

12-Deoxyphorbol 13-acetate

PKC activator

A non-tumor promoting phorbol ester, which acts as a PKC partial agonist. Protects T-lymphoblastoid cells from the killing effects of HIV-1. Upregulates latent HIV-1 provirus expression and inhibits HIV infection and viral spread at the entry/fusion step of viral life cycle. Does not induce tumor formation.

Citations: 7

[View Online »](#)

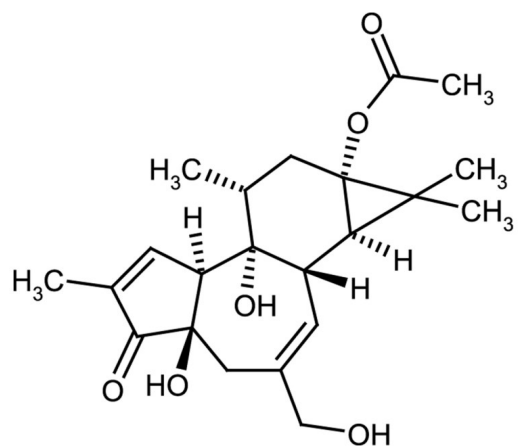
Ordering Information

[Order Online »](#)

BML-PE187-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Prostratin
Appearance	White to off white crystalline powder.
CAS	60857-08-1
Couple Target	PKC
Couple Type	Activator
Formula	$C_{22}H_{30}O_6$
MW	390.5
Purity	≥99% (HPLC, TLC)
Solubility	Soluble in 100% ethanol (25mg/ml) or DMSO (25mg/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