



Acetylcholine Assay Kit

Catalog Number KA1624

100 assays

Version: 04

Intended for research use only

www.abnova.com

Table of Contents

Introduction	3
Intended Use	3
Background	3
General Information	4
Materials Supplied	4
Storage Instruction	4
Materials Required but Not Supplied	4
Precautions for Use	4
Assay Protocol	5
Reagent Preparation	5
Sample Preparation	5
Assay Procedure	5
Data Analysis	6
Calculation of Results	6
Resources	7
References	7

Introduction

Intended Use

Applications:

- ✓ Assays: acetylcholine in biological samples such as serum, plasma, urine, saliva, milk, tissue, and cell culture.
- ✓ Drug Discovery/Pharmacology: effects of drugs on acetylcholine metabolism.

Features:

Use 20 μ L samples. Linear detection range: colorimetric assay 10 to 200 μ M, fluorimetric assay 0.4 to 10 μ M acetylcholine.

Background

ACETYLCHOLINE is a neurotransmitter produced in acetylcholinergic neurons. It plays important roles in skeletal muscle movement, regulation of smooth and cardiac muscles, as well as in learning, memory and mood. Acetylcholine Assay Kit provides a simple, direct and high-throughput assay for measuring acetylcholine in biological samples. In this assay, acetylcholine is hydrolyzed by acetylcholinesterase to choline which is oxidized by choline oxidase to betaine and H_2O_2 . The resulting H_2O_2 reacts with a specific dye to form a pink colored product. The color intensity at 570nm or fluorescence intensity (530/585 nm) is directly proportional to the acetylcholine concentration in the sample.

General Information

Materials Supplied

List of component

Component	Amount
Assay Buffer	10 mL
ACHE Enzyme	120 μ L
Enzyme Mix: Dried	1 vial
Dye Reagent	120 μ L
Standard: 2 mM acetylcholine	400 μ L

Storage Instruction

Store all components at -20°C. Shelf life of six months after receipt.

Materials Required but Not Supplied

- ✓ Pipetting devices
- ✓ Centrifuge tubes
- ✓ Clear flat-bottom uncoated 96-well plates
- ✓ Optical density plate reader
- ✓ Black flat-bottom uncoated 96-well plates
- ✓ Fluorescence plate reader.

Precautions for Use

Reagents are for research use only. Normal precautions for laboratory reagents should be exercised while using the reagents.

Assay Protocol

Reagent Preparation

Equilibrate all components to room temperature. Briefly centrifuge the tubes before opening. Keep thawed tubes on ice during assay. Keep thawed tubes on ice during assay.

- Reconstitute Enzyme mix with 120 μL Assay Buffer. Reconstituted Enzyme mix is stable for 1 month when stored at -20°C .
- Mix 24 μL 2 mM Standard with 216 μL dH_2O (final 200 μM). Dilute standard in dH_2O as follows.

No	200 μM STD + H_2O	Vol (μL)	Acetylcholine (μM)
1	100 μL + 0 μL	100	200
2	60 μL + 40 μL	100	120
3	30 μL + 70 μL	100	60
4	0 μL + 100 μL	100	0

Sample Preparation

- ✓ Liquid samples such as serum and plasma can be assayed directly. Tissue and cell lysates can be prepared by homogenization in cold 1 x PBS and centrifugation (5 min at 14,000 rpm). Use clear supernatants for assay. Milk samples should be cleared by mixing 600 μL milk with 100 μL 6 N HCl. Centrifuge 5 min at 14,000 rpm. Transfer 300 μL supernatant into a clean tube and neutralize with 50 μL 6 N NaOH. The neutralized supernatant is ready for assay (dilution factor $n = 1.36$).

