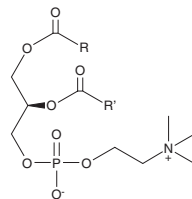


PRODUCT INFORMATION

Phosphatidylcholine (hydrogenated) (soy)

Item No. 45786

CAS Registry No.: 97281-48-6
Synonyms: HSPC, Hydro Soy PC
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 (hydrogenated) (soy) (HSPC) is supplied as a solid. A stock solution may be made by dissolving the HSPC in the solvent of choice, which should be purged with an inert gas. HSPC is slightly soluble (0.1-1 mg/ml) in chloroform.

Description

HSPC is a hydrogenated glycerophospholipid derived from natural soy lecithin.¹ HSPC has been widely used in the generation of liposomes.¹⁻³ HSPC (1.25 wt% in the diet) decreases hepatic total lipid and cholesterol levels in mice fed a high-fat diet.⁴ Formulations containing HSPC have been used in cosmetic products as skin-conditioning agents and in the generation of liposomes for the delivery of anticancer and antifungal agents. This product contains phosphatidylcholine molecular species with primarily C16:0 and C18:0 fatty acyl chains.

References

1. Rao, P., Rudrakshula, M., Potham, R., *et al.* Unified synthesis of DSPC and PSPC: Chemical entities of hydrogenated soy L- α -phosphatidylcholine (HSPC), a key component of liposomal drug formulations. *Tetrahedron Lett.* **155**, 155424 (2025).
2. Sha, X., Guo, J., Chen, Y., *et al.* Effect of phospholipid composition on pharmacokinetics and biodistribution of epirubicin liposomes. *J. Liposome Res.* **22**(1), 80-88 (2012).
3. Son, K., Ueda, M., Taguchi, K., *et al.* Evasion of the accelerated blood clearance phenomenon by polysarcosine coating of liposomes. *J. Control Release* **322**, 209-216 (2020).
4. Tandy, S., Chung, R.W., Kamili, A., *et al.* Hydrogenated phosphatidylcholine supplementation reduces hepatic lipid levels in mice fed a high-fat diet. *Atherosclerosis* **213**(1), 142-147 (2010).

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, 05/22/2026

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