

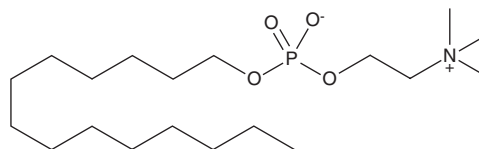
# PRODUCT INFORMATION



## Tetradecylphosphocholine

Item No. 25700

**CAS Registry No.:** 77733-28-9  
**Formal Name:** 2-[[hydroxy(tetradecyloxy)phosphinyl]oxy]-N,N,N-trimethylethanaminium, inner salt  
**Synonym:** Fos-choline 14  
**MF:**  $C_{19}H_{42}NO_4P$   
**FW:** 379.5  
**Purity:**  $\geq 95\%$   
**Supplied as:** A crystalline solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 4$  years



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

### Laboratory Procedures

Tetradecylphosphocholine is supplied as a crystalline solid. A stock solution may be made by dissolving the tetradecylphosphocholine in the solvent of choice. Tetradecylphosphocholine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of tetradecylphosphocholine in ethanol is approximately 15 mg/ml and approximately 16 mg/ml in DMSO and DMF.

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 tetradecylphosphocholine can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of tetradecylphosphocholine in PBS, pH 7.2, is approximately 25 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Tetradecylphosphocholine is a zwitterionic surfactant. It has been used in the solubilization and purification of G protein-coupled receptors (GPCRs).<sup>1,2</sup> It has also been used as the stationary phase in ion chromatography for the determination of inorganic acids.<sup>3</sup> Tetradecylphosphocholine has a critical micelle concentration (CMC) of approximately 0.12 mM.

### References

1. Cook, B.L., Ernberg, K.E., Chung, H., *et al.* Study of a synthetic human olfactory receptor 17-4: Expression and purification from an inducible mammalian cell line. *PLoS One* **3(8)**, e2920 (2008).
2. Ren, H., Yu, D., Ge, B., *et al.* High-level production, solubilization and purification of synthetic human GPCR chemokine receptors CCR5, CCR3, CXCR4, and CX3CR1. *PLoS One* **4(2)**, e4509 (2009).
3. Hu, W., Hasebe, K., Tanaka, K., *et al.* Determination of inorganic acids by ion chromatography with n-tetradecylphosphocholine (zwitterionic surfactant) as the stationary phase and pure water as the mobile phase. *Fresenius J. Anal. Chem.* **370(4)**, 399-402 (2001).

#### 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, 12/19/2022

#### 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](http://WWW.CAYMANCHEM.COM)