

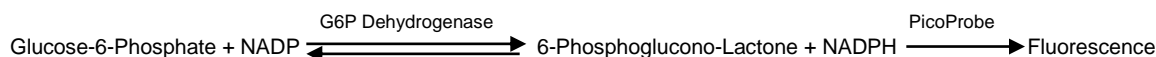
## PicoProbe™ Glucose-6-Phosphate Dehydrogenase Activity Assay Kit (Fluorometric)

(Catalog # K751-100; 100 assays; Store at -20°C)

7/18

### I. Introduction:

Glucose-6-Phosphate Dehydrogenase (G6PDH: EC 1.1.1.49) is a cytosolic enzyme in the pentose phosphate pathway, a metabolic pathway that supplies reducing energy to cells (such as erythrocytes) by maintaining the level of the co-enzyme nicotinamide adenine dinucleotide phosphate (NADPH). The NADPH in turn maintains the level of glutathione in these cells that helps protect the red blood cells against oxidative damage. Of greater quantitative importance is the production of NADPH for tissues actively engaged in biosynthesis of fatty acids and/or isoprenoids, such as liver, mammary gland, adipose tissue, and adrenal gland. BioVision's PicoProbe™ Glucose-6-Phosphate Dehydrogenase Assay kit provides a quick and easy method for monitoring G6PDH activity in a wide variety of samples. In this assay, G6PDH converts G6P into pyruvate and NADPH, which further reduces PicoProbe™ to generate an intense fluorescence product (Ex/Em = 535/587 nm). This kit is simple, sensitive and high-throughput adaptable and can detect as low as 1  $\mu$ U of G6PDH activity.



### II. Applications:

- Measurement of G6PDH activity in various tissues and cells
- Evaluation of pentose phosphate pathway

### III. Sample Type:

- Animal tissues: muscle, liver, heart, kidney, etc.
- Cell culture: adherent or suspension cells
- Plant tissues

### IV. Kit Contents:

| Components                | K751-100 | Cap Code | Part Number |
|---------------------------|----------|----------|-------------|
| G6PDH Assay Buffer        | 25 ml    | WM       | K751-100-1  |
| PicoProbe™ (in DMSO)      | 0.4 ml   | Blue     | K751-100-2  |
| G6PDH Substrate           | 1 Vial   | Orange   | K751-100-3  |
| G6PDH Developer           | 1 Vial   | Red      | K751-100-4  |
| G6PDH Positive Control    | 1 Vial   | Green    | K751-100-5  |
| NADPH Standard (200 nmol) | 1 Vial   | Yellow   | K751-100-6  |

### V. User Supplied Reagents and Equipment:

- 96-well white plate with flat bottom
- Multi-well spectrophotometer (ELISA reader)

### 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.

- **G6PDH Assay Buffer:** Bring to room temperature before use. Store at 4°C or -20°C.
- **PicoProbe™:** Before use, thaw at room temperature. Store at -20°C. Use within two months.
- **G6PDH Substrate and Developer:** Reconstitute with 220  $\mu$ l Assay Buffer. Pipette up and down to dissolve completely. Store at -20°C. Use within two months.
- **G6PDH Positive Control:** Reconstitute with 100  $\mu$ l Assay Buffer and mix thoroughly. Aliquot and store at -70°C. Avoid freeze/thaw. Use within two months. Keep on ice while in use.
- **NADPH Standard:** Reconstitute with 200  $\mu$ l dH<sub>2</sub>O to generate 1 mM (1 nmol/ $\mu$ l) NADPH Standard solution. Aliquot and store at -20°C. Use within two months. Keep on ice while in use.

### VII. G6PDH Activity Assay Protocol:

1. **Sample Preparation:** Homogenize tissue (~10 mg) or cells ( $1 \times 10^6$ ) with 100  $\mu$ l ice cold G6PDH Assay Buffer. Keep on ice for 10 min. Centrifuge at 10,000 X g, 4°C for 5 min. and collect supernatant. Dilute the supernatant ~10 fold in Assay Buffer and add 1-50  $\mu$ l into desired well(s) in a 96-well plate. For Positive Control, dilute G6PDH Positive Control 200 times with G6PDH Assay Buffer just before use and add 2-20  $\mu$ l of diluted G6PDH Positive Control into desired well(s). Adjust the volume of Positive Control and sample wells to 50  $\mu$ l/well with G6PDH Assay Buffer.

#### Notes:

- a. For unknown samples, we suggest doing pilot experiment and testing several amounts of G6PDH to ensure the readings are within the Standard Curve range.
  - b. If sample has high background, prepare parallel sample well(s) as sample background control.
  - c. Don't store the diluted G6PDH Positive Control.
2. **NADPH Standard Curve:** Dilute NADPH Standard to 40  $\mu$ M (40 pmol/ $\mu$ l) by adding 40  $\mu$ l of 1 mM NADPH Standard to 960  $\mu$ l of dH<sub>2</sub>O. Add 0, 2, 4, 6, 8, and 10  $\mu$ l of 40  $\mu$ M NADH Standard into a series of wells in a 96-well plate to generate 0, 80, 160, 240, 320 and 400 pmol/well of NADPH Standard. Adjust the volume to 50  $\mu$ l/well with G6PDH Assay Buffer.
  3. **Reaction Mix:** Mix enough reagents for the number of assays to be performed. For each well, prepare 50  $\mu$ l Mix containing

|                    | Reaction Mix | *Background Control Mix |
|--------------------|--------------|-------------------------|
| G6PDH Assay Buffer | 44 $\mu$ l   | 46 $\mu$ l              |
| PicoProbe™         | 2 $\mu$ l    | 2 $\mu$ l               |
| G6PDH Developer    | 2 $\mu$ l    | 2 $\mu$ l               |
| G6PDH Substrate    | 2 $\mu$ l    | ---                     |

Mix. Add 50  $\mu$ l of Reaction Mix to each well containing Standards, Positive Control, and samples. Mix well.

