

AACOCF₃

Inhibitor of PLA₂ and FAAH

AACOCF₃ is a trifluoromethylketone derivative of arachidonic acid. It is a selective inhibitor of human cytosolic PLA₂ and Ca²⁺-independent PLA₂ (iPLA₂, IC₅₀=15 µM). It is active in intact platelets and U937 cells (IC₅₀=8 µM). It also inhibits fatty acid amide hydrolase (100% at 7.5 µM).

Citations: 11

[View Online »](#)

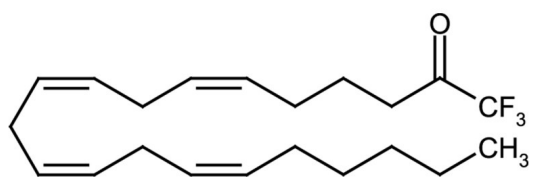
Ordering Information

[Order Online »](#)

BML-ST335-0010	10mg
BML-ST335-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions are stable for up to 1 month when stored under nitrogen or argon in tightly sealed container at -80°C.
Handling	Protect from air. Dissolve in solvent purged with nitrogen or argon. Store under nitrogen or argon in tightly sealed container.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance	Yellow oil.
CAS	149301-79-1
Couple Target	FAAH, Phospholipase, PLA
Couple Type	Inhibitor
Formula	$C_{21}H_{31}F_3O$
Identity	Determined by NMR.
MW	356.5
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (>25mg/ml) or 100% ethanol (>25mg/ml).

Last modified: May 29, 2024



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-
fr@enzolifesciences.com](mailto:info-fr@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)