

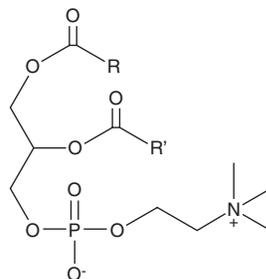
PRODUCT INFORMATION



Phosphatidylcholine (chicken egg)

Item No. 24343

CAS Registry No.:	8002-43-5
Formal Name:	Lecithins
Synonyms:	E 322, Lecithin (egg), Lecithinon, Phospholutein, PtdCho, 3-fluoro UB
MF:	C ₄₄ H ₈₄ NO ₈ P (for oleoyl)
FW:	786.1
Purity:	≥98%
Supplied as:	A solution in chloroform
Storage:	-20°C
Stability:	≥2 years



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

Laboratory Procedures

Phosphatidylcholine (chicken egg) is supplied as a solution in chloroform. To change the solvent, simply evaporate the chloroform under a gentle stream of nitrogen and immediately add the solvent of choice. Phosphatidylcholine (chicken egg) is soluble in the organic solvent ethanol.

Description

Phosphatidylcholine is the most abundant phospholipid in mammalian, plant, and yeast cells.¹ It is found mainly in the outer leaflet of cell membranes and can make up approximately half of the total phospholipids.² 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 glucosylceramide molecular species with primarily C16:0, C22:0, and C24:0 fatty acyl chain lengths. As this product is derived from a natural source, there may be variations in the sphingoid backbone.

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. *Biochemistry Journal* **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, 08/24/2022

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