

Glucose Uptake Colorimetric Assay Kit

rev.07/20

(Catalog # K676-100; 100 assays; Store at -20°C)

I. Introduction:

Glucose uptake is an important biological tool for studying cell signaling and glucose metabolism. Of many different methods available for measuring glucose uptake, 2-deoxyglucose (2-DG) has been most widely used because of its structural similarity to glucose. As with glucose, 2-DG can be taken up by glucose transporters and metabolized to 2-DG-6-phosphate (2-DG6P). 2-DG6P, however, cannot be further metabolized, and thus accumulates in the cells. The accumulated 2-DG6P is directly proportional to 2-DG (or glucose) uptake by cells. In BioVision's Glucose Uptake Colorimetric Assay Kit, 2-DG6P is oxidized to generate NADPH, which is determined by an enzymatic recycling amplification reaction. This easy to use non-radioactive kit is highly sensitive and can detect glucose uptake as low as 10 pmol/sample.

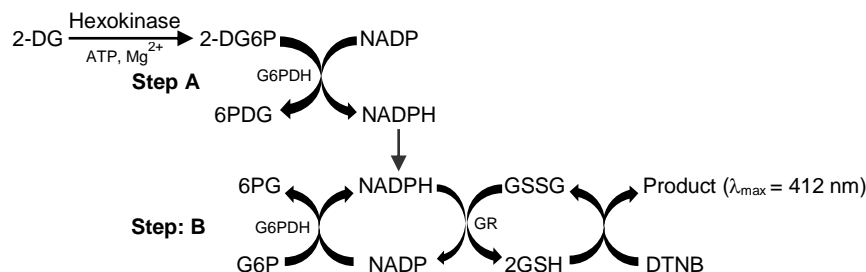


Figure 1. Assay Procedure. Step A: 2-DG oxidation to generate NADPH; Step B: NADPH recycling amplification reaction

II. Applications:

- Measurement of glucose uptake in response to insulin, growth factors, cytokines, mitogens, and nutrients, etc.
- Analysis of glucose metabolism and cell signaling in various cell types.
- Screening anti-diabetic drugs.

III. Sample Type:

- Adherent or suspension cells cultured in 96-well microtiter plate.

IV. Kit Contents:

Components	K676-100	Cap Code	Part Number
Extraction Buffer	17 ml	NM	K676-100-1
Neutralization Buffer	2.5 ml	Clear	K676-100-2
2-Deoxyglucose (2-DG, 10 mM)	1 ml	Purple	K676-100-3
Assay Buffer	25 ml	WM	K676-100-4
Enzyme Mix (Lyophilized)	1 vial	Orange	K676-100-5
Recycling Mix (Lyophilized)	1 vial	Blue	K676-100-6
2-DG6P Standard (Lyophilized)	1 vial	Yellow	K676-100-7
Glutathione Reductase	2 X 25 µl	Green	K676-100-8
Substrate-DTNB (Lyophilized)	2 vials	Red	K676-100-9

V. User Supplied Reagents and Equipment:

- 96-well clear plate with flat bottom
- Plate sealing tape
- Multi-well spectrophotometer (ELISA reader)
- Multi-channel pipette
- KRPH Buffer: 20 mM HEPES, 5 mM KH₂PO₄, 1 mM MgSO₄, 1 mM CaCl₂, 136 mM NaCl, 4.7 mM KCl, pH 7.4.

VI. Storage and Handling:

Store kit at -20°C, protected from light. Warm all Buffers to room temperature (RT) before use. Briefly centrifuge all small vials prior to opening.

