

# N-(p-Amylcinnamoyl)anthranilic acid

## Inhibitor of PLA<sub>2</sub> and TRP channels

ACA is a close structural analog of the PLA<sub>2</sub> inhibitor, ONO-RS-082 and possesses similar but less potent inhibitory activity. Inhibits TRP channels independently of PLA<sub>2</sub>, including TRPM2, TRPC6 and TRPM8. It is a useful tool for studying the physiological role of TRP channels.

Citations: 6

[View Online »](#)

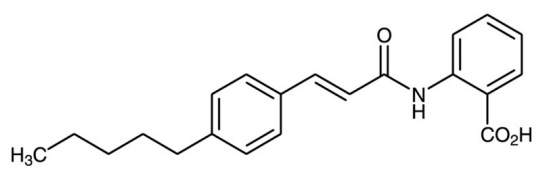
## Ordering Information

[Order Online »](#)

BML-EI178-0250	250mg
----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to 1 year. Make solutions fresh daily.
Long Term Storage	Ambient
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	ACA
Appearance	White solid.
CAS	110683-10-8
Couple Target	Phospholipase, TRP channel
Couple Type	Blocker, Inhibitor
Formula	$C_{21}H_{23}NO_3$
MW	337.4
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO (50 mg/ml) or ethanol (50 mg/ml @ 60°C).

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)