

Ingenol 3,20-dibenzoate

PKC activator

Ingenol is a diterpenoid related to phorbol. Activates protein kinase C (PKC) and competes with [³H]phorbol dibutyrate for binding to PKCα (K_i for binding=240 nM). Induces apoptosis. Anticancer agent. Shows potent antileukemic activity in the P388 mouse assay. Displays unusually high mouse ear inflammatory activity, even though the 20-hydroxy group is blocked with a benzoate ester. It appears likely that this high potency is due to metabolic cleavage of the 20-benzoate group, which releases the more potent ingenol 3-benzoate.

Citations: 7

[View Online »](#)

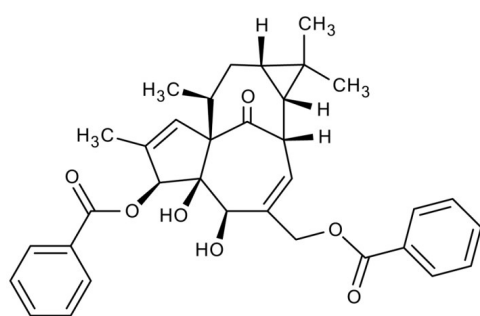
Ordering Information

[Order Online »](#)

BML-PE186-0001	1mg
----------------	-----

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, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	IDB
Appearance	White solid.
CAS	59086-90-7
Couple Target	PKC
Couple Type	Activator
Formula	$C_{34}H_{36}O_7$
MW	556.7
Purity	≥97%
Solubility	Soluble in DMSO (20mg/ml) or ethanol (20mg/ml).

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