |
|-------------------------------------|-------------|
| Glucose Assay Buffer | 25 mL |
| Glucose Substrate Mix (Lyophilized) | 1 vial |
| Glucose Enzyme Mix (Lyophilized) | 1 vial |
| Glucose Standard (100 mM) | 100 μ L |

* Store kit at -20°C , protect from light.

- Allow Assay Buffer to warm to room temperature before use.
- Briefly centrifuge all small vials prior to opening.
- Read the entire protocol before performing the assay.

GLUCOSE SUBSTRATE MIX AND ENZYME MIX: Dissolve separately in 220 μ L Glucose Assay Buffer. Aliquot and store at -20°C , protect from light and moisture. 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:

Prepare test samples in 50 μl /well with Glucose Assay Buffer in a 96-well plate. If using serum sample, serum (0.5 - 2 μl /assay. Normal serum contains ~ 5 nmol/ μl glucose) can be directly diluted in the Glucose Assay Buffer.

Note:

It is recommended to de-proteinize samples by centrifugation using a 10 kDa spin column (ab93349) to remove enzymes and interfering proteins.

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

2. Standard Curve Preparation:

Dilute the Glucose Standard to 1 nmol/ μl by adding 10 μl of the Glucose Standard to 990 μl of Glucose Assay Buffer, mix well. Add 0, 2, 4, 6, 8, 10 μl into a series of wells of a 96 well plate. Adjust volume of all wells to 50 μl with Glucose Assay Buffer to generate 0, 2, 4, 6, 8, 10 nmol/well of Glucose Standard.

3. Glucose Reaction Mix: Mix enough reagent for the number of assays to be performed. For each well, prepare a total 50 μ l Reaction Mix containing:

| | |
|-----------------------|------------|
| Glucose Assay Buffer | 46 μ l |
| Glucose Enzyme Mix | 2 μ l |
| Glucose Substrate Mix | 2 μ l |

Mix well. Add 50 μ l of the Reaction Mix to each well containing the Glucose Standard and test samples, mix well. Incubate the reaction for 30 min, protect from light.

4. Measurement: Measure absorbance at 450 nm in a microplate reader.

5. Data Analysis

Correct background by subtracting the value derived from the zero glucose control from all readings. The background reading can be significant and must be subtracted from sample readings.

Plot the standard curve. Apply the sample readings to the standard curve. Glucose concentrations of the test samples can then be calculated:

$$\text{Concentration} = \text{Sa} / \text{Sv} \text{ (nmol/}\mu\text{l or mM)}$$

Where:

Sa is sample amount (in nmol) calculated from standard curve.

Sv is sample volume (μl) added to the wells.

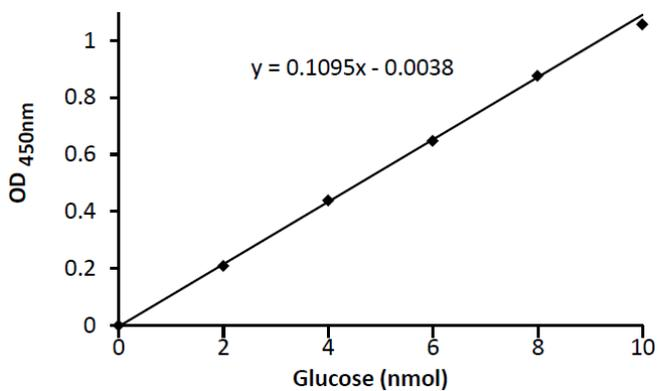
Note:

If sample was pre-diluted before added to reaction well you must correct for this dilution factor

Glucose Molecular Weight: 180.16.

Normal serum glucose range: 3-7 mM.

Normal urine glucose range: 0-0.8 mM.



Standard Curve for Glucose carried out using the 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 |

| Problem | Reason | Solution |
|---|---|---|
| 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 |

UK, EU and ROW

Email: technical@abcam.com | Tel: +44-(0)1223-696000

Austria

Email: wissenschaftlicherdienst@abcam.com | Tel: 019-288-259

France

Email: supportscientifique@abcam.com | Tel: 01-46-94-62-96

Germany

Email: wissenschaftlicherdienst@abcam.com | Tel: 030-896-779-154

Spain

Email: soportecientifico@abcam.com | Tel: 911-146-554

Switzerland

Email: technical@abcam.com
Tel (Deutsch): 0435-016-424 | Tel (Français): 0615-000-530

US and Latin America

Email: us.technical@abcam.com | Tel: 888-77-ABCAM (22226)

Canada

Email: ca.technical@abcam.com | Tel: 877-749-8807

China and Asia Pacific

Email: hk.technical@abcam.com | Tel: 400 921 0189 / +86 21 2070 0500

Japan

Email: technical@abcam.co.jp | Tel: +81-(0)3-6231-0940

www.abcam.com | www.abcam.cn | www.abcam.co.jp