

HBDDE

PKC inhibitor

Inhibits PKC α (IC₅₀=43 μ M) and PKC γ (IC₅₀=50 μ M) without inhibiting the δ , β I and β II isotypes.

Citations: 1

[View Online »](#)

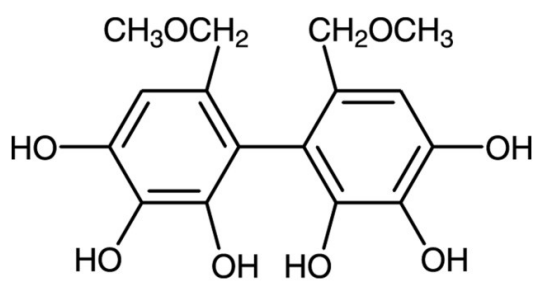
Ordering Information

[Order Online »](#)

BML-EI273-0005	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	Product is subject to oxidation. Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Orange solid.
CAS	154675-18-0
Couple Target	PKC
Couple Type	Inhibitor
Formula	$C_{16}H_{18}O_8$
MW	338.3
MeltingPoint	165-168°C
Purity	≥95% (TLC)
Solubility	Soluble in DMSO (25 mg/ml), ethanol (15 mg/ml warm) or water (10 mg/ml warm).

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)