A convenient, easy-to-follow shortened protocol is now being provided with this assay. For a detailed protocol go to www.caymanchem.com/pdfs/10009582.pdf

## Glucose Colorimetric Assay Kit Short Protocol Item No. 10009582

#### **REAGENT PREPARATION**

- 1. Glucose Assay Standard (1,000 mg/dL; Item No. 10010098) This standard is ready to use as supplied.
- 2. Sodium Phosphate Assay Buffer (Item No. 700003) Dilute the contents with 40 ml of HPLC-grade water (final formulation is 50 mM sodium phosphate, pH 7.2); stable for three months at 4°C.
- **3.** Glucose Colorimetric Enzyme Mixture (Item No. 10010100) Reconstitute vial (1 vial sufficient for 60 wells) with 6 ml of diluted Assay Buffer and mix well; stable for one hour at 4°C.

## STANDARD PREPARATION

Dilute 50  $\mu$ l of the 1,000 mg/dL Glucose Standard with 450  $\mu$ l of diluted Assay Buffer to make a 100 mg/dL stock. Prepare Assay Standards as shown in Table 1. Standards are stable for 2 hrs at room temperature.

Tube	Glucose Stock (µl) (100 mg/dL)	Assay Buffer (µl)	Glucose Concentration (mg/dL)
А	0	200	0
В	5	195	2.5
С	10	190	5
D	15	185	7.5
E	20	180	10
F	30	170	15
G	40	160	20
Н	50	150	25



Short Protocol Item 10009582

Tech Support 888.526.5351 techserv@caymanchem.com

For Research Use Only www.caymanchem.com

## Glucose Colorimetric Assay Kit Short Protocol Item No. 10009582

#### PERFORMING THE ASSAY

- 1. Glucose Standard wells Add 85 µl of diluted Assay Buffer and 15 µl of each Standard (tubes A-H) to duplicate wells.
- 2. Sample wells Add 85 µl of diluted Assay Buffer and 15 µl of sample to duplicate wells.
- 3. Initiate the reaction Add 100 µl of Enzyme Mixture to all standard and sample wells.
- 4. Cover plate and Incubate for 10 minutes at 37°C.
- 5. Remove the plate cover and read the absorbance at 500-520 nm.

# CALCULATIONS

- 1. Calculate the average absorbance of each standard and sample.
- 2. Subtract the absorbance value of the standard A (0 mg/dL) from itself and all other values (both standards and samples). This is the corrected absorbance.
- 3. Plot the corrected absorbance values of each standard as a function of the concentration of glucose and fit the data to a linear regression.
- 4. Calculate the concentration of glucose for each sample from the standard curve using the equation below.

$$Glucose (mg/dL) = \left[\frac{(Corrected absorbance) - (y-intercept)}{Slope}\right] x dilution$$



www.caymanchem.com

03/19/2013