

Phorbol 12-myristate 13-acetate

PKC activator

PKC activator (10-100 ng/ml). Most commonly used phorbol ester. Causes an extremely wide range of effects in cells and tissues. Very potent mouse skin tumor promoter. Inhibits apoptosis induced by Fas, but induces apoptosis in HL-60 promyelocytic leukemia cells.

Citations: 61

[View Online »](#)

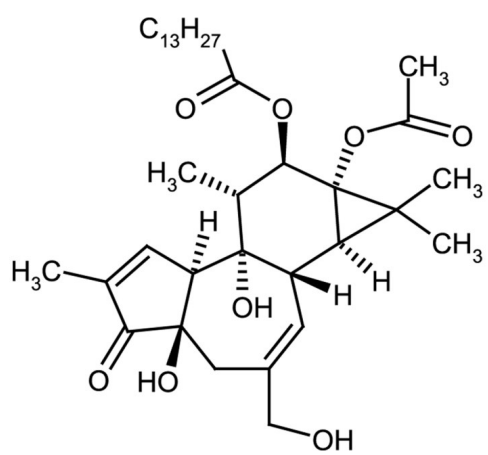
Ordering Information

[Order Online »](#)

| | |
|----------------|-----|
| BML-PE160-0001 | 1mg |
| BML-PE160-0005 | 5mg |

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

| | |
|-------------------|--|
| Use/Stability | As indicated on product label or CoA when stored as recommended. Dissolve in an organic solvent prior to dilution into aqueous buffers. Add calculated volume of organic solvent to vial to prepare concentrated stock solution. Intermediate dilutions into the same organic solvent are recommended prior to final dilution into aqueous buffer. 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

| | |
|------------------|---|
| Alternative Name | PMA, TPA |
| Appearance | Colorless amorphous film or foam. |
| CAS | 16561-29-8 |
| Couple Target | PKC |
| Couple Type | Activator |
| Formula | $C_{36}H_{56}O_8$ |
| MW | 616.8 |
| Purity | ≥98% (HPLC) |
| Solubility | Soluble in DMSO (25mg/ml), 100% ethanol (25mg/ml), acetone, ether or dimethyl formamide; almost insoluble in aqueous buffers. |

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-
be@enzolifesciences.com](mailto:info-be@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)