

Choline/Acetylcholine Assay Kit ab65345

56 References **画像数 5**

製品の概要

製品名	Choline/Acetylcholine Assay Kit
検出方法	Colorimetric/Fluorometric
サンプルの種類	Cell culture supernatant, Urine, Serum, Plasma, Cell culture extracts, Other biological fluids, Tissue Extracts
アッセイタイプ	Quantitative
検出感度	> 0.01 nmol/well
検出範囲	0.01 nmol/well - 5 nmol/well
全工程の試験時間	0h 40m
種交差性	交差種: Mammals, Other species
製品の概要	Choline/Acetylcholine Assay Kit (ab65345) provides a simple and sensitive means for quantifying Choline and Acetylcholine by either a colorimetric or fluorometric method in samples such as blood, cells, culture media, fermentation media, etc.

In the choline / acetylcholine assay protocol, free choline is oxidized to betaine, via the intermediate betaine aldehyde. The reaction generates products which react with the Choline Probe to generate color ($\lambda = 570$ nm), and fluorescence (Ex/Em 535/587 nm).

Acetylcholine can be converted to choline by adding acetylcholinesterase to the reaction to measure total choline (choline + acetyl choline). The amount of acetyl choline is calculated by subtracting choline from total choline.

There is no need for pretreatment or purification of samples. The kit can detect ~10 pmol-5 nmol of choline or acetylcholine.

Cells should be cultured in choline chloride free medium.

Choline / acetylcholine assay protocol summary:

- add samples and standards to wells
- add choline reaction mix with, or without, acetylcholinesterase, and incubate for 30 min at room temp
- analyze with microplate reader

特記事項 This product is manufactured by BioVision, an Abcam company and was previously called K615

Choline/Acetylcholine Quantification Colorimetric/Fluorometric Kit. K615-100 is the same size as the 100 test size of ab65345.

If you need to purchase additional Choline Probe for this product, please see [ab139505](#).

試験プラットフォーム

Microplate reader

製品の特性

保存方法

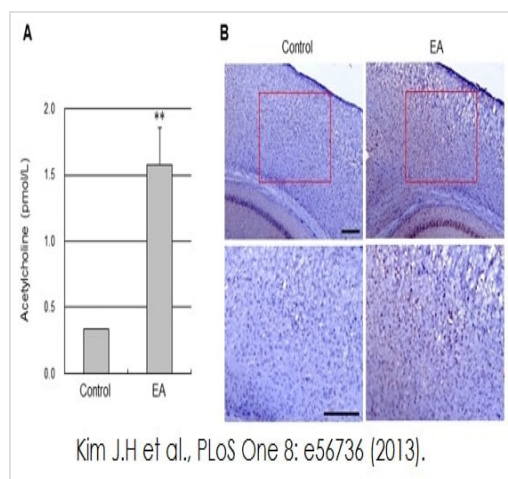
Store at -20°C. Please refer to protocols.

内容	100 tests	2000 tests
Acetylcholinesterase Enzyme	1 vial	20 vials
Assay Buffer VI	1 x 25ml	20 x 25ml
Choline Standard	1 vial	20 vials
Enzyme Mix IV	1 vial	20 vials
OxiRed Probe	1 x 200µl	20 x 200µl

関連性

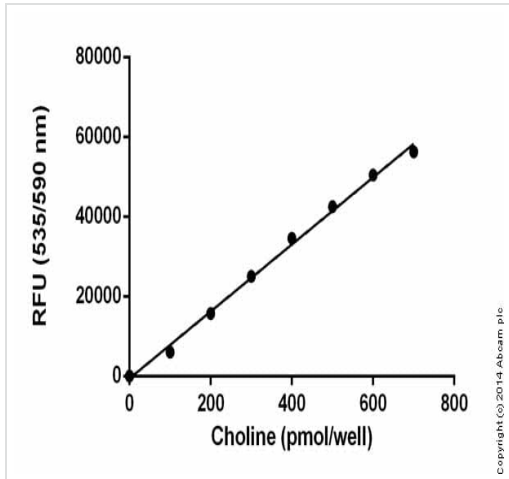
Choline and acetylcholine (often abbreviated ACh) play important roles in many biological processes. Choline is a natural amine, and is generally found in phosphatidylcholine and sphingomyelin, phospholipids abundant in cell membranes. Choline is also the precursor molecule for the neurotransmitter acetylcholine which is involved in many functions both in the peripheral nervous system (PNS) and in the central nervous system (CNS). In the PNS, ACh activates muscles and is a major neurotransmitter in the autonomic nervous system. In the CNS, ACh and the associated neurons form the cholinergic system, which tends to cause anti-excitatory actions.

