

# PACOCF3

## Phospholipase inhibitor

Inhibitor of both  $\text{Ca}^{2+}$ -dependent cytosolic  $\text{cPLA}_2$  ( $\text{IC}_{50}=45\text{ }\mu\text{M}$ ) and  $\text{Ca}^{2+}$ -independent phospholipases  $\text{A}_2$   $\text{iPLA}_2$  ( $\text{IC}_{50}=3.8\text{ }\mu\text{M}$ ).

Citations: 2

[View Online »](#)

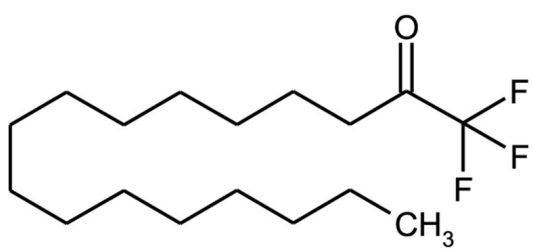
## Ordering Information

[Order Online »](#)

BML-ST336-0010	10mg
BML-ST336-0050	50mg

## 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	Palmityl 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](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)