

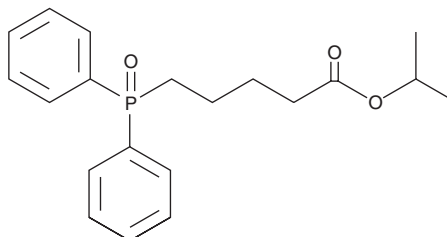
PRODUCT INFORMATION



Isopropyl 5-(Diphenylphosphoryl)pentanoate

Item No. 10010978

CAS Registry No.: 2088449-88-9
Formal Name: 5-(diphenylphosphoryl)-pentanoic acid, 1-methylethyl ester
Synonym: DPP isopropyl ester
MF: C₂₀H₂₅O₃P
FW: 344.4
Purity: ≥98%
UV/Vis.: λ_{max}: 222, 266, 272 nm
Supplied as: A crystalline 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

Isopropyl 5-(diphenylphosphoryl)pentanoate is supplied as a crystalline solid. A stock solution may be made by dissolving the isopropyl 5-(diphenylphosphoryl)pentanoate in the solvent of choice, which should be purged with an inert gas. Isopropyl 5-(diphenylphosphoryl)pentanoate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of isopropyl 5-(diphenylphosphoryl)pentanoate in ethanol is approximately 50 mg/ml and approximately 25 mg/ml in DMSO and DMF.

Isopropyl 5-(diphenylphosphoryl)pentanoate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, isopropyl 5-(diphenylphosphoryl)pentanoate should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Isopropyl 5-(diphenylphosphoryl)pentanoate has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Latanoprost is the isopropyl ester of 17-phenyl-13,14-dihydro prostaglandin F2α. It is a prodrug of the free acid, which is a potent agonist of the FP receptor in the eye. Isopropyl 5-(diphenylphosphoryl)pentanoate is a potential trace impurity in commercial preparations of latanoprost.

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, 01/31/2024

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