

## 2-NBDG Glucose Uptake Assay Kit (Cell-Based)

8/15

(Catalog # K682-50; 50 assays; Store at -20°C)

### I. Introduction:

Glucose is a ubiquitous energy source in most organisms and plays a pivotal role in cellular metabolisms and homeostasis. Cancer cells exhibit increased glucose uptake to support their high proliferation rate. 2-NBDG (2-deoxy-2-[(7-nitro-2,1,3-benzoxadiazol-4-yl) amino]-D-glucose) is a fluorescent deoxyglucose analog that can be taken up by cells through glucose transporters. However, 2-NBDG cannot be fully utilized in glycolysis because of its modification and thus accumulates inside the cells. Fluorescence generated by this fluorescent glucose analog is proportional to glucose uptake by the cells and can be used to measure glucose uptake using fluorescent microscopy and flow cytometry. To validate the assay, the kit includes phloretin, a natural phenol that inhibits glucose uptake. This easy to use non-radioactive kit allows imaging and accurate measurement of glucose uptake in cultured cells in response to insulin, growth factors etc.

### 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 compounds

### III. Sample Type:

- Adherent or suspension cells

### IV. Kit Contents:

Components	K682-50	Cap Code	Part Number
Analysis Buffer (50X)	1.8 ml	Brown	K682-50-1
2-NBDG Reagent (100X)	200 µl	Red	K682-50-2
Glucose Uptake Enhancer	1 ml	Blue	K682-50-3
Phloretin (100X)	75 µl	Yellow	K682-50-4

### V. User Supplied Reagents and Equipment:

- Cell culture medium
- PBS
- 24 or 12-well tissue culture plate
- Fluorescence microscope
- Flow cytometer with excitation filter at 488 nm wavelength

### VI. Storage Conditions and Reagent Preparation:

Store kit at -20°C, protected from light. Briefly centrifuge small vials prior to opening. Read entire protocol before performing the assay.

- **Analysis Buffer (50X):** Dilute Analysis Buffer with 1X PBS (not supplied) to make 1X Analysis Buffer. Keep on ice while in use.
- **2-NBDG Reagent (100X):** Aliquot and store at -20°C. For consistent results, avoid repeated freeze/thaw.

### VII. Glucose Uptake Assay Protocol:

The protocol described below is for 24-well tissue culture plate. Reagents, buffer, and the number of cells to be seeded should be adjusted for different size culture plates.

**1. Sample Preparation:** Seed adherent cells ( $2-5 \times 10^4$  cells/well) one day before starting the assay. After 8-12 hrs, remove regular culture medium (10% FBS) and treat cells with test compound or vehicle control in 400 µl cell culture medium with 0.5% FBS. Incubate cells at 37°C with 5% CO<sub>2</sub> for 1 hr or desired time period depending upon the test compound. To use Phloretin as a control, treat cells with 4 µl Phloretin (final concentration 1X) in 400 µl of cell culture medium with 0.5% FBS at 37°C with 5% CO<sub>2</sub> for 1 hr.

**Note:** Cell seeding is not required for suspension cells. Use up to  $1-2 \times 10^5$  suspension cells/well in 400 µl of cell culture medium with 0.5% FBS to treat with test compound or vehicle control.

**2. Glucose Uptake:** Prepare 400 µl glucose uptake mix for each well as following:

Cell culture medium (0.5% FBS)	376 µl
2-NBDG Reagent	4 µl
Glucose Uptake Enhancer	20 µl
Test compound or 1X Phloretin	same conc. as in step 1

