

## QuantiChrom™ Glucose Dehydrogenase Kit (DGDH-100)

### Quantitative Colorimetric Kinetic Glucose Dehydrogenase Activity Determination

#### DESCRIPTION

GLUCOSE DEHYDROGENASE (GDH) belongs to the family of oxidoreductases, specifically those acting on the CH-OH group of donor with other acceptors. GDH participates in the pentose phosphate pathway. BioAssay Systems' non-radioactive, colorimetric GDH assay is based on the reduction of the tetrazolium salt MTT in a NADH-coupled enzymatic reaction to a reduced form of MTT which exhibits an absorption maximum at 565 nm. The increase in absorbance at 565 nm is proportional to the enzyme activity.

#### KEY FEATURES

**Fast and sensitive.** Linear detection range (20 µL sample): 0.5 to 200 U/L for 15 min reaction.

**Convenient and high-throughput.** Homogeneous "mix-incubate-measure" type assay. Can be readily automated on HTS liquid handling systems for processing thousands of samples per day.

#### APPLICATIONS

GDH activity determination in biological samples (e.g. plasma, serum, tissue and culture media).

#### KIT CONTENTS (100 TESTS IN 96-WELL PLATES)

<b>Assay Buffer:</b>	10 mL	<b>Diaphorase:</b>	120 µL
<b>NAD/MTT:</b>	1 mL	<b>Calibrator:</b>	1.5 mL
<b>Substrate:</b>	1 mL		

**Storage conditions.** The kit is shipped at ambient temperature. Store all components at -20°C upon receiving. Shelf life: 6 months after receipt.

**Precautions:** reagents are for research use only. Normal precautions for laboratory reagents should be exercised while using the reagents. Please refer to Material Safety Data Sheet for detailed information.

#### PROCEDURES

This assay is based on a kinetic reaction. To ensure identical incubation time, addition of Working Reagent to samples should be quick and mixing should be brief but thorough. Use of a multi-channel pipettor is recommended. Assays can be executed at any desired temperature (e.g. 25°C or 37°C).

**Sample Preparation:** Serum and plasma are assayed directly.

**Tissue:** prior to dissection, rinse tissue in phosphate buffered saline (pH 7.4) to remove blood. Homogenize tissue (50 mg) in ~200 µL buffer containing 50 mM potassium phosphate (pH 7.5). Centrifuge at 10,000 × g for 15 min at 4°C. Remove supernatant for assay.

**Cell Lysate:** collect cells by centrifugation at 2,000 × g for 5 min at 4°C. For adherent cells, do not harvest cells using proteolytic enzymes; rather use a rubber policeman. Homogenize or sonicate cells in an appropriate volume of cold buffer containing 50 mM potassium phosphate (pH 7.5). Centrifuge at 10,000 × g for 15 min at 4°C. Remove supernatant for assay.

All samples can be stored at -20 to -80°C for at least one month.

**Reagent Preparation:** Equilibrate reagents to desired reaction temperature (e.g. 25°C or 37°C). Briefly centrifuge tubes before use.

Prepare enough Working Reagent (WR) for all assay wells by mixing, for each 96-well assay: 8 µL Substrate, 8 µL NAD/MTT Solution, 1 µL Diaphorase and 70 µL Assay Buffer.

#### Reaction Preparation:

1. Transfer 100 µL H<sub>2</sub>O (OD<sub>H2O</sub>) and 100 µL Calibrator (OD<sub>CAL</sub>) solution into wells of a clear flat bottom 96-well plate.
2. Transfer 20 µL of each sample into separate wells and then add 80 µL WR to each sample well. Tap plate briefly to mix.
3. Read OD<sub>565nm</sub> (OD<sub>0</sub>), and again after 15 min (OD<sub>15</sub>) on a plate reader.

#### CALCULATION

Subtract the OD<sub>0</sub> from OD<sub>15</sub> for each sample to compute the ΔOD<sub>S</sub> values. GDH activity can then be calculated as follows:

$$\text{GDH Activity} = \frac{\Delta\text{OD}_S}{\epsilon_{\text{mtt}} \cdot l} \times \frac{\text{Reaction Vol } (\mu\text{L})}{t \text{ (min)} \cdot \text{Sample Vol } (\mu\text{L})} \times n$$

$$= \frac{\Delta\text{OD}_S}{\text{OD}_{\text{CAL}} - \text{OD}_{\text{H}_2\text{O}}} \times \frac{273}{t \text{ (min)}} \times n \text{ (U/L)}$$

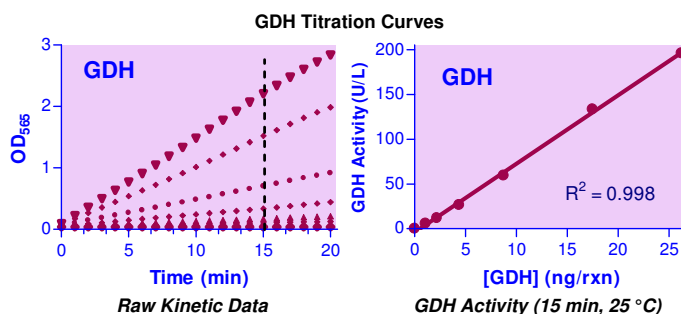
where  $\epsilon_{\text{mtt}}$  is the molar absorption coefficient of reduced MTT.  $l$  is the light pathlength which is calculated from the calibrator. OD<sub>CAL</sub> and OD<sub>H2O</sub> are OD<sub>565nm</sub> (OD<sub>0</sub>) values of the Calibrator and water.  $t$  is the reaction time (15 min is the recommended time). Reaction Vol and Sample Vol are 100 µL and 20 µL, respectively.  $n$  is the dilution factor.

**Unit definition:** 1 Unit (U) of GDH will catalyze the conversion of 1 µmole of NAD to NADH per min at pH 8.2.

**Note:** If sample GDH activity exceeds 200 U/L, either use a shorter reaction time or dilute samples in water and repeat the assay. For samples with GDH activity < 5 U/L, the incubation time can be extended up to 2 hours.

#### MATERIALS REQUIRED, BUT NOT PROVIDED

Pipetting devices and accessories (e.g. multi-channel pipettor), clear flat-bottom 96-well plates (e.g. VWR cat# 82050-760), centrifuge tubes and plate reader.



#### LITERATURE

1. Bak, TG (1967) "Studies on glucose dehydrogenase of *Aspergillus oryzae*. II Purification and physical and chemical properties". *Biochim. Biophys. Acta*. 139: 277-93.
2. Brink, NG, et al. (1953) "Beef liver glucose dehydrogenase. 1. Purification and properties". *Acta Chem. Scand.* 7: 1081-1089.
3. Thompson RE, Carper WR (1970) "Glucose dehydrogenase from pig liver. I. Isolation and purification". *Biochim. Biophys. Acta* 198 (3): 397-406.

