

EnzyChrom™ Glutamine Assay Kit (EGLN-100)

Quantitative Colorimetric Determination of Glutamine at 565 nm

DESCRIPTION

Glutamine is an amino acid synthesized in the muscle that plays major roles in protein synthesis, acid-base balance, anabolic processes and is utilized for cellular energy and as a carbon source. It is used in treatment of injury, trauma, burns, and also as a supplement for muscle growth and post-surgery healing.

Simple, direct and automation-ready procedures for measuring glutamine concentration are very desirable. BioAssay Systems' EnzyChrom™ glutamine assay kit is based on hydrolysis of glutamine to glutamate and colorimetric determination of the product. The intensity of the product color, measured at 565 nm, is proportional to the glutamine concentration in the sample.

APPLICATIONS

Direct Assays: glutamine in serum, plasma, urine, tissue extracts and cell culture samples.

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

KEY FEATURES

Sensitive and accurate. Use 20 μL sample. Linear detection range 0.023 - 2 mM glutamine in 96-well plate assay.

Convenient. The procedure involves adding a single working reagent, incubation for 40 min at room temperature, adding a stop reagent and reading the optical density. No 37°C heater is needed.

High-throughput. Can be readily automated as a high-throughput 96-well plate assay for thousands of samples per day.

KIT CONTENTS (100 tests in 96-well plates)

Assay Buffer: 15 mL **NAD Solution:** 1 mL
Enzyme A: 120 μL **MTT Solution:** 2 x 1.5 mL
Enzyme B: 220 μL **Stop Reagent:** 25 mL
Standard: 400 μL 100 mM Glutamine

Storage conditions. The kit is shipped on ice. Store all reagents 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.

PROCEDURE

Note: (1). this assay is based on an enzyme-catalyzed kinetic reaction. Addition of Working Reagent should be quick and mixing should be brief but thorough. Use of multi-channel pipettor is recommended. (2). the following substances interfere and should be avoided in sample preparation: ascorbic acid, SDS (>0.2%), sodium azide, NP-40 (>1%) and Tween-20 (>1%).

- Standard Curve.** Prepare 2.0 mM glutamine Premix by mixing 5 μL 100 mM Standard and 245 μL distilled water. Dilute standard as follows. Transfer 20 μL standards into wells of a clear flat-bottom 96-well plate.

No	Premix + H ₂ O	Vol (μL)	Glutamine (mM)
1	100 μL + 0 μL	100	2.0
2	60 μL + 40 μL	100	1.2
3	30 μL + 70 μL	100	0.6
4	0 μL + 100 μL	100	0.0

Samples: add 20 μL sample per well in separate wells.

IMPORTANT: if a sample is known to contain glutamate, a sample blank control is required. In this case, transfer an additional 20 μL sample into a separate well.

- Reaction.** Spin the enzyme and reagent tubes briefly before pipetting. Fresh reconstitution is recommended.

For each standard and sample well, prepare Working Reagent by mixing 65 μL Assay Buffer, 1 μL Enzyme A, 1 μL Enzyme B, 2.5 μL NAD and 14 μL MTT.

Where a sample blank is required, prepare a Blank Working Reagent by mixing 65 μL Assay Buffer, 1 μL Enzyme B, 2.5 μL NAD and 14 μL MTT (i.e. No Enzyme A).

Add 80 μL Working Reagent per well to standards and sample wells.

Where appropriate, add 80 μL Blank Working Reagent to the Sample Blank wells.

Tap plate to mix briefly and thoroughly.

- Incubate 40 min at room temperature. Add 100 μL Stop Reagent to each well. Read OD at 565 nm (520-600 nm).

CALCULATION

Subtract water (#4) blank OD from OD values for the standards. Plot ΔOD against standard concentrations. Determine the slope and calculate sample glutamine concentration,

$$[\text{Glutamine}] = \frac{\text{OD}_{\text{SAMPLE}} - \text{OD}_{\text{BLANK}}}{\text{Slope (mM}^{-1})} \times n \quad (\text{mM})$$

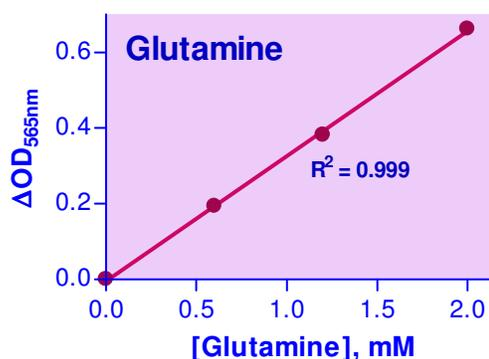
where $\text{OD}_{\text{SAMPLE}}$ and OD_{BLANK} are the OD values of the sample and water (if sample does not contain glutamate) or sample blank (if sample contains glutamate).

Note: if the calculated glutamine concentration is higher than 2 mM, dilute sample in distilled water and repeat the assay. Multiply the results by the dilution factor n .

Conversions: 1 mM glutamine = 14.6 mg/dL or 146 ppm.

MATERIALS REQUIRED, BUT NOT PROVIDED

Pipeting (multi-channel) devices. Clear flat-bottom 96-well plates (e.g. Corning Costar) and plate reader.



Standard Curve in 96-well plate assay

PUBLICATIONS

- Velickovic, K., et al. (2020). Targeting glutamine synthesis inhibits stem cell adipogenesis in vitro. Cellular Physiology and Biochemistry, 54(5), 917-927.
- Tran, T. Q., et al. (2020). α -Ketoglutarate attenuates Wnt signaling and drives differentiation in colorectal cancer. Nature Cancer. 1(3): 345-358.
- Ma, Huanrong, et al (2018). Inhibition of SLC1A 5 sensitizes colorectal cancer to cetuximab. International Journal of Cancer 142.12: 2578-2588.