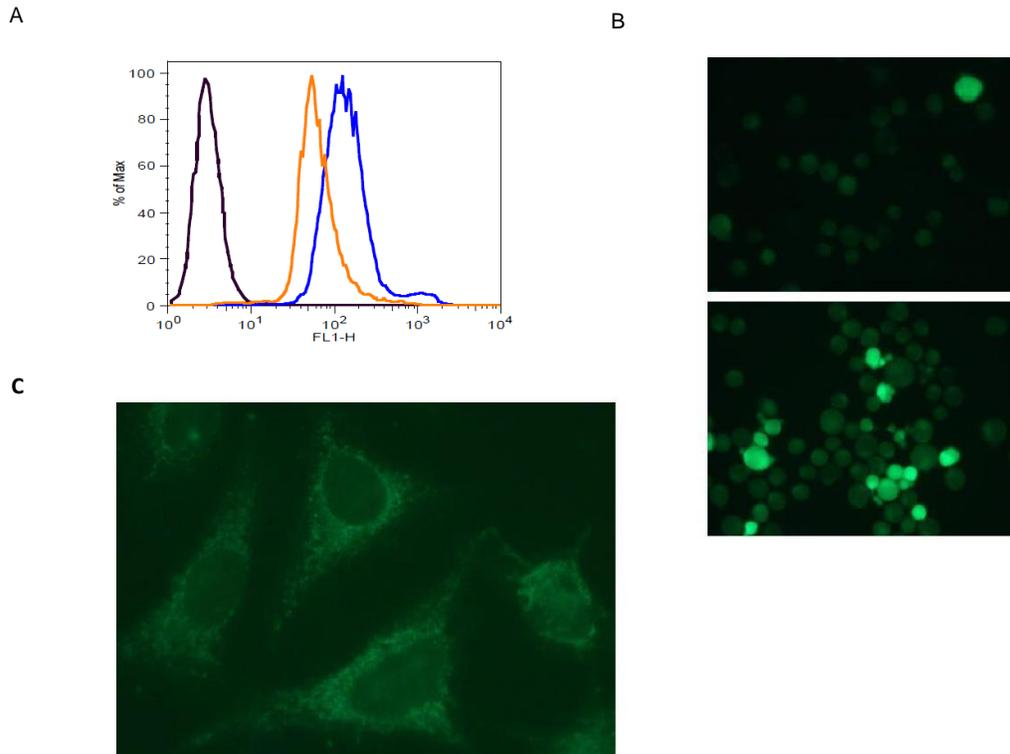
Mix well. After the incubation in Step 1, spin down the plate at 400 x g for 5 min. and carefully remove the medium without disturbing cells. Gently add glucose uptake mix to each well and incubate cells at 37°C with 5% CO<sub>2</sub> for 30 min.

**3. Measurement:** After incubation, collect cells from the plate and keep on ice. Wash once with 1 ml ice-cold 1X Analysis Buffer. Spin down at 400 x g for 5 min. and resuspend cell pellet in 400 µl of 1X Analysis Buffer. Cells are ready to be analyzed on flow cytometer (488 nm excitation laser). For flow acquisition and analysis, select the main cell population in the FSC vs SSC plot to exclude dead cells and cellular debris. Within the main cell population, mean fluorescence intensity in FL1 can be quantified and compared between cells treated with test compounds and untreated control cells.

#### Notes:

- Trypsin can be used to collect the adherent cells for performing this assay.

- b. The assay can be used to measure and compare glucose uptake levels in various cell types.
- c. Optional: To visualize the level of glucose uptake under fluorescence microscope, centrifuge plate at 400 x g for 5 min. Wash cells once with 500  $\mu$ l ice-cold 1X Analysis Buffer, and replace with fresh 200  $\mu$ l of 1X Analysis Buffer. Observe cells under fluorescence microscope using blue excitation fluorescence filter (excitation range 420 nm-495 nm).



**Figure: Glucose uptake in Jurkat and HeLa cells.**  $2.5 \times 10^5$  Jurkat cells were pretreated with or without 4  $\mu$ l phloretin (1X concentration) for 45 min. After pretreatment, cells were washed and incubated with 2-NBDG Reagent, Glucose Uptake Enhancer, and the same concentration of phloretin for another 30 min. according to kit's protocol. (A) Comparison of histograms from flow analysis showing the inhibition of glucose uptake by phloretin in Jurkat cells (Black: negative control cells; orange: in the presence of phloretin; blue: without phloretin). (B) Images of Jurkat cells obtained using fluorescence microscope (Top: treated with phloretin; Bottom: without phloretin treatment). (C) Glucose Uptake in HeLa cells: HeLa cells showing the uptake of 2-NBDG Reagent in the cytoplasm. Cells were stained with 2-NBDG Reagent for 30 min. and fixed. Image was taken using a fluorescent microscope with a 60X objective lens.

#### VIII. RELATED PRODUCTS:

Glucose Uptake Colorimetric Assay Kit (K676)	Glucose Uptake Fluorometric Assay Kit (K666)
GluTracker™ Glucose Uptake Assay Kit (Cell-Based) (K681)	Glucose-6-Phosphate Colorimetric Assay Kit (K657)
Glucose Colorimetric Assay Kit II (K686)	Glucose Colorimetric/Fluorometric Assay Kit (K606)
Glucose and Sucrose Colorimetric/Fluorometric Assay Kit (K616)	Maltose and Glucose Colorimetric/Fluorometric Assay Kit (K618)
PicoProbe™ Glucose-6-Phosphate Fluorometric Assay Kit (K687)	PicoProbe™ Glucose Fluorometric Assay Kit (K688)
Glucose Dehydrogenase Activity Colorimetric Assay Kit (K786)	Glucose Oxidase Activity Colorimetric/Fluorometric Assay Kit (K788)
Glucose-1-Phosphate (G1P) Colorimetric Assay Kit (K697)	GluTracker™ Glucose Uptake Assay Kit (Cell-Based) (K681)
2-NBDG (2760)	Glucose AGE-BSA-II (2223)
G6PD, Human Recombinant (7532)	SGLT-2 Antibody (3690)
GPI, human recombinant (7822)	GLUT4 Antibody (3945)
SGLT-1 Antibody (3711)	

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