

## QuantiChrom™ Glyoxalase I Assay Kit (DGLO-100)

### Colorimetric Determination of Glyoxalase I Activity

#### DESCRIPTION

**GLYOXALASE I (GLO-1)**, a lactoylglutathione lyase also known as methylglyoxalase, aldoketomutase, ketone-aldehyde mutase, and (R)-S-lactoylglutathione methylglyoxal-lyase, is an enzyme that catalyzes the isomerization of hemithioacetal adducts which are formed in spontaneous reactions between glutathionyl groups and aldehydes. The primary physiological function of glyoxalase I is the detoxification of methylglyoxal, a reactive 2-oxoaldehyde that is cytostatic at low concentrations and cytotoxic at millimolar concentrations. Glyoxalase I is a target for the development of pharmaceuticals against bacteria, protozoans and human cancer.

Simple, direct and automation-ready procedures for measuring GLO-1 activity in biological samples are highly desirable in research and drug discovery. BioAssay Systems' Glyoxalase I assay kit provides a sensitive and convenient method for GLO-1 activity determination. The method involves monitoring the increase in the product of the GLO-1 reaction, S-lactoylglutathione, by measuring the change in absorbance at 240 nm.

#### KEY FEATURES

**Sensitive and accurate.** Detection limit 4 U/L GLO-1 activity.

**Simple and high-throughput.** The procedure involves incubation of the provided substrate with the sample in a microplate. Can be readily automated as a high-throughput assay for thousands of samples per day.

#### APPLICATIONS

**Direct Assays:** GLO-1 activity in enzyme preparations or biological samples.

**Drug Discovery/Pharmacology:** effects of drugs on GLO-1 activity.

#### KIT CONTENTS

**Assay Buffer (pH 6.6):** 20 mL    **Substrate:** 1 mL  
**96 well UV Titer Plate** 1 Plate    **Cosubstrate:** 1 mL

**Storage conditions.** Kit is shipped at room temperature. Store the plate at room temperature and other components at -20°C. 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.

#### PROCEDURE FOR PURE ENZYME PREPARATIONS

- Bring all reagents to room temperature prior to assay.
- Add 40 µL of each sample to separate wells of the 96 well UV titer plate. Also, include one blank (sample buffer without GLO-1) well per assay run.
- Prepare Working Reagent for all wells by mixing per well: 160 µL Assay Buffer with 8 µL Substrate and 8 µL Cosubstrate. Add 160 µL Working Reagent to each well.
- Read the optical density at 240 nm at t=0 min and again at t=10 min. If measuring low GLO-1 activities, longer reaction times can be used.

#### PROCEDURE FOR PROTEINOUS SAMPLES

- Each sample requires 2 tubes: one for the GLO-1 Reaction and one for the Sample Blank. Add 40 µL of each sample (serum samples should be diluted at least 2x with assay buffer) to 2 separate Eppendorf tubes.
- Prepare Working Reagent for all tubes by mixing per tube: 160 µL Assay Buffer with 8 µL Substrate and 8 µL Cosubstrate.
- GLO-1 Reaction.** Add 160 µL Working Reagent to the GLO-1 Reaction tubes. Incubate for 20 min at room temperature.
- Protein Precipitation of GLO-1 Reaction.** After 20 min incubation, add 70 µL 4 M Perchloric Acid to each GLO-1 reaction tube, vortex to mix and chill for 15 min on ice. (Note: *It is important that the protein precipitation is carried out on ice.*) After 15 min centrifuge samples for 5 min at 14,000 rpm and transfer 200 µL of each clear supernatant to separate wells of the 96 well UV titer plate.

- Sample Blank.** Add 70 µL 4 M Perchloric Acid to each Sample Blank tube, vortex to mix and *chill* for 15 min on ice. After 15 min, add 160 µL Working Reagent to each Sample Blank tube. Vortex to mix and *chill* on ice for 15 min. (Note: *It is important that the sample is deproteinated prior to adding the Working Reagent.*) After the second chill, centrifuge the sample blanks for 5 min at 14,000 rpm and transfer 200 µL of each clear supernatant to separate wells of the 96 well UV titer plate.

- Read optical density at 240 nm.

#### CALCULATION

For pure samples glyoxalase activity is calculated as follows:

$$\text{GLO-1} = \frac{\text{OD}_{10} - \text{OD}_0}{\epsilon \times l} \times \frac{V_T}{t} \times \frac{1}{V_S} = 350 \times (\text{OD}_{10} - \text{OD}_0) \text{ (U/L)}$$

where  $\text{OD}_{10}$  and  $\text{OD}_0$  are the optical density values of the sample taken at 10 min and 0 min respectively.  $V_T$  is the total reaction volume (0.2 mL),  $V_S$  is the sample volume (40 µL),  $\epsilon$  is the S-lactoylglutathione extinction coefficient ( $3.37 \text{ mM}^{-1}\text{cm}^{-1}$ ),  $l$  is the path length (0.425 cm for 0.2 mL in provided plate) and  $t$  is the reaction time (10 min).

For proteinous samples glyoxalase activity is calculated as follows:

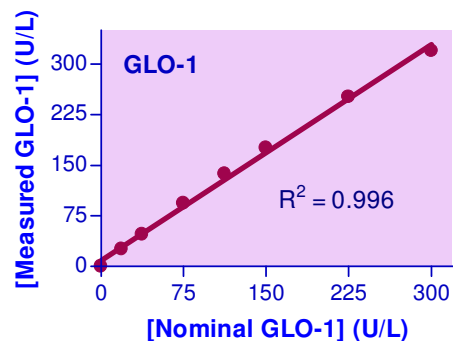
$$\begin{aligned} \text{GLO-1} &= \frac{\text{OD}_{\text{SAMPLE}} - \text{OD}_{\text{BLANK}}}{\epsilon \times l} \times \frac{V_T}{t} \times \frac{1}{V_S} \times 1.35 \times n \\ &= 175 \times (\text{OD}_{\text{SAMPLE}} - \text{OD}_{\text{BLANK}}) \times 1.35 \times n \text{ (U/L)} \end{aligned}$$

where  $\text{OD}_{\text{SAMPLE}}$  and  $\text{OD}_{\text{BLANK}}$  are the optical density values of the sample and sample blank respectively.  $t$  is the reaction time (20 min),  $1.35$  is the dilution factor for the deproteination step and  $n$  is the dilution factor if a sample dilution is required.

**Unit definition:** 1 unit of Glyoxalase-1 forms 1 µmole of S-lactoylglutathione from methylglyoxal and reduced glutathione per minute at pH 6.6 and 25°C.

#### MATERIALS REQUIRED, BUT NOT PROVIDED

Perchloric Acid (e.g. Sigma Cat. No. 311413), pipetting devices and accessories and plate reader.



#### REFERENCES

- Davis, KA and Williams, GR. (1969) Glyoxalase I, a lyase or an oxidoreductive isomerase? *Can. J. Biochem* 47: 553-6.
- Ditzen, C et al. (2006) Protein biomarkers in a mouse model of extremes in trait anxiety. *Mol Cell Proteomics* 5: 1914-20.
- Strzinek, RA. et al. (1972) The purification and characterization of liver glyoxalase I from normal mice and from mice bearing a lymphosarcoma. *Cancer Res* 32:2359-64.

