

Methyl Arachidonyl Fluorophosphonate

Phospholipase inhibitor

Selective, active-site directed, irreversible inhibitor of cPLA₂ (IC₅₀=5μM) and Ca²⁺-independent iPLA₂ (IC₅₀=0.5μM). Potent, irreversible inhibitor of fatty acid amide hydrolase (FAAH) (IC₅₀=2.5nM), 800 -fold more potent than AACOCF₃ (Prod. No. BML-ST335). Irreversible CB₁ receptor antagonist in rat brain membrane (IC₅₀=20nM).

Citations: 4

[View Online »](#)

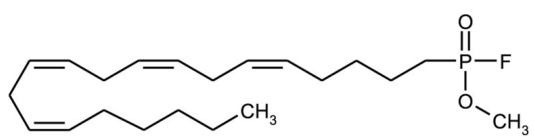
Ordering Information

[Order Online »](#)

BML-ST360-0005	5mg
BML-ST360-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at -80°C. Stock solutions are stable for up to 2 months at -20°C.
Handling	Protect from air.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	MAFP, Methylphosphonofluoridic acid 5,8,11,14-eicosatetraenyl ester
Appearance	Colorless to yellow oil.
CAS	188404-10-6
Couple Target	Cannabinoid receptor, FAAH, Phospholipase, PLA
Couple Type	Inhibitor
Formula	$C_{21}H_{36}FO_2P$
MW	370.5
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol (50mg/ml) or DMSO (50mg/ml).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)