

Glutamate Assay Kit ab83389

★★★★★ [2 Abreviews](#) [43 References](#) [画像数 2](#)

製品の概要

製品名	Glutamate Assay Kit
検出方法	Colorimetric
サンプルの種類	Urine, Serum, Plasma, Other biological fluids, Cell Lysate
アッセイタイプ	Quantitative
全工程の試験時間	0h 40m
製品の概要	Glutamate Assay Kit ab83389 provides a sensitive detection method of the glutamate in a variety of samples.

The glutamate assay kit will only measure free glutamate levels but not glutamic acid found in the backbone of peptides or proteins.

In the glutamate assay protocol, the enzyme mix recognizes glutamate as a specific substrate leading to proportional color development. The amount of glutamate can therefore be easily quantified by colorimetric spectrophotometry at OD = 450 nm.

Glutamate assay protocol summary:

- add samples and standards to wells
- add reaction mix and incubate for 30 min at 37°C
- analyze with microplate reader

特記事項

This product is manufactured by BioVision, an Abcam company and was previously called K629 Glutamate Colorimetric Assay Kit. K629-100 is the same size as the 100 test size of ab83389.

EDTA plasma may not be used with the ab83389 Glutamate Assay Kit. EDTA can interfere with the assay.

Review our [Metabolism Assay Guide](#) to learn about assays for metabolites, metabolic enzymes, mitochondrial function, and oxidative stress, and also about how to assay metabolic function in live cells using your plate reader.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

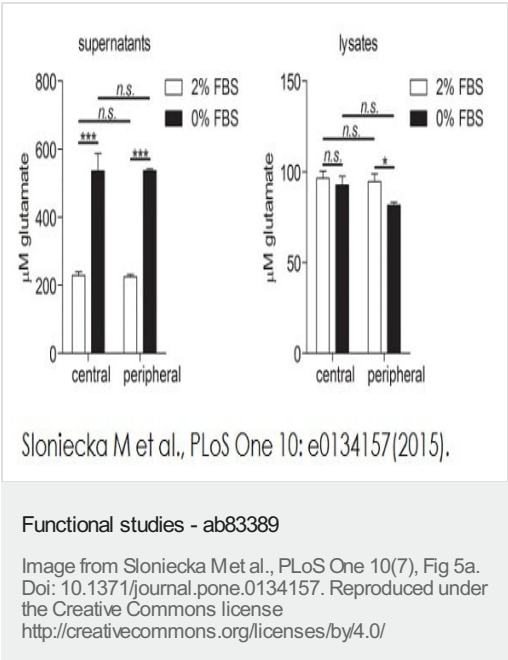
製品の特性

保存方法Store at -20°C. Please refer to protocols.

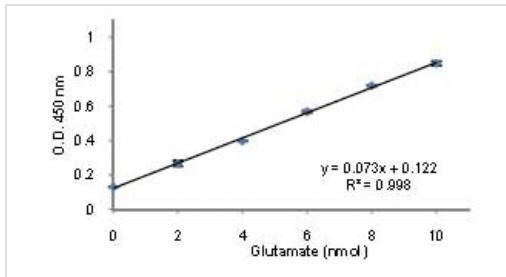
内容	100 tests
Assay Buffer XXXI	1 x 25ml
Developer Solution III	1 vial
Enzyme Mix VIII	1 vial
Glutamate Standard	1 x 100µl

関連性Glutamate, one of the two acidic proteinogenic amino acids, is also a key molecule in cellular metabolism. In humans, glutamate plays an important role both in amino acid degradation and disposal of excess or waste nitrogen. Glutamate is the most abundant swift excitatory neurotransmitter in the mammalian nervous system. It is believed to be involved in learning and memory and has appeared to be involved in diseases like amyotrophic lateral sclerosis, lathyrism, autism, some forms of mental retardation and Alzheimer's disease. Glutamic acid is also present in a wide variety of foods, and has been used as a flavor enhancer in food industry.

画像



Glutamate levels of central and peripheral keratocytes were determined using Glutamate assay kit (ab83389).



Example of Glutamate standard curve obtained using ab83389

Functional Studies - Glutamate Assay Kit (ab83389)

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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