

PRODUCT INFORMATION



Acetylcholinesterase (eel)

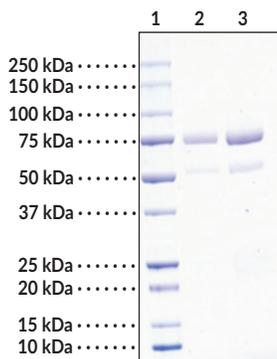
Item No. 44687

Overview and Properties

Synonym:	AChE
Source:	Native AChE purified from electric eel
Amino Acids:	1-633 (full length)
Uniprot No.:	O42275
Molecular Weight:	71.81 kDa
Storage:	-80°C (as supplied)
Stability:	≥1 year
Purity:	≥70% estimated by SDS-PAGE
Supplied in:	20 mM potassium phosphate, pH 7.4
Protein	
Concentration:	<i>batch specific</i> mg/ml
Activity:	<i>batch specific</i> U/ml
Specific Activity:	<i>batch specific</i> U/mg
Unit Definition:	One unit is defined as the amount of AChE required to produce 1 μmol of thiocholine per minute at room temperature in 1X PBS buffer, pH 7.4, containing 1 mM acetylthiocholine as a substrate.

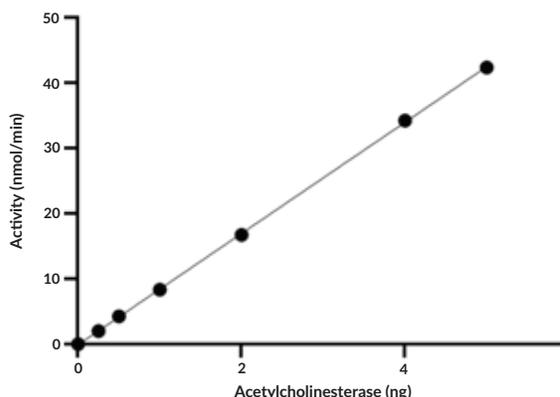
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: Acetylcholinesterase (2 μg)
Lane 3: Acetylcholinesterase (4 μg)

SDS-PAGE Analysis of Acetylcholinesterase.



Native acetylcholinesterase activity was measured using 1 mM acetylthiocholine (Item No. 34313) as a substrate in 1X PBS, pH 7.4 at room temperature.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/20/2025

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Acetylcholinesterase (AChE) is a highly conserved homotetrameric carboxylesterase that is composed of four ~70 kDa subunits, each containing a single active site.¹⁻³ In the electric eel (*E. electricus*), AChE is highly expressed in the electric organs at the posterior innervated membrane, where it hydrolyzes acetylcholine to acetate and choline to terminate synaptic transmission between electromotor neurons and electrocytes.⁴ AChE has commonly been used in biomedical research in assays, such as ELISA, for the detection or quantification of target molecules.⁵ It has also been used in the detection of organophosphate and carbamate pesticides due to their ability to inhibit AChE activity. Cayman's Acetylcholinesterase (eel) protein can be used for ELISA, enzyme activity assay, and Western blot applications.

References

1. Ha, Z.H., Matthew, S., and Yeong, K.Y. Butyrylcholinesterase: A multifaceted pharmacological target and tool. *Curr. Protein Pept. Sci.* **21(1)**, 99-109 (2020).
2. Cartaud, J., Rieger, F., Bon, S., *et al.* Fine structure of electric eel acetylcholinesterase. *Brain Res.* **88(1)**, 127-130 (1975).
3. Rosenberry, T.L., Chen, Y.T., and Bock, E. Structure of 11S acetylcholinesterase. Subunit composition. *Biochemistry* **13(15)**, 3068-79 (1974).
4. Gotter, A.L., Kaetzel, M.A., and Dedman, J.R. *Electrophorus electricus* as a model system for the study of membrane excitability. *Comp. Biochem. Physiol. A Mol. Integr. Physiol.* **119(1)**, 225-241 (1998).
5. Liu, D.-M., Xu, B., and Dong, C. Recent advances in colorimetric strategies for acetylcholinesterase assay and their applications. *Trends Analyt. Chem.* **142**, 116320 (2021).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM