

## QuantiChrom™ β-Glucuronidase Assay Kit (DGCD-100)

### Quantitative Fluorimetric Kinetic β-Glucuronidase Activity Determination

#### DESCRIPTION

**BETA-GLUCURONIDASE (BG)** is an enzyme which catalyzes the hydrolysis of β-glucuronide bonds. In humans, β-glucuronidase catalyzes the hydrolysis of β-D-glucuronic acids from glycoproteins and mucopolysaccharides such as heparan sulfate. β-glucuronidase activity is important for drug metabolism as drugs are often conjugated to glucuronic acid to make for a more water soluble delivery system, relying on subsequent β-glucuronidase activity to release the active drug. Bilirubin clearance is also facilitated by glucuronidation. GUSB is often used as a reporter gene to track gene expression through β-glucuronidase activity. BioAssay Systems' fluorimetric BG assay uses a β-D-glucuronide substrate which fluoresces upon hydrolysis with β-glucuronidase and is read at 365/450 nm.

#### KEY FEATURES

**Highly Sensitive.** The fluorimetric nature of the kit detects  $1 \times 10^{-4}$  U/L to 8 U/L BG activity in 96 well plate with only 10 μL sample within 30 min.

**Versatile.** This assay has been validated to work with both animal and *E. coli* derived β-glucuronidase.

**High-throughput.** Homogeneous "mix-incubate-measure" type assay. Can be readily automated to assay thousands of samples per day.

#### APPLICATIONS

For quantitative determination of beta glucuronidase enzyme activity in tissue and bacterial samples.

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

**Substrate:** 4 mL                      **Standard:** 100 μL  
**Stop Reagent:** 1.5 mL

**Storage conditions.** The kit is shipped on ice. Store all kit components at -20 °C. Shelf life of 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 substrate to samples should be quick and mixing should be brief but thorough. Use of a multi-channel pipettor is recommended. Assays can be run at room temperature or 37°C.

**Sample Preparation:** we recommend to avoid surfactants such as SDS or Triton 100-X in sample preparation.

**Tissue:** Prior to dissection, rinse tissue in phosphate buffered saline (pH 7.4) to remove blood. Homogenize tissue (50 mg) in ~200 μL lysis buffer.

**Cell Lysate:** Collect cells by centrifugation at 2,000 x 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 lysis buffer. All samples can be stored at -20 to -80°C for at least one month.

**Reagent Preparation:** Equilibrate all kit reagents to room temperature. Invert or gently vortex tubes and bottles to ensure reagents are mixed.

**Standard Preparation:** Prepare 100 μM Premix by mixing 5 μL of the provided Standard with 995 μL dH<sub>2</sub>O. Prepare the standards as shown in the following table.

No	Premix + H <sub>2</sub> O	Vol (μL)	Standard (μM)
1	100μL + 0μL	100	100
2	60μL + 40μL	100	60
3	30μL + 70μL	100	30
4	0μL + 100μL	100	0

**Sample Preparation:** For tissue lysates we recommend you dilute samples 5-10 fold with deionized water prior to assay. For serum samples a 15-fold dilution may be required.

1. Transfer 10 μL Standards and 10 μL samples into wells of a black flat bottom 96 well plate.
2. Add 40 μL substrate to each well. Tap plate to mix. Incubate plate at RT or 37°C for 30 minutes.
3. Add 15 μL stop reagent to each well. Tap plate to mix. Incubate for 15 minutes.
4. Read Fluorescence at λ<sub>ex</sub>/em = 365/450nm.

This assay procedure can be readily miniaturized for 384- or 1536-wells.

#### CALCULATION

BG activity can then be calculated as follows:

$$\text{BG Activity} = \frac{F_{\text{SAMPLE}} - F_{\text{BLANK}}}{\text{Time} \cdot \text{Slope}} \times n \text{ (U/L)}$$

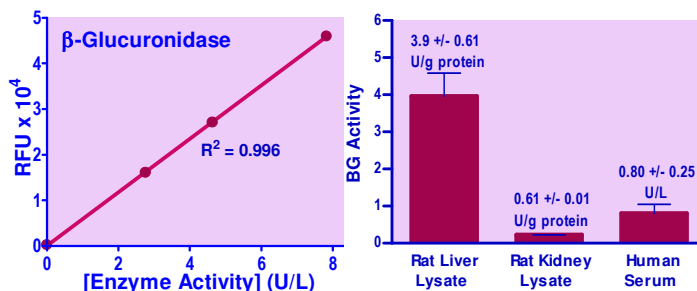
where F<sub>SAMPLE</sub> is the RFU value for each sample and F<sub>BLANK</sub> is the RFU value of the water (standard #4). Slope is the slope of the standard curve and Time is the incubation time (30 min). n is the sample dilution factor.

**Unit definition:** 1 Unit (U) of BG will catalyze the conversion of 1 μmole of the fluorescent glucuronide at 37°C and pH 5.0.

*Note: If sample exceeds 8 U/L calculated BG activity, either use a shorter incubation time or dilute samples in water and repeat the assay. For samples with calculated BG activity < 0.2 U/L, the incubation time can be extended up to 2 hours for greater sensitivity.*

#### MATERIALS REQUIRED, BUT NOT PROVIDED

Pipetting devices and accessories (e.g. multi-channel pipettor), black flat-bottom 96-well plates (e.g. VWR cat# 89089-582), centrifuge tubes and plate reader.



#### LITERATURE

1. Matsushita-Oikawa, H, et al. (2006) Novel carbohydrate-binding activity of bovine liver β-glucuronidase toward lactose/N-acetylglucosamine sequences. *Glycobiology*, 16(10), 891-901.
2. Marathe, S. V., & McEwen, J. E. (1995). Vectors with the gus reporter gene for identifying and quantitating promoter regions in *Saccharomyces cerevisiae*. *Gene*, 154(1), 105-107.

