

PACOFC3

Phospholipase inhibitor

Inhibitor of both Ca^{2+} -dependent cytosolic cPLA_2 ($\text{IC}_{50}=45 \mu\text{M}$) and Ca^{2+} -independent phospholipases A_2 iPLA_2 ($\text{IC}_{50}=3.8 \mu\text{M}$).

Citations: 2

[View Online »](#)

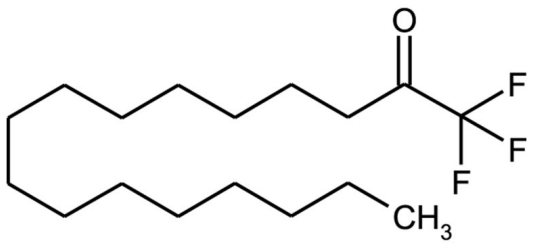
Ordering Information

[Order Online »](#)

BML-ST336-0010	10mg
----------------	------

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 -20°C. Stock solutions are stable for up to 3 months at -20°C. Not stable in basic solutions (pH > 7.4). We do not recommend storing aqueous solutions for more than one day.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Palmitoyl trifluoromethylketone, PTK
Appearance	White to off-white solid.
CAS	141022-99-3
Couple Target	Phospholipase
Couple Type	Inhibitor
Formula	$C_{17}H_{31}F_3O$
MW	308.4
Purity	≥98% (TLC)
Solubility	Soluble in 100% ethanol, dimethyl formamide or DMSO (warm); sparingly soluble in neutral buffers.
Technical Info / Product Notes	Note: Product is not sterile.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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