

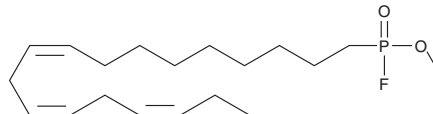
Product Information



Methyl α -Linolenyl Fluorophosphonate

Item No. 70662

Formal Name: 9Z,12Z,15Z-octadecatrienyl-phosphonofluoridic acid, methyl ester
Synonym: MLnFP
MF: C₁₉H₃₄FO₂P
FW: 344.4
Purity: \geq 98%
Stability: \geq 1 year at -80°C
Supplied as: A solution in methyl acetate



Laboratory Procedures

For long term storage, we suggest that methyl α -linolenyl fluorophosphonate (MLnFP) be stored as supplied at -80°C. It should be stable for at least one year.

MLnFP is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of MLnFP in these solvents is approximately 5 mg/ml.

Further dilutions of the organic solvent solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Store aqueous solutions of MLnFP on ice and use within 12 hours of preparation. Although the aqueous solutions of MLnFP may be stable for more than 12 hours, we strongly recommend using a fresh preparation each day.

MLnFP is an analog of methyl arachidonyl fluorophosphonate (MAFP; Item No. 70660), which has been widely studied as an inhibitor of phospholipases, FAAH, and as a CB receptor ligand.¹⁻³ The pharmacology of the α -linolenyl analog of MAFP has not been completely investigated.

References

- Huang, Z., Liu, S., Street, I., *et al.* Methyl arachidonyl fluorophosphonate, a potent irreversible cPLA₂ inhibitor, blocks the mobilization of arachidonic acid in human platelets and neutrophils. *Mediators Inflamm.* **3**, 307-308 (1994).
- Deutsch, D.G., Omeir, R., Arreaza, G., *et al.* Methyl arachidonyl fluorophosphonate: A potent irreversible inhibitor of anandamide amidase. *Biochem. Pharmacol.* **53**, 255-260 (1997).
- Lio, Y.-C., Reynolds, L.J., Balsinde, J., *et al.* Irreversible inhibition of Ca²⁺-independent phospholipase A₂ by methyl arachidonyl fluorophosphonate. *Biochim. Biophys. Acta* **1302**, 55-60 (1996).

Related Products

URB597 - Item No. 10046 • URB937 - Item No. 10674 • Methyl Arachidonyl Fluorophosphonate - Item No. 70660 • Methyl γ -Linolenyl Fluorophosphonate - Item No. 70664 • Methyl Nonadecadienyl Fluorophosphonate - Item No. 70667 • α -Linolenic Acid - Item No. 90210

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any **obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 03/09/2011