

Calphostin C

PKC