

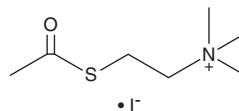
# PRODUCT INFORMATION



## Acetylthiocholine (iodide)

Item No. 34313

<b>CAS Registry No.:</b>	1866-15-5
<b>Formal Name:</b>	2-(acetylthio)-N,N,N-trimethyl-ethanaminium, monoiodide
<b>Synonyms:</b>	ASCh, ATCh
<b>MF:</b>	C <sub>7</sub> H <sub>16</sub> NOS • I
<b>FW:</b>	289.2
<b>Purity:</b>	≥95%
<b>Supplied as:</b>	A solid
<b>Storage:</b>	-20°C
<b>Stability:</b>	≥4 years



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

### Laboratory Procedures

Acetylthiocholine (iodide) is supplied as a solid. A stock solution may be made by dissolving the acetylthiocholine (iodide) in the solvent of choice, which should be purged with an inert gas. Acetylthiocholine (iodide) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of acetylthiocholine (iodide) in these solvents is approximately 30 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of acetylthiocholine (iodide) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of acetylthiocholine (iodide) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Acetylthiocholine is a substrate for acetylcholinesterase (AChE) and an agonist of  $\alpha 4\beta 2$  subunit-containing nicotinic acetylcholine receptors (nAChRs).<sup>1,2</sup> It binds to  $\alpha 4\beta 2$  subunit-containing nAChRs in SH-EP1 cells ( $IC_{50} = 7.1 \mu M$  for the human receptor) and induces rubidium efflux in an agonist-induced rubidium efflux assay ( $EC_{50} = 100 \mu M$ ).<sup>2</sup> Acetylthiocholine has been used as part of a thiocholine sensor for the detection of pesticides in complex sample types, such as soil, groundwater, and spinach.<sup>1</sup>

### References

1. Chen, L., Tian, X., Li, Y., *et al.* Broad-spectrum pesticide screening by multiple cholinesterases and thiocholine sensors assembled high-throughput optical array system. *J. Hazard. Mater.* **402**, 123830 (2021).
2. Eaton, J.B., Peng, J.-H., Schroeder, K.M., *et al.* Characterization of human  $\alpha 4\beta 2$ -nicotinic acetylcholine receptors stably and heterologously expressed in native nicotinic receptor-null SH-EP1 human epithelial cells. *Mol. Pharmacol.* **64**(6), 1283-1294 (2003).

#### 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, 10/17/2022

#### 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