

PRODUCT INFORMATION



10-Pyrene-PC

Item No. 62245

CAS Registry No.: 95864-17-8
Formal Name: 1-O-hexadecanoyl-2-O-(10-(1-pyrenyl) decanoyl)-sn-glycerol-3-phosphorylcholine
Synonym: 1-Palmitoyl-2-pyrenedecanoyl Phosphatidylcholine

MF: C₅₀H₇₆NO₈P

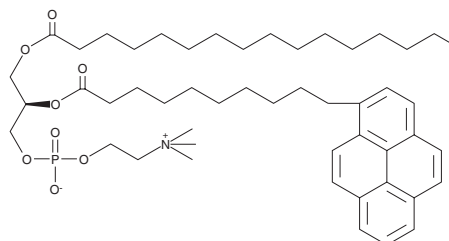
FW: 850.1

Purity: ≥98%

Supplied as: A solution in chloroform

Storage: -20°C

Stability: ≥1 year



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

Laboratory Procedures

10-Pyrene-PC is supplied as a solution in chloroform. The solubility of 10-pyrene-PC in water is approximately 1 mg/ml. In water, 10-pyrene-PC phospholipids form micelles. Avoid dissolving 10-pyrene-PC in basic solutions, since base treatment promotes hydrolysis of the thioester.

Description

10-Pyrene-PC is a substrate for all PLA₂s with the exception of cPLA₂ and PAF-AH.¹ Upon phospholipid hydrolysis, 10-pyrenyldecanoic acid is produced from 10-pyrene-PC. The monomeric 10-pyrenyldecanoic acid exhibits fluorescence (excitation 345 nm, emission 395 nm), allowing quantitation of phospholipase activity.²

References

1. Roberts, M.F. Phospholipases: Structural and functional motifs for working at an interface. *FASEB J.* 10(10), 1159-1172 (1996).
2. Lusa, S., Myllärniemi, M., Volmonen, K., *et al.* Degradation of pyrene-labelled phospholipids by lysosomal phospholipases in vitro. Dependence of degradation on the length and position of the labelled and unlabelled acyl chains. *Biochem. J.* 315(Pt 3), 947-952 (1996).

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, 11/11/2021

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