

Phenylalanine Assay Kit ab83376

画像数 3

製品の概要

製品名	Phenylalanine Assay Kit
検出方法	Fluorescent
サンプルの種類	Urine, Serum, Plasma, Other biological fluids, Tissue Extracts
アッセイタイプ	Quantitative
検出感度	> 2 μ M
検出範囲	2 μ M - 20 μ M
全工程の試験時間	1h 00m
製品の概要	Phenylalanine Assay Kit ab83376 provides a quick, simple, accurate method for quantifying phenylalanine in biological samples.

In the phenylalanine assay protocol, phenylalanine is reductively deaminated with the simultaneous formation of NADH which reacts with our fluorescent probe to generate fluorescence at Ex/Em=535/587 nm. The assay is linear in the range from 0.1 to 1.0 nmol (2-20 μ M) of Phenylalanine.

Phenylalanine assay protocol summary:

- add samples and standards to wells
- add reaction mix and incubate for 60 min
- analyze with a microplate reader

特記事項

This product is manufactured by BioVision, an Abcam company and was previously called K572 Phenylalanine Fluorometric Assay Kit. K572-100 is the same size as the 100 test size of ab83376.

L-Phenylalanine (PHE) is an electrically-neutral amino acid, one of the twenty common and one of the three aromatic amino acids used to biochemically form proteins. Phenylalanine uses the same active transport channel as tryptophan to cross the blood-brain barrier, and, in large quantities, interferes with the production of serotonin.

試験プラットフォーム

Microplate reader

製品の特性

保存方法

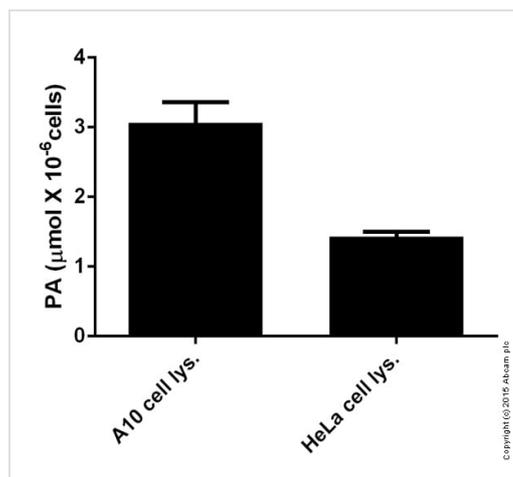
Store at -20°C. Please refer to protocols.

内容	100 tests
Phenylalanine Standard	1 vial
Fructose Assay Buffer	1 x 25ml
Phenylalanine Developer Mix	1 vial
Phenylalanine Enzyme Mix	1 vial
Tyrosinase	1 vial

関連性

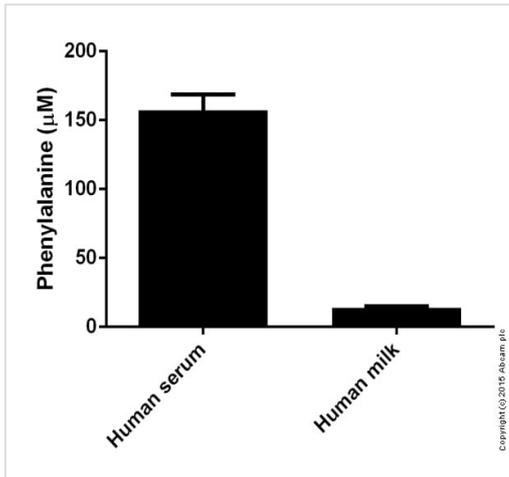
L-Phenylalanine (PHE) is an electrically-neutral amino acid, one of the twenty common and one of the three aromatic amino acids used to biochemically form proteins. Phenylalanine uses the same active transport channel as tryptophan to cross the blood-brain barrier, and, in large quantities, interferes with the production of serotonin. Errors in PHE metabolism lead to phenylketonuria or PKU which can have dire consequences.

画像



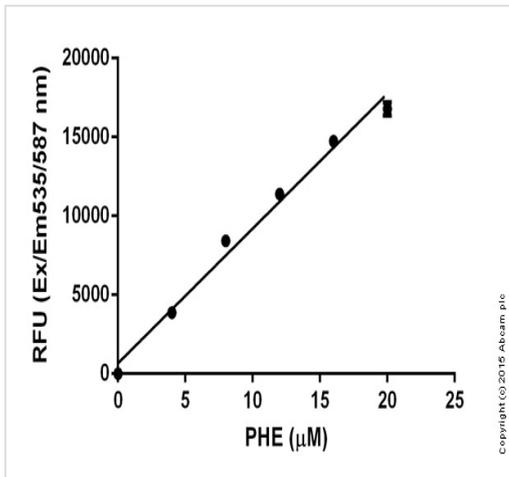
Phenylalanine measured in cell lysates showing quantities (micromoles) per million of cells tested.

OEM testing Phase VI ; Kits & Assays Initiative
Assay Testing Jan/Feb 2015



Phenylalanine measured in biofluids showing concentration in micromolar.

Functional assay - ab83376 Phenylalanine Assay Kit



Phenylalanine fluorimetric standard curve: mean of duplicates (+/- SD) with background readings subtracted

Functional assay - ab83376 Phenylalanine Assay Kit

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