*Note: (1). SH-containing reagents (e.g. *b*-mercaptoethanol, dithiothreitol, > 5 μM) are known to interfere in this assay and should be avoided in sample preparation. (2). This assay is based on an enzyme-catalyzed kinetic reaction. Addition of Working Reagent should be quick and mixing should be brief but thorough.*

Assay Procedure

- ✓ Colorimetric assay
- 1. Transfer 20 μL of diluted standards and samples into separate wells of a clear flatbottom 96-well plate.
Note: if a sample is known to contain choline, prepare an extra sample blank well with 20 μL of the sample.
- 2. Color reaction. Prepare enough Working Reagent by mixing, for each well, 85 μL Assay Buffer, 1 μL ACHE Enzyme, 1 μL Enzyme Mix and 1 μL Dye Reagent. Add 80 μL Working Reagent to each well. Immediately tap plate to mix. Incubate 30 min at room temperature.
Note: for samples that contain choline, prepare a blank control reagent with no ACHE Enzyme (i.e., 85 μL

Assay Buffer, 1 μL Enzyme Mix and 1 μL Dye Reagent). Add 80 μL of the control Reagent to each Sample Blank well.

3. Read optical density at 570 nm (550-585 nm).

✓ Fluorimetric assay

The fluorimetric assay procedure is similar to the colorimetric procedure except that (1) 0, 3, 6 and 10 μM acetylcholine standards and (2) a black 96-well plate are used. Read fluorescence intensity at $\lambda_{\text{ex}} = 530$ nm and $\lambda_{\text{em}} = 585$ nm.

Note: if the calculated acetylcholine concentration of a sample is higher than 200 μM in the Colorimetric Assay or 10 μM in the Fluorimetric Assay, dilute sample in water and repeat the assay. Multiply result by the dilution factor n.

Data Analysis

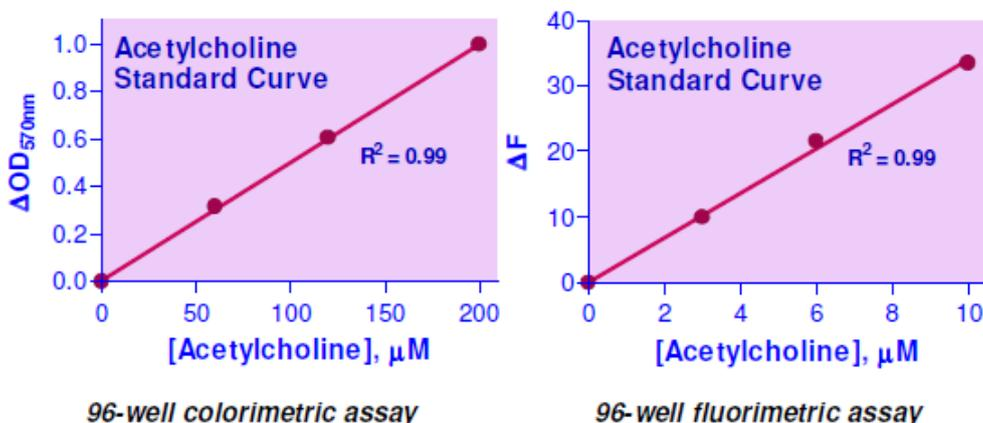
Calculation of Results

Subtract blank value (#4) from the standard values and plot the ΔOD or ΔF against standard concentrations. Determine the slope and calculate the acetylcholine concentration of Sample,

$$[\text{Acetylcholine}] = \frac{R_{\text{SAMPLE}} - R_{\text{BLANK}}}{\text{Slope}(\mu\text{M}^{-1})} \times n (\mu\text{M})$$

R_{SAMPLE} and R_{BLANK} are optical density or fluorescence intensity readings of the Sample and H₂O Blank (or Sample Blank if sample contains choline), respectively. n is the sample dilution factor.

Conversions: 1 mM acetylcholine equals 14.6 mg/dL, 0.015% or 146 ppm.



Resources

References

1. Vizi, E.S. et al (1985). A simple and sensitive method of acetylcholine identification and assay. Bioassay combined with minicolumn gel filtration or high-performance liquid chromatography. J Pharmacol Methods. 13:201-211.
2. Gilberstadt, M.L. and Russell, J.A. (1984). Determination of picomole quantities of acetylcholine and acetylcholine in physiologic salt solutions. Anal Biochem. 138:78-85.
3. Israel, M. and Lesbats, B. (1982). Application to mammalian tissues of the chemiluminescent method for detecting acetylcholine. J Neurochem. 39:248-250.