

## QuantiFluo™ Glutaminase Assay Kit (DGLN-100)

### Quantitative Fluorimetric Assay for Glutaminase Activity

#### DESCRIPTION

**GLUTAMINASE** (EC 3.5.1.2) is a key enzyme in the metabolism of glutamine. Glutamine is an important amino acid in the metabolism of cancer cells, and increased glutaminolysis has been linked to cancer. BioAssay Systems' DGLN-100 Kit provides a convenient fluorimetric method to measure glutaminase activity in biological samples. In this assay, o-phthalaldehyde reacts with liberated ammonia where the increase in fluorescence at  $\lambda_{\text{ex/em}} = 415/475$  nm is directly proportional to enzyme activity.

#### KEY FEATURES

**Fast and Safe.** Assay can be completed within 50 minutes. Non-radioactive assay.

**Sensitive and accurate.** Linear detection range is 0.66-500 U/L glutaminase in a 96-well plate assay.

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

#### APPLICATIONS

For quantitative determination of glutaminase activity in biological samples.

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

**Assay Buffer:** 5 mL      **20 mM Standard:** 50  $\mu$ L  
**20 mM Substrate:** 500  $\mu$ L      **Detection Reagent:** 5 mL

**Storage conditions.** The kit is shipped at room temperature. Store all components at  $-20^{\circ}\text{C}$  upon receipt. 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 the Material Safety Data Sheet for detailed information.

#### PROCEDURES

This assay is based on a kinetic reaction. To ensure identical incubation time, addition of the Working Reagent (*WR*) to samples should be quick and mixing should be brief but thorough. Use of a multi-channel pipettor is recommended. The assay can be run at room temperature.

**Reagent Preparation:** Prior to the assay, equilibrate all components to room temperature and briefly centrifuge tubes before opening. The *WR* should be prepared fresh for each assay run.

**Enzyme Preparation:** Enzyme should be prepared in an enzyme buffer, e.g. 20 mM Potassium Phosphate, pH 7.4. The following protocol is optimized for glutaminase from *Bacillus amyloliquefaciens*. If using a different enzyme, we recommend that you experimentally determine the optimal amount of enzyme to use per well.

#### Sample Preparation:

*Serum* and *Plasma* samples must be diluted at least 1:10.

*Urine* must be diluted 1:50 in water prior to the assay run.

**Standard Preparation:** Prepare a 2 mM Standard Premix by combining 20  $\mu$ L 20 mM Standard and 180  $\mu$ L Assay Buffer. Prepare all standards according to table below. Transfer 10  $\mu$ L of Standard to each well plus 40  $\mu$ L of Assay Buffer.

No.	2 mM Std Premix + Assay Buffer	Total Volume ( $\mu$ L)	Std (mM)
1	100 $\mu$ L + 0 $\mu$ L	100 $\mu$ L	2 mM
2	50 $\mu$ L + 50 $\mu$ L	100 $\mu$ L	1 mM
3	25 $\mu$ L + 75 $\mu$ L	100 $\mu$ L	0.5 mM
4	0 $\mu$ L + 100 $\mu$ L	100 $\mu$ L	0 mM

#### Reaction Preparation:

- Transfer 10  $\mu$ L of each sample to separate wells of the plate.
- Prepare enough *WR* for all sample wells by mixing 5  $\mu$ L of 20 mM Substrate and 45  $\mu$ L of Assay Buffer for each well.
- Initiate the reaction by addition of 40  $\mu$ L of *WR* to all sample wells. Incubate the reaction for 30 minutes at RT.
- Add 50  $\mu$ L *Detection Reagent* to all wells. Tap plate to mix and read for 20 min. Measure fluorescence intensity at  $\lambda_{\text{ex/em}} = 415/475$  nm.

#### CALCULATION

Subtract the blank value (Standard #4) from the standard values and plot  $\Delta F$  against the standard concentrations. Determine the slope ( $\text{mM}^{-1}$ ) and calculate the glutaminase activity in each Sample as follows,

$$\text{Glutaminase Activity} = \frac{(F_{\text{Sample}} - F_{\text{Blank}})}{\text{Slope (mM}^{-1})} \times \frac{\text{Reaction Vol}}{t \text{ (min)} \times \text{Enzyme Vol (}\mu\text{L)}} \times n \text{ (U/L)}$$

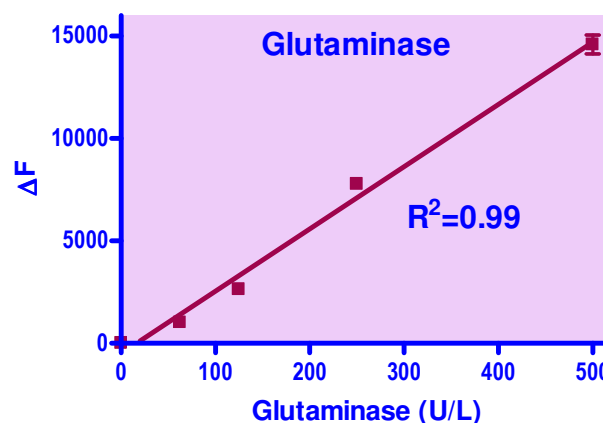
Where  $F_{\text{Sample}}$  and  $F_{\text{Blank}}$  are the measured fluorescence values of the sample and blank,  $t$  is the reaction time (30 min), Reaction Vol and Enzyme Vol are 50  $\mu$ L and 10  $\mu$ L, respectively, and  $n$  is the sample dilution factor.

**Unit definition:** 1 Unit (U) of glutaminase will catalyze the conversion of 1 micromole L-glutamine per min at room temperature and pH 7.4.

**Note:** If sample glutaminase activity exceeds 500 U/L, dilute samples in enzyme buffer and repeat the assay.

#### MATERIALS REQUIRED, BUT NOT PROVIDED

Pipetting devices and accessories (e.g. multi-channel pipettor), black, flat-bottom 96-well plates (e.g. Corning Costar), centrifuge tubes and a plate reader.



Varying amounts of Glutaminase were assayed in the presence of 2 mM Glutamine for 30 minutes at RT.

#### LITERATURE

- Milano, S. K., et al (2022). New insights into the molecular mechanisms of glutaminase C inhibitors in cancer cells using serial room temperature crystallography. *Journal of Biological Chemistry*, 298(2), 101535.
- Sugawara, K, Oyama, F. (1981). Fluorogenic reaction and specific microdetermination of ammonia. *J. Biochem.* 89, 771–774.