画像

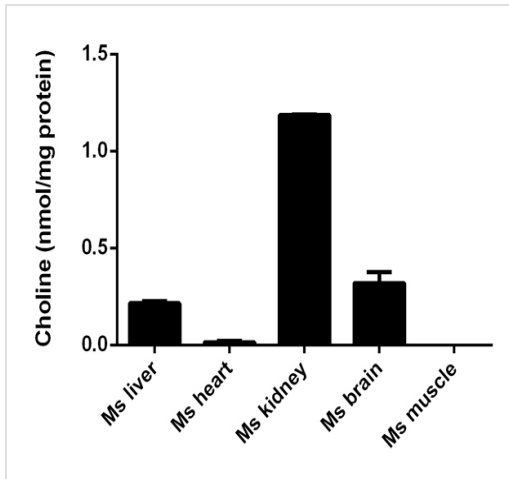


Functional studies- ab65345

Image from Kim J.H et al., PLoS One 8(2), Fig 4. doi: 10.1371/journal.pone.0056736. Reproduced under the Creative Commons license <http://creativecommons.org/licenses/by/4.0/>

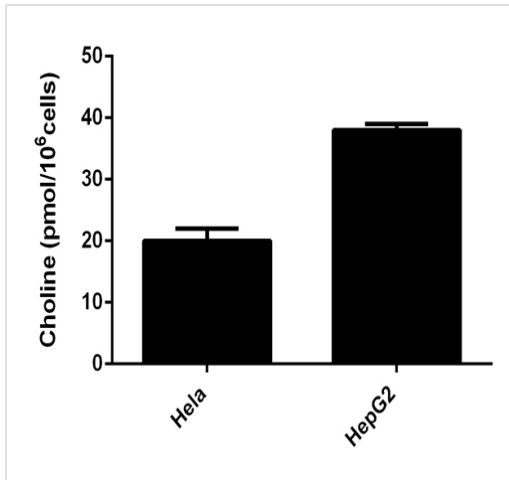


Functional Studies - Choline/Acetylcholine Assay
Kit (ab65345)



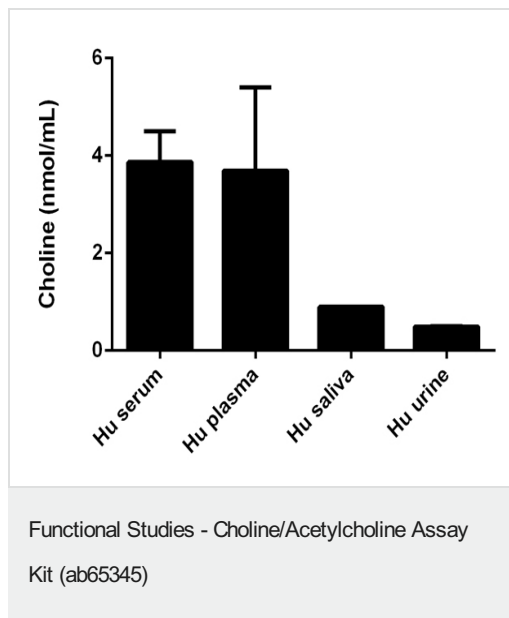
Choline measured fluorometrically in mouse tissue lysates showing quantity (nmol) per mg protein of tested sample

Functional Studies - Choline/Acetylcholine Assay
Kit (ab65345)



Choline measured fluorometrically in cell lysates showing quantity (pmol) per 1 mln of tested cells

Functional Studies - Choline/Acetylcholine Assay
Kit (ab65345)



Choline measured fluorometrically in biological fluids showing quantity (nmol) per mL of tested sample

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