

Branched Chain Amino Acid Assay Kit / BCAA Assay Kit ab83374

21 References **画像数 4**

製品の概要

製品名	Branched Chain Amino Acid Assay Kit / BCAA Assay Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Urine, Serum, Plasma, Other biological fluids
アッセイタイプ	Quantitative
検出感度	> 0.2 nmol/well
全工程の試験時間	0h 40m
製品の概要	<p>Branched Chain Amino Acid Assay Kit / BCAA Assay Kit ab83374 provides a simple convenient means of measuring the BCAA's in biological samples.</p> <p>The BCAA assay protocol uses an enzyme reaction in which BCAA is oxidatively deaminated, producing NADH which reduces a probe, generating a colored product ($\lambda_{\text{max}} = 450 \text{ nm}$).</p> <p>The Branched Chain Amino Acid Assay Kit measures BCAA's in the range of 0 to 10 nmol per sample with a detection limit of ~0.2 nmol (~10 μM BCAA in sample). BCAA's are present in serum ~ 0.1-0.4 mM each (~0.125-1.5 mM combined).</p> <p>BCAA assay protocol summary:</p> <ul style="list-style-type: none"> - add samples and standards to wells - add reaction mix and incubate for 30 min - analyze with a microplate reader
特記事項	<p>This product is manufactured by BioVision, an Abcam company and was previously called K564 Branched Chain Amino Acid (Leu/Ile/Val) Colorimetric Assay Kit. K564-100 is the same size as the 100 test size of ab83374.</p> <p>The branched-chain amino acids or BCAA's, refer to the amino acids with non-linear aliphatic side-chains, namely leucine, isoleucine and valine. These three essential amino acids make up approximately 1/3 of skeletal muscle in the human body.</p>
試験プラットフォーム	Microplate reader

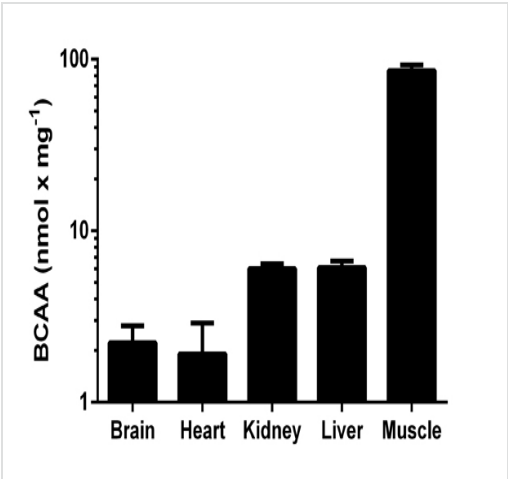
製品の特性

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

内容	100 tests
Assay Buffer IV	1 x 25ml
BCAA Enzyme Mix	1 vial
Developer Solution III	1 vial
Leucine Standard	1 x 100µl

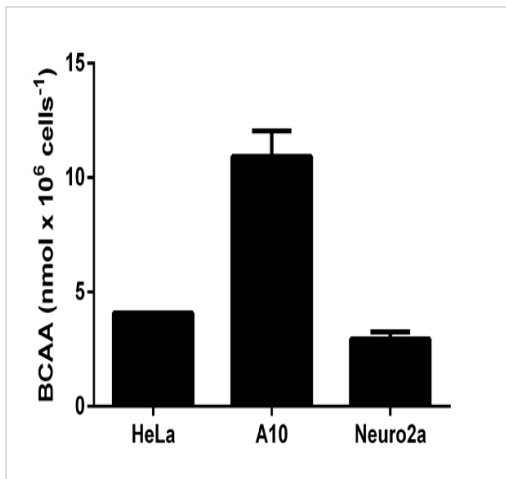
関連性The branched-chain amino acids or BCAA's, refer to the amino acids with non-linear aliphatic side-chains, namely leucine, isoleucine and valine. These three essential amino acids make up approximately 1/3 of skeletal muscle in the human body. BCAA's are currently used clinically to aid in the recovery of burn victims, as well as for strength supplementation for athletes. BCAA's, primarily Leu, can stimulate insulin secretion. The BCAA's have also been implicated in a wide range of other physiological effects.

画像



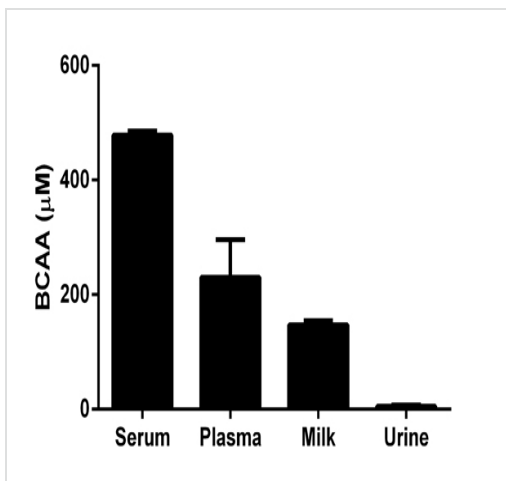
BRCAA levels measured in mouse tissue lysates (mg of extracted protein; background signal subtracted, mean of duplicates; +/- SD).

Functional Studies - Branched Chain Amino Acid Assay Kit (ab83374)



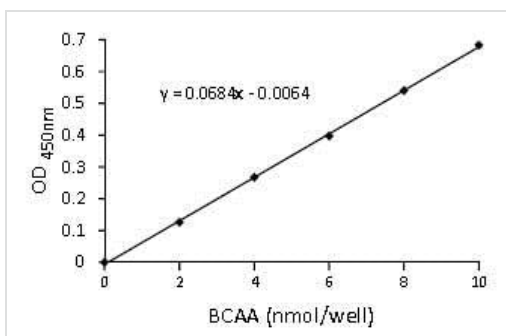
BCAA levels measured in cell lysates (background signal subtracted, mean of duplicates; +/- SD).

Functional Studies - Branched Chain Amino Acid Assay Kit (ab83374)



BCAA levels measured in human biological fluids (background signal subtracted, mean of duplicates; +/- SD).

Functional Studies - Branched Chain Amino Acid Assay Kit (ab83374)



Leucine Standard Curve performed according to the attached protocol.

Functional Studies - Branched Chain Amino Acid Assay Kit (ab83374)

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