VII. Reagent Preparation and Storage Conditions:

- **Enzyme Mix (Lyophilized):** Reconstitute with 220 µl Assay Buffer. Pipette up and down to dissolve completely. Divide into aliquots and store at -20°C. Avoid repeated freeze/thaw cycles.
- **Recycling Mix (Lyophilized):** Reconstitute with 220 µl Assay Buffer. Pipette up and down to dissolve completely. Divide into aliquots and store at -20°C. Avoid repeated freeze/thaw cycles.
- **2-DG6P Standard (Lyophilized):** Reconstitute the vial with 100 µl dH₂O to generate a 10 mM (10 nmol/µl) 2-DG6P Standard solution. Keep on ice while in use. Store at -20°C. Use within two months.
- **Glutathione Reductase:** Add 1.1 ml Assay Buffer. Mix well, aliquot and store at -20°C. Avoid repeated freeze/thaw cycles.
- **Substrate-DTNB (Lyophilized):** Reconstitute each vial with 1 ml Assay Buffer. Dissolve completely by pipetting up and down. Store at -20°C. Reconstituted substrate is stable for 2 months.

VIII. Glucose Uptake Assay Protocol:

1. **2-DG6P Standard Curve:** Dilute 2-DG6P Standard to 0.1 mM (100 pmol/µl) by adding 10 µl of 10 mM 2-DG6P to 990 µl Assay Buffer and mix well. Dilute further to 0.01 mM (10 pmol/µl) by adding 50 µl of 0.1 mM 2-DG6P to 450 µl Assay Buffer. Add 0, 2, 4, 6, 8 and 10 µl into a series of well on a 96 well plate in duplicate to generate 0, 20, 40, 60, 80 and 100 pmol/well of 2-DG6P Standard. Adjust volume to 50 µl/well with Assay Buffer. **Note:** Make fresh dilution with the 10mM 2-DG6P standard stock solution each time.

2. **Sample Preparation:** Treat cells with desired method. Example: For 3T3-L1 adipocytes, cells were seeded at a density of ~1500 cells per well in a 96-well plate, differentiated, then maintained for another 4 days prior to use. To assay glucose uptake, adipocytes were washed twice with PBS and starved in 100 μ l serum free adipocyte medium overnight (to increase glucose uptake), then rewashed 3X with PBS. The cells were starved for glucose by preincubating with 100 μ l Krebs-Ringer-Phosphate-HEPES (KRPH) buffer containing 2 % BSA for 40 min. then stimulated or not with 1 μ M insulin for 20 min to activate the glucose transporter. 10 μ l of 10 mM 2-DG was added and the cells incubated for 20 min. Wash cells 3X with PBS to remove exogenous 2-DG. To degrade endogenous NAD(P) and to denature enzymes, lyse cells with 80 μ l of Extraction Buffer, freeze/thaw once and heat at 85°C for 40 min. Cool the cell lysate on ice for 5 min and neutralize by adding 10 μ l of Neutralization Buffer. Briefly spin & dilute 1:10 times by adding 45 μ l of Assay Buffer to 5 μ l sample. Add 1-50 μ l sample per well. Adjust final volume to 50 μ l with Assay Buffer.

Notes:

1) As a control, prepare a parallel sample well not treated with insulin and 2-DG.

2) For other cell types, optimal incubation times and treatment protocols may vary from these conditions. We suggest testing several doses of cell lysate to ensure the readings are within the standard curve range. We suggest starving cells overnight in serum free medium to increase 2-DG uptake.

3. Reaction:

a. **NADPH Generation:** Mix enough reagents for the number of assays (Control, samples and Standards) to be performed. For each well, prepare 10 μ l Reaction Mix A:

<u>Reaction Mix A</u>	
Assay Buffer	8 μ l
Enzyme Mix	2 μ l

Add 10 μ l Reaction Mix A into each well. Mix and incubate at 37°C for 1 hr.

b. **NADP Degradation:** To degrade unused NADP (Figure 1, step A), add 90 μ l of Extraction Buffer to each well, seal with aluminum sealing tape and heat at 90°C for 40 min. Cool on ice for 5 min and add 12 μ l of Neutralization Buffer.

c. **Recycling Amplification Reaction:** For each well (control, Standards and samples), prepare 38 μ l recycling Reaction Mix B:

<u>Reaction Mix B</u>	
Glutathione Reductase	20 μ l
Substrate (DTNB)	16 μ l
Recycling Mix	2 μ l

Mix. Add 38 μ l Reaction Mix B into each well. Mix well.

