

N-(p-Amylcinnamoyl)anthranilic acid

Inhibitor of PLA₂ and TRP channels

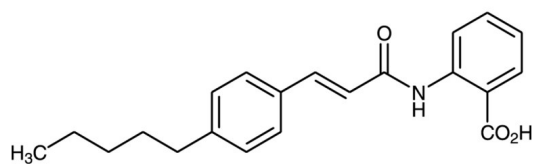
ACA is a close structural analog of the PLA₂ inhibitor, ONO-RS-082 and possesses similar but less potent inhibitory activity. Inhibits TRP channels independently of PLA₂, including TRPM2, TRPC6 and TRPM8. It is a useful tool for studying the physiological role of TRP channels.

Citations: 6

[View Online »](#)

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

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