

Phorbol 12,13-dibutyrate

PKC activator

Widely used PKC activator (10-100ng/ml). Less potent than phorbol 12-myristate 13-acetate (PMA; TPA) (Prod. No. BML-PE160), but also less hydrophobic.

Citations: 12

[View Online »](#)

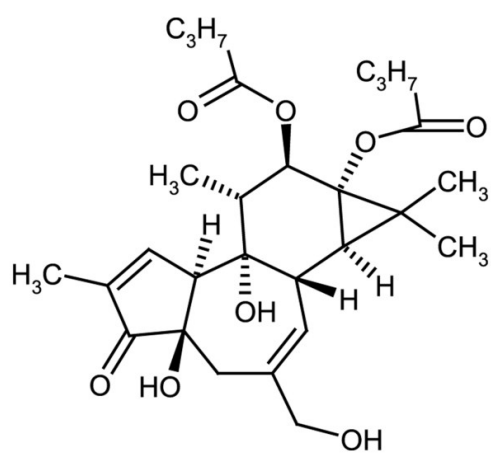
Ordering Information

[Order Online »](#)

BML-PE135-0005	5mg
----------------	-----

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.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Colorless oil.
CAS	37558-16-0
Couple Target	PKC
Couple Type	Activator
Formula	$C_{28}H_{40}O_8$
MW	504.6
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO (25mg/ml) or 100% ethanol (20mg/ml).

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)