

Acetylcholinesterase Assay Kit (Fluorometric -Red) ab138873

★★★★★ [1 Abreviews](#) [3 References](#) [画像数 3](#)

製品の概要

製品名	Acetylcholinesterase Assay Kit (Fluorometric -Red)
検出方法	Colorimetric/Fluorometric
サンプルの種類	Serum, Plasma, Cell Lysate
アッセイタイプ	Enzyme activity
検出感度	0.1 mU/ml
全工程の試験時間	0h 20m
種交差性	交差種: Mammals, Other species
製品の概要	<p>Acetylcholinesterase Assay Kit (Fluorometric -Red) (ab138873) uses a red indicator to quantify the choline produced from the hydrolysis of acetylcholine by AChE through choline oxidase-mediated enzyme coupling reactions.</p> <p>The assay can be used for monitoring and quantifying AChE activity in blood, cell extracts or other solutions. The fluorescence intensity of the indicator is proportional to the amount of choline formed, which is proportional to the AChE activity.</p> <p>The assay is an optimized “mix and read” assay using a simple one-step fluorometric protocol to detect as little as 0.01 mU AChE in a 100 µl assay volume (0.1 mU/ml). The assay can be performed in a 96-well or 384-well microtiter-plate format, and can be easily read with a fluorescence microplate reader at Ex/Em = 540/590 nm or with an absorbance microplate reader at OD = 575 nm.</p> <p>Acetylcholinesterase assay protocol summary:</p> <ul style="list-style-type: none">- add samples and standards to wells- add assay mixture and incubate for 10-30 min- analyze with microplate reader
特記事項	<p>Please note this product does not differentiate between acetylcholinesterase (AChE) or butyrylcholinesterase (BChE) activity as both enzymes can hydrolyze acetylcholine.</p> <p>If you are looking for a colorimetric kit, we recommend Acetylcholinesterase Assay Kit (Colorimetric) (ab138871). If your plate reader has a filter to measure green fluorescence (Ex/Em = 490/520 nm), you can use Acetylcholinesterase Assay Kit (Fluorometric - Green) (ab138872).</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	200 tests
AbRed Indicator	1 vial
Acetylcholine	1 vial
Acetylcholinesterase Probe	2 units
Acetylcholinesterase Standard	1 vial
Assay Buffer	1 x 10ml
Dilution Buffer	1 x 10ml
DMSO	1 x 100µl

機能

Terminates signal transduction at the neuromuscular junction by rapid hydrolysis of the acetylcholine released into the synaptic cleft. Role in neuronal apoptosis.

組織特異性

Isoform H is highly expressed in erythrocytes.

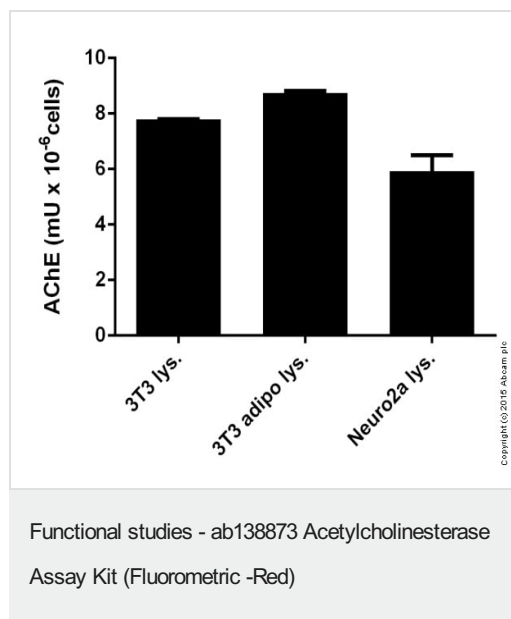
配列類似性

Belongs to the type-B carboxylesterase/lipase family.

細胞内局在

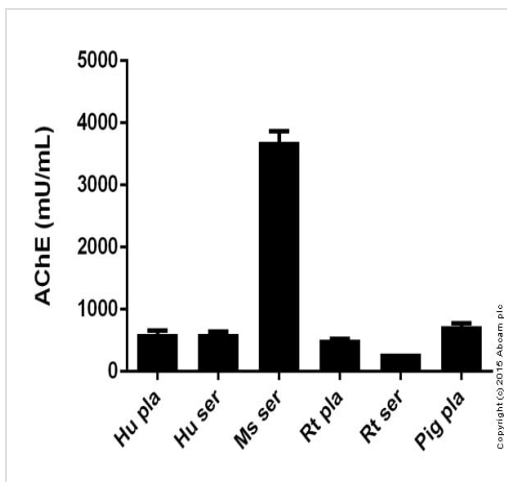
Cell membrane; Cell junction > synapse. Secreted. Cell membrane and Nucleus. Only observed in apoptotic nuclei.

画像



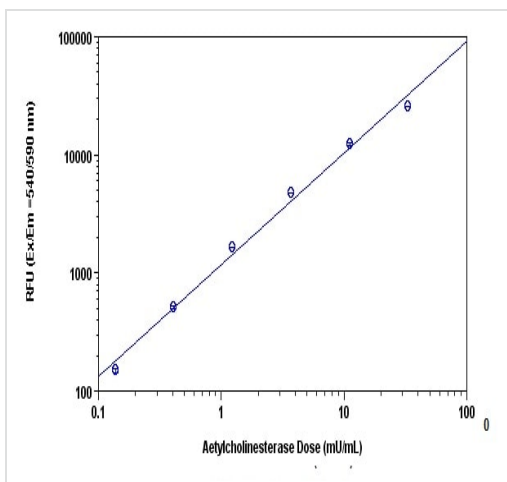
AChE measured in cell lysates showing activity per 1 million cells

Functional studies - ab138873 Acetylcholinesterase Assay Kit (Fluorometric -Red)



AChE measured in various species showing activity per mL sample tested

Functional studies - ab138873 Acetylcholinesterase Assay Kit (Fluorometric -Red)



Acetylcholinesterase dose response was measured in a solid black 96-well plate with ab138873 using a fluorescence microplate reader. As low as 0.01 mU/well (0.1 mU/ml) acetylcholinesterase can be detected with 20 minutes incubation (n=3).

Sample Standard Curve for Acetylcholinesterase

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