\* For samples having background, add 50  $\mu$ l of Background Control Mix to sample background control well(s).

**4. Measurement:** Measure fluorescence (Ex/Em = 535/587 nm) immediately in kinetic mode for 10-40 min. at 37°C.

**Note:** Incubation time depends on the G6PDH activity in the samples. We recommend measuring fluorescence in kinetic mode, and choosing two time points ( $T_1$  and  $T_2$ ) in the linear range to calculate the G6PDH activity of the samples. The NADPH Standard Curve can be read in endpoint mode (i.e. at the end of incubation time).

**5. Calculation:** Subtract 0 Standard reading from all readings. Plot the NADPH Standard curve. If sample background control reading is significant, subtract the sample background control reading from sample reading. Calculate the G6PDH activity of the test sample:  $\Delta RFU = RFU_2 - RFU_1$ . Apply  $\Delta RFU$  to NADPH Standard Curve to get B pmol of NADPH generated by G6PDH during the reaction time ( $\Delta T = T_2 - T_1$ ).

$$\text{Sample G6PDH Activity} = B / (\Delta T \times V) \times D = \text{pmol/min}/\mu\text{l} = \mu\text{U}/\mu\text{l} = \text{mU/ml}$$

Where: **B** is NADPH amount in the sample well from Standard Curve (pmol)

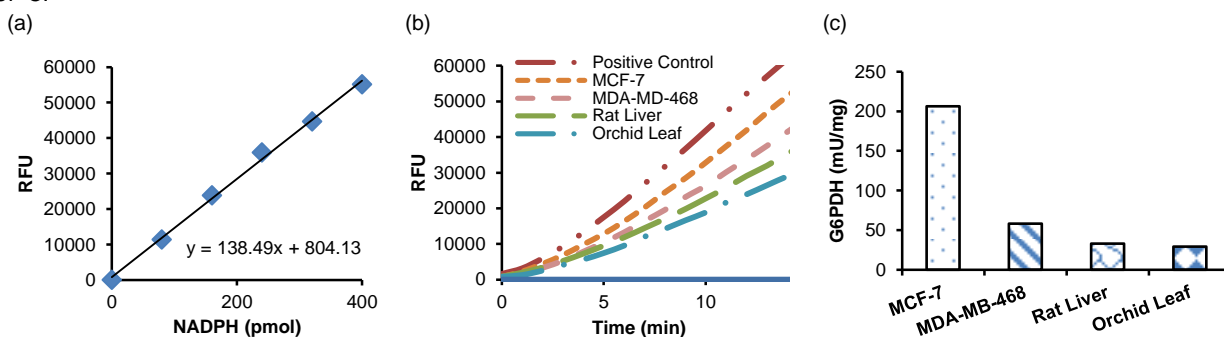
$\Delta T$  is reaction time (min.)

**V** is sample volume added into the reaction well ( $\mu$ l)

**D** is dilution factor

G6PDH Activity in samples can also be expressed in mU/mg of protein.

**Unit Definition:** One unit of G6P Dehydrogenase is the amount of enzyme that generates 1.0  $\mu$ mol of NADPH per min. at pH 8.0 at 37°C.



**Figure:** (a) NADPH Standard Curve. (b) Kinetic measurement of G6PDH activity in various samples. (c) G6PDH specific activity was calculated in lysates prepared from MCF-7 (0.29  $\mu$ g), MDA-MB-468 (0.41  $\mu$ g), rat liver (0.6  $\mu$ g) and orchid leaf (0.56  $\mu$ g). Assays were performed following the kit protocol.

**VIII. RELATED PRODUCTS:**

Glucose-6-Phosphate Dehydrogenase Activity Assay Kit (K757)  
 Pyruvate Colorimetric/Fluorometric Assay Kit (K609)  
 Triose Phosphate Isomerase Assay Kit (K670)  
 Lactate Colorimetric Assay Kit II (K627)  
 Phosphoglucumutase Assay Kit (K774)  
 PicoProbe™ D-Lactate Fluorometric Assay Kit (K668)  
 Glucose-6-Phosphate Assay Kit (K657)

Hexokinase Colorimetric Assay Kit (K789)  
 Pyruvate Dehydrogenase Activity Assay Kit (K679)  
 PicoProbe™ NADH Assay Kit (K338)  
 PicoProbe™ NADPH Assay Kit (K349)  
 Phosphoglucose Isomerase Assay Kit (K775)  
 Glucose-1-Phosphate Assay Kit (K697)  
 Glucose Dehydrogenase Activity Assay Kit (K786)

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