

Glutamine Assay Kit (Colorimetric) ab197011

★★★★★ 1 Abreviews 31 References 3 Images

Overview

Product name	Glutamine Assay Kit (Colorimetric)
Detection method	Colorimetric
Sample type	Urine, Serum, Plasma, Other biological fluids, Tissue
Assay type	Quantitative
Sensitivity	25 µM
Assay time	0h 90m
Species reactivity	<b>Reacts with:</b> Mammals, Other species
Product overview	<p>Glutamine Assay kit (Colorimetric) ab197011 is a simple and sensitive assay that detects the biologically relevant concentrations of glutamine in biological fluids and tissues.</p> <p>The glutamine assay protocol is based on the hydrolysis of Glutamine to Glutamate producing a stable colored signal, which is directly proportional to the amount of glutamine in the sample.</p> <p>The assay can detect as little as 25 µM of glutamine in a variety of biological samples.</p> <p>Glutamine assay protocol summary:</p> <ul style="list-style-type: none"><li>- add samples and standards to wells</li><li>- add hydrolysis mix and incubate for 30 min</li><li>- add reaction mix and incubate for 60 min</li><li>- analyze with a microplate reader.</li></ul>
Notes	<p>This product is manufactured by BioVision, an Abcam company and was previously called K556 Glutamine Colorimetric Assay Kit. K556-100 is the same size as the 100 test size of ab197011.</p>
Platform	Microplate reader

Properties

Storage instructions Store at -20°C. Please refer to protocols.

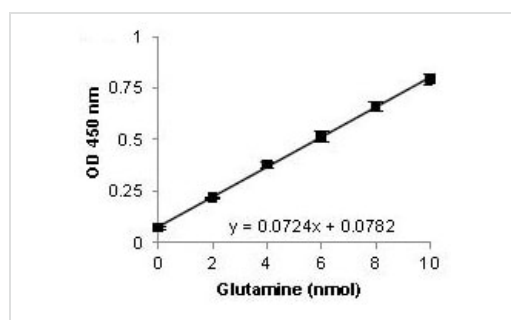
Components	100 tests
Assay Buffer XXIX	1 x 25ml

Components	100 tests
Assay Buffer XXXI	1 x 25ml
Developer Solution III	1 vial
Enzyme Mix VIII	1 vial
Glutamine Standard	1 vial
Hydrolysis Enzyme Mix II	1 vial

## Relevance

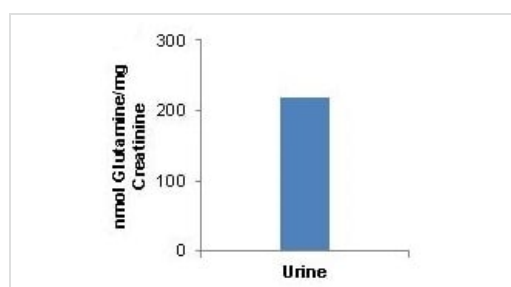
Glutamine (abbreviated as Gln or Q; the abbreviation Glx or Z represents either glutamine or glutamic acid) is one of the 20 amino acids encoded by the standard genetic code. Its side chain is an amide formed by replacing the side-chain hydroxyl of glutamic acid with an amine functional group. It can therefore be considered the amide of glutamic acid. Its codons are CAA and CAG. Recent studies have proven its effectiveness in anabolic muscle growth from prolonged consumption.

## Images



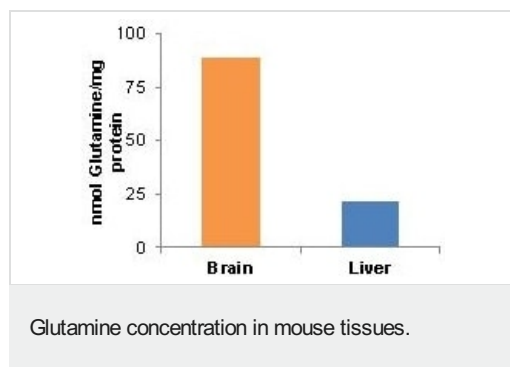
Glutamine Standard Curve. For demonstrative purposes only.

Glutamine Standard Curve.



Measurement of Glutamine concentration in Human urine (5  $\mu$ l).

Glutamine concentration in Human urine.



Measurement of Glutamine concentration in Mouse tissues (brain [15 µg], liver [50 µg]). Samples were deproteinized using 10K Spin Column and spiked with known amount of Glutamine (6 nmol).

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### Our Abpromise to you: Quality guaranteed and expert technical support

---

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

### Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors