Glucose Assay Kit-WST

Technical Manual

Technical Manual (Japanese version) is available at http://www.dojindo.co.jp/manual/g264.pdf

General Information

Glucose, the most important biological energy source, is a known marker for energy metabolism. It is used not only as a marker for diabetes and obesity but also for monitoring changes in intracellular metabolism, such as lactate production. In addition, a recent study suggested that inhibition of the enzymatic activities involved in both glucose and lipid metabolism effectively inhibits cancer cell proliferation¹⁾.

Glucose Assay Kit-WST enables quantitation of glucose as a substrate in energy metabolism. This kit has been optimized to quantify glucose in cell culture supernatants by measuring absorbance in a colorimetric WST reaction. This kit is formatted for 96-well microplates and the assay has a detection limit of 0.02 mmol/L glucose.



	lactone					
		Figure 1. Principle	of Glucose	Assay Kit-	WST	
Kit Contents		50 tests	200 tests			
	Dye Mixture	×1	×1			
	Glucose Standard (10 mmol/L) (Red cap)	150 μL × 1	600 µL × 1			
	Enzyme (Green cap)	×1	×1			
	Assay Buffer	3.5 mL × 1	14 mL × 1			
	Reconstitution Buffer (Blue cap)	350 μL × 1	1.4 mL × 1			
Storage Conditions	Store at 0−5 °C					
Required Equipment and Materials	- Microplate reader (450 nm filter)- 96-well microplate- Incubator (37 °C)- 20-200 μL multichannel pipette- 100-1000 μL and 20-200 μL micropipettes- Phosphate buffered saline (PBS)- Conical tube- 0.1% Triton X-100					
Precautions	 Equilibrate reagents to room temperature prior to use. Centrifuge the tube with Enzyme (green cap) briefly before opening to remove all content from walls of the tube and inside the cap. Analyzing samples and standards in triplicate is recommended for accuracy. Use a multichannel pipette to dispense working solution, to minimize time lag in pipetting, because the enzymatic reaction is initiated immediately upon its addition. Prepare samples at different dilutions to determine the most suitable concentration, ranging from 0 to 0.5 mmol/L glucose. Dye Mixture is stored in a glass bottle with an aluminum cap. Handle with caution, wearing gloves. The kit was designed for analyzing cell culture supernatants. To instead measure intracellular glucose, prepare cell lysates and glucose standards in a 0.1% Triton X-100 solution. 					
	 Preparation of Dye Mixture stock solution Add the entire volume of Reconstitution Buffer to the Dye Mixture vial and dissolve contents completely. Transfer the resulting Dye Mixture stock solution to the vial from the Reconstitution Buffer and store at 0-5 °C, protected from light. The Dye Mixture stock solution is stable for 2 months under these conditions. 					
	Preparation of Enzyme stock solution Add PBS to the Enzyme tube and dissolve with pipetting. ※ Refer to Table 1. ※ Centrifuge the tube briefly each time before opening to remove all content from the tube wall or inside the cap. ※ Enzyme stock solution should be maintained in an ice bath during use. ※ The Enzyme stock solution is stable for 2 months when stored at 0-5 °C.					
	 Preparation of working solution (1) Add the Dye Mixture stock solution to a conical tube and dilute with Assay Buffer. (2) Add the Enzyme stock solution to the solution prepared in step (1). ※ Refer to Table 2. ※ The working solution is light-sensitive. Prepare it immediately prior to use and protect it from light by covering the tube with aluminum foil. Prepare working solution fresh each day. 					
	F	action outer duy.	for 24 well	for 48 well	for 96 well	
	l –	Due Mixture steek colution				
	-	Dye Mixture stock solution	150 µL	300 µL	600 µL	
	-	Assay Buffer	1.35 mL	2.7 mL	5.4 mL	
	L	Enzyme stock solution	27 µL	54 µL	108 µL	
		Table 2. Examples	s of working	solution p	reparation	

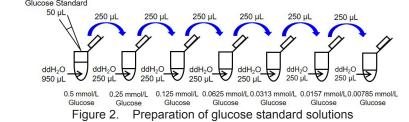
General Protocol 1. Sample preparation

- Prepare cell culture supernatants (samples).
- % Prepare samples at different dilutions to determine the most suitable concentration. Glucose concentrations should range from 0 to 0.5 mmol/L. Use double-deionized H₂O (ddH₂O) for dilutions.
- % Sample volume in the assay is 50 μL per well.

2. Preparation of glucose standard solution

Mix 50 μ L Glucose Standard (10 mmol/L) and 950 μ L ddH₂O in a microtube to prepare a 0.5 mmol/L glucose standard solution. Serially dilute with ddH₂O to prepare 0.5, 0.25, 0.125, 0.0625, 0.0313, 0.0157, 0.00785, and 0 mmol/L (Figure 2).

* To measure intracellular glucose concentrations, prepare glucose standard solutions in 0.1% Triton X-100 instead of ddH₂O. 10 mmol/L Glucose Standard



3. Measurements

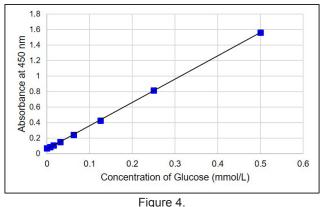
- (1) Add 50 µL glucose standard and sample solutions to separate wells in the 96-well microplate (Figure 3).
- * To obtain accurate data, triplicate samples are recommended.
- (2) Add 50 µL working solution to each well.
- * Because the enzymatic reaction starts immediately after working solution is added, it should be dispensed with a multichannel pipette to minimize time lags in pipetting.
- (3) Incubate the microplate at 37 °C for 30 minutes.
- ※ Seal the microplate during incubation to prevent evaporation.

(4) Measure absorbance at 450 nm with a microplate reader.

(5) Calculate the glucose concentration in each sample from a calibration curve generated from the glucose stan dards.

% Multiply the results by the dilution factor to determine glucose concentrations in the original samoles.

	1 2 3	4 5 6				
Α	0 mmol/L Glucose	Sample 1				
в	0.00785 mmol/L Glucose	Sample 2				
С	0.0157 mmol/L Glucose	Sample 3				
D	0.0313 mmol/L Glucose	Sample 4				
Е	0.0625 mmol/L Glucose	Sample 5				
F	0.125 mmol/L Glucose	Sample 6				
G	0.25 mmol/L Glucose	Sample 7				
н	0.5 mmol/L Glucose	Sample 8				



Typical calibration curve generated with glucose standards

Figure 3. Example of an assay plate format (n=3)

Experimental Inhibition of glucose uptake by phloretin

Example

- (1) Jurkat cells (5×10⁵ cells/mL, in RPMI containing 10% fetal bovine serum and 1% penicillin-streptomycin) were suspended in medium containing phloretin.
- (2) The cells prepared in step (1) were seeded in a 6-well microplate (1×10⁶ cells/well) and cultured overnight at 37 °C in a 37 °C incubator equilibrated with 95% air, 5% CO₂.
- (3) After the incubation, the cell suspension was transferred to a conical tube and centrifuged at 300×g for 5 minutes.
- (4) The supernatant (100 μL) was transferred to a 1.5 mL microtube and diluted 30-fold with ddH₂O (sample solution). Sample or glucose standard solutions, -50 μL each, were added to individual wells of 96-well micro plate.
- (5) Working solution (50 µL) was added to each well.
- (6) The assay plate was incubated at 37 °C for 30 minutes.
- (7) Absorbance at 450 nm was measured with a microplate reader and the concentrations of glucose in the samples were calculated from the calibration curve.

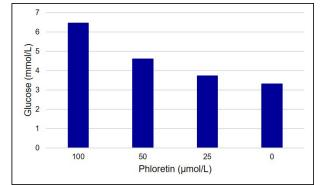


Figure 5. Inhibition of glucose uptake by phloretin Decreased glucose consumption in accordance with the concentration of phloretin, a glucose transport inhibitor, was confirmed.

Reference 1) Xiaoping, Z. et al., J. Biol. Chem., 2018, 293, 6623-6634.

For research use only. Not for use in diagnostic procedures. If you need more information, please contact Dojindo technical service.

Dojindo Laboratories 2025-5 Tabaru, Mashiki-machi, Kamimashi

2025-5 Tabaru, Mashiki-machi, Kamimashiki-gun, Kumamoto 861-2202, Japan Phone: +81-96-286-1515 Fax: +81-96-286-1525 E-mail: Info@dbjirdo.co.jp Web: www.dbjirdo.co.jp Dojindo Molecular Technologies,Inc. Tel: +1-301-987-2667 Web:http://www.dojindo.com/ Dojindo EU GmbH Tel: +49-89-3540-4805 Web: http://www.dojindo.eu.com/ Dojindo China Co., Ltd Tel: +86-21-6427-2302 Web:http://www.dojindo.cn/