

QuantiChrom™ Glucose Assay Kit (DIGL-100)

Quantitative Colorimetric Glucose Determination at 630nm

DESCRIPTION

Glucose (C₆H₁₂O₆) is a ubiquitous fuel molecule in biology. It is oxidized through a series of enzyme-catalyzed reactions to form carbon dioxide and water, yielding the universal energy molecule ATP. Due to its importance in metabolism, glucose level is a key diagnostic parameter for many metabolic disorders. Increased glucose levels have been associated with diabetes mellitus, hyperactivity of thyroid, pituitary and adrenal glands. Decreased levels are found in insulin secreting tumors, myxedema, hypopituitarism and hypoadrenalism.

Simple, direct and automation-ready procedures for measuring glucose concentrations find wide applications in research and drug discovery. BioAssay Systems' glucose assay kit is designed to measure glucose directly in serum or plasma without any pretreatment. The improved o-toluidine method utilizes a specific color reaction with glucose. The absorbance at 630nm is directly proportional to glucose concentration in the sample.

KEY FEATURES

Sensitive and accurate. Use as little as 5 µL samples. Linear detection range 0.7 mg/dL (39 µM) to 300 mg/dL (16.6 mM) glucose in 96-well plate.

Simple and convenient. The procedure involves addition of a single working reagent and incubation for 8 min in a boiling water bath.

Improved reagent stability. The optimized formulation has greatly enhanced the reagent and signal stability.

Low interference in biological samples. No pretreatments are needed. Assays can be directly performed on serum and plasma samples.

APPLICATIONS:

Direct Assays: glucose in biological samples (e.g. serum and plasma).

Drug Discovery/Pharmacology: effects of drugs on glucose metabolism.

Food and Beverages: glucose in food, beverages etc.

KIT CONTENTS

Catalog #	Tests (96-well)	Reagent	Standard
DIGL-100	100	50 mL	1 mL 300 mg/dL

Storage conditions. The kit is shipped at room temperature. Store the reagent at room temperature and standard at -20°C, respectively. Shelf life: 12 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

IMPORTANT: THE REAGENT CONTAINS ACETIC ACID. THIS ASSAY IS PREFERABLY CARRIED OUT IN A CHEMICAL FUME HOOD.

Procedure using 96-well plate:

1. Dilute standard in distilled water as follows.

No	STD + H ₂ O	Vol (µL)	Glucose (mg/dL)
1	150µL + 0µL	150	300
2	100µL + 50µL	150	200
3	50µL + 100µL	150	100
4	25µL + 125µL	150	50
5	0µL + 150µL	150	0

Set up 1.5-mL centrifuge tubes. Transfer 5 µL diluted standards and samples to appropriately labeled tubes. Transfer 500 µL Reagent to each tube. Close the tubes tightly and mix. Store diluted standards at -20°C for future use.

- Place the tubes in a tube holder and heat in a boiling water bath or heat block for 8 min. Cool down in cold water bath for 4 min.
- Transfer 200 µL in duplicate into a clear bottom 96-well plate. Careful: avoid bubble formation. Read optical density at 620-650nm (peak absorbance at 630nm).

Procedure using cuvette:

1. Dilute standards and transfer 12 µL water blank, Standards and

samples to appropriately labeled tubes. Transfer 1200 µL Reagent to each tube. Close the tubes tightly and mix.

2. Place the tubes in a tube holder and heat in a boiling water bath for 8 min. Cool down in cold- water bath for 4 min.

3. Transfer 1000 µL reaction mixture into cuvet. Read optical density at 620-650nm (peak absorbance at 630nm) against blank.

Note: 1. if the Sample OD is higher than the Standard OD at 300mg/dL, dilute sample in water and repeat the assay. 2. To determine low glucose concentrations, use 50 µL sample and standards (instead of 5 µL) per 500 µL Reagent.

CALCULATION

Subtract blank OD (water, #5) from the standard OD values and plot the OD against standard concentrations. Determine the slope using linear regression fitting. The glucose concentration of Sample is calculated as

$$= \frac{OD_{\text{SAMPLE}} - OD_{\text{BLANK}}}{\text{Slope}} \quad (\text{mg/dL})$$

OD_{SAMPLE} and OD_{BLANK} are optical density values of the sample and sample "Blank" (water or buffer in which the sample was diluted). Typical serum/plasma glucose values: 70 - 110 mg/dL.

Conversions: 1mg/dL glucose equals 55.5 µM, 0.001% or 10 ppm.

MATERIALS REQUIRED, BUT NOT PROVIDED

Pipeting devices, centrifuge tubes, boiling water bath, tube holder.

Procedure using 96-well plate:

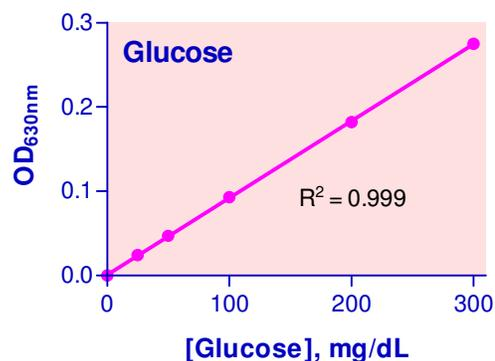
Clear bottom 96-well plates (e.g. Corning Costar) and plate reader.

Procedure using cuvette:

Spectrophotometer and Cuvets for measuring OD at 620-650nm.

EXAMPLES:

Rat plasma, rat serum, goat serum and human plasma were assayed using the 96-well plate assay protocol. The glucose concentrations were 128 ± 2 (n = 4), 72.5 ± 0.8 (n = 4), 78.6 ± 0.6, 69.3 ± 0.7 mg/dL (n = 4), respectively. Coefficient of variance < 3%.



Standard Curve in 96-well plate assay

PUBLICATIONS

1. Jeong, S., et al. (2021). High fructose drives the serine synthesis pathway in acute myeloid leukemic cells. *Cell Metabolism* 33(1): 145-159.e6.

2. Alghadir, A. H., et al. (2021). Moderate aerobic training modulates cytokines and cortisol profiles in older adults with cognitive abilities. *Cytokine* 138: 155373.

3. Hedbacker, K et al (2010). Antidiabetic effects of IGFBP2, a leptin-regulated gene. *Cell Metab.* 11(1):11-22.

