

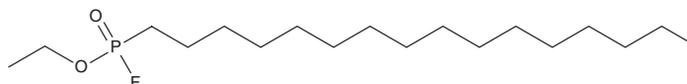
PRODUCT INFORMATION



Ethyl HDFP

Item No. 38209

CAS Registry No.: 1407160-81-9
Formal Name: p-hexadecyl-phosphonofluoridic acid, ethyl ester
Synonyms: Ethyl Hexadecylfluorophosphonate, Ethyl P-Hexadecylphosphonofluoridate
MF: C₁₈H₃₈FO₂P
FW: 336.5
Purity: ≥98%
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

Ethyl HDFP is supplied as a solid. A stock solution may be made by dissolving the Ethyl HDFP in the solvent of choice, which should be purged with an inert gas. Ethyl HDFP is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of Ethyl HDFP in DMSO is approximately 20 mg/ml and approximately 30 mg/ml in ethanol and DMF.

Description

Ethyl HDFP is a lipase inhibitor.¹ It inhibits 12 lipases, including fatty acid amide hydrolase (FAAH), fatty acid synthase (FASN), and α/β -hydrolase domain-containing protein 4 (ABHD4), by greater than 75% in mouse T cell hybridoma lysates when used at a concentration of 2 μ M. Ethyl HDFP inhibits depalmitoylation of H-Ras in HEK293 cells.

Reference

1. Martin, B.R., Wang, C., Adibekian, A., *et al.* Global profiling of dynamic protein palmitoylation. *Nat. Methods* **9**(1), 84-89 (2012).

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, 10/20/2025

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