

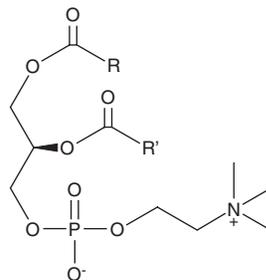
# PRODUCT INFORMATION



## Phosphatidylcholine (soy)

Item No. 34800

**CAS Registry No.:** 97281-47-5  
**Synonyms:** E 322, Lecithin (soy), Lecithinon, Phospholutein, PtdCho, 3-fluoro UB  
**MF:** C<sub>44</sub>H<sub>80</sub>NO<sub>8</sub>P (for linoleoyl)  
**FW:** 782.1  
**Purity:** ≥95% (mixture)  
**Supplied as:** A 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

Phosphatidylcholine (soy) is supplied as a solid. A stock solution may be made by dissolving the phosphatidylcholine (soy) in the solvent of choice, which should be purged with an inert gas. Phosphatidylcholine (soy) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of phosphatidylcholine (soy) in these solvents is approximately 10 mg/ml.

### Description

Phosphatidylcholine is the most abundant phospholipid in mammalian, plant, and yeast cells.<sup>1</sup> It is found mainly in the outer leaflet of cell membranes and can make up approximately half of the total phospholipids.<sup>2</sup> In mammalian tissues, phosphatidylcholine commonly contains a saturated and unsaturated fatty acid at the C-1 and C-2 positions of glycerol, respectively. It is a substrate for various enzymes in cell signaling pathways that is cleaved by phospholipases into diacylglycerol and phosphocholine or phosphatidic acid and choline. This product contains phosphatidylcholine molecular species with variable fatty acyl chain lengths at the *sn*-1 and *sn*-2 positions.

### References

1. Vance, J.E. Phosphatidylserine and phosphatidylethanolamine in mammalian cells: Two metabolically related aminophospholipids. *J. Lipid Res.* **49(7)**, 1377-1387 (2008).
2. Billah, M.M. and Anthes, J.C. The regulation and cellular functions of phosphatidylcholine hydrolysis. *Biochem. J.* **269(2)**, 281-291 (1990).

#### 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, 01/26/2024

#### 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