

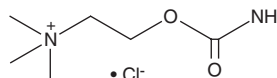
# PRODUCT INFORMATION



## Carbamoylcholine (chloride)

Item No. 14486

**CAS Registry No.:** 51-83-2  
**Formal Name:** 2-[(aminocarbonyl)oxy]-N,N,N-trimethyl-ethanaminium, monochloride  
**Synonyms:** Carbachol  
**MF:** C<sub>6</sub>H<sub>15</sub>N<sub>2</sub>O<sub>2</sub> • Cl  
**FW:** 182.7  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Carbamoylcholine (chloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the carbamoylcholine (chloride) in the solvent of choice, which should be purged with an inert gas. Carbamoylcholine (chloride) is soluble in the organic solvent DMSO at a concentration of approximately 0.1 mg/ml 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 carbamoylcholine (chloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of carbamoylcholine (chloride) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Carbamoylcholine, also known as carbachol, is an analog of acetylcholine that activates acetylcholine receptors (AChR). Carbamoylcholine is an agonist of both nicotinic (nAChR) and muscarinic (mAChR) receptors, with reported K<sub>i</sub> values ranging from 10 to 10,000 nM for different receptors and different preparations.<sup>1-5</sup> Carbamoylcholine (chloride) is used to study responses mediated by nAChR and mAChR, including smooth muscle contraction, gut motility, and neuronal signaling.<sup>6-8</sup>

### References

1. Dávila-García, M.I., Musachio, J.L., Perry, D.C., et al. *J. Pharmacol. Exp. Ther.* **282(1)**, 445-451 (1997).
2. Eaton, J.B., Peng, J.-H., Schroeder, K.M., et al. *Mol. Pharmacol.* **64(6)**, 1283-1294 (2003).
3. Grinevich, V.P., Letchworth, S.R., Lindenberger, K.A., et al. *J. Pharmacol. Exp. Ther.* **312(2)**, 619-626 (2005).
4. Jensen, A.A., Mikkelsen, I., Frølund, B., et al. *Mol. Pharmacol.* **64(4)**, 865-875 (2003).
5. Xiao, Y. and Kellar, K.J. *J. Pharmacol. Exp. Ther.* **310(1)**, 98-107 (2004).
6. Aldini, R., Micucci, M., Cevenini, M., et al. *PLoS One* **9(9)**, 1-16 (2014).
7. Lee, H.S., Ghetti, A., Pinto-Duarte, A., et al. *Proc. Natl. Acad. Sci. USA* **111(32)**, E3343-E3352 (2014).
8. Mori, D., Watanabe, N., Kaminuma, O., et al. *J. Pharmacol. Sci.* **125(4)**, 394-405 (2014).

#### 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, 11/10/2022

#### 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