

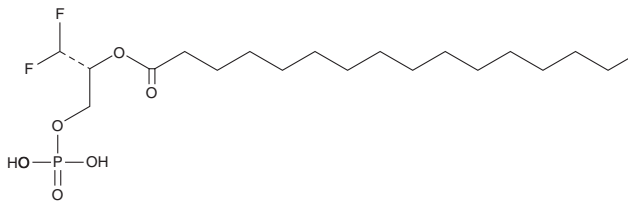
PRODUCT INFORMATION



XY-4

Item No. 39273

CAS Registry No.: 474329-47-0
Formal Name: hexadecanoic acid, (1R)-2,2-difluoro-1-[(phosphonoxy)methyl]ethyl ester
MF: C₁₉H₃₇F₂O₆P
FW: 430.5
Purity: ≥95%
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

XY-4 is supplied as a solid. A stock solution may be made by dissolving the XY-4 in the solvent of choice, which should be purged with an inert gas. XY-4 is soluble in methanol.

Description

XY-4 is a derivative of 1-palmitoyl-2-hydroxy-*sn*-glycero-3-PA (1-palmitoyl LPA; Item No. 10010094) and an agonist of peroxisome proliferator-activated receptor γ (PPAR γ).¹ It induces reporter gene expression in RAW 264.7 macrophages expressing a peroxisome proliferator-response element (PPRE) when used at a concentration of 5 μ M. XY-4 does not act as an agonist at lysophosphatidic acid receptor 1 (LPA₁), LPA₂, or LPA₃. It induces platelet aggregation *in vitro* when used at a concentration of 1 μ M.² XY-4 induces carotid artery neointima formation in rats.³

References

- McIntyre, T.M., Pontsler, A.V., Silva, A.R., *et al.* Identification of an intracellular receptor for lysophosphatidic acid (LPA): LPA is a transcellular PPAR γ agonist. *Proc. Natl. Acad. Sci. USA* **100**(1), 131-136 (2003).
- Xu, Y., Qian, L., and Prestwich, G.D. Synthesis of monofluorinated analogues of lysophosphatidic acid. *J. Org. Chem.* **68**(13), 5320-5330 (2003).
- Zhang, C., Baker, D.L., Yasuda, S., *et al.* Lysophosphatidic acid induces neointima formation through PPAR γ activation. *J. Exp. Med.* **199**(6), 763-774 (2004).

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, 01/05/2024

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