4. **Measurement:** Measure absorbance at 412 nm in microplate reader at 37°C every 5 min until the 100 pmol standard reaches 1.5-2.0 OD. Take an endpoint reading of all samples and Standards.
5. **Calculation:** Subtract the 0 pmol Standard from all Standard readings. Plot the 2-DG6P Standard Curve. Correct the sample background by subtracting the value derived from the untreated cells (Control i.e. not treated with insulin and 2-DG) from all sample readings (Note: The background reading can be significant and must be subtracted from all sample readings). 2-DG concentration of the test samples, which is proportional to accumulated 2-DG6P can then, be calculated.

$$2\text{-DG uptake} = \frac{Sa}{Sv} \text{ (pmol/}\mu\text{l or nmol/ml or } \mu\text{M)}$$

Where: **Sa** is the amount of 2-DG6P (in pmol) in sample well calculated from the Standard Curve.

Sv is sample volume (in μ l) added into the sample well.

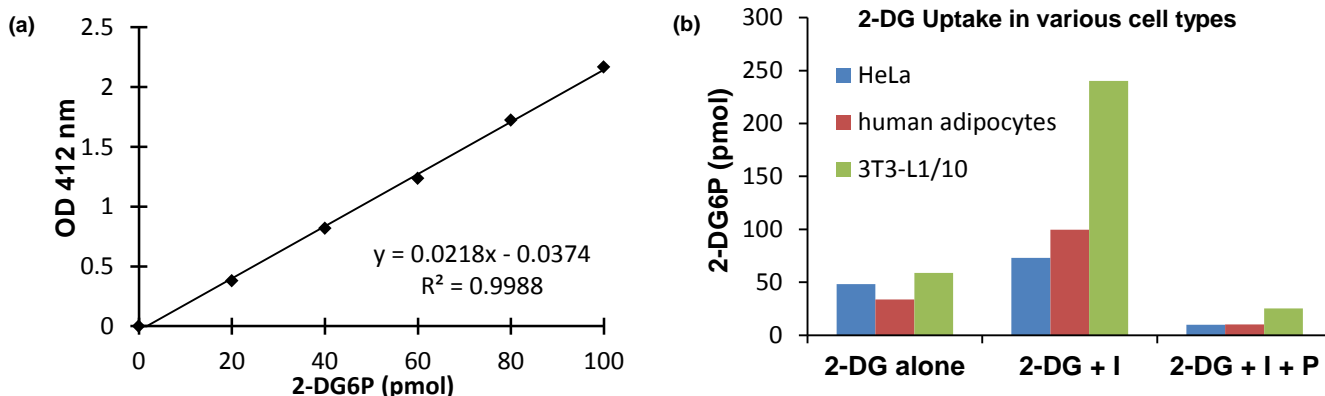


Figure 1. (a) 2-DG6P Standard Curve **(b)** 2-DG uptake in 3T3-L1, human adipocyte and HeLa cells. To scale on the same graph, data from 3T3-L1 cells is plotted at 10% of true value. Assays were performed following kit protocol. 2-DG = 2-deoxyglucose, I = Insulin; P = Phloretin.

IX. Related Products:

Glucose uptake Fluorometric Assay Kit

Glucose and Sucrose Assay Kit

Glucose Assay kit

NAD/NADH Quantification Kit

NADP/NADPH Quantification Kit

Glucose Dehydrogenase Activity Assay Kit

Glucose-6-Phosphate Dehydrogenase Assay Kit

Maltose and Glucose Assay Kit

PicoProbe™ Glucose-6-Phosphate Assay Kit

Adipogenesis Assay Kit

FOR RESEARCH USE ONLY! Not to be used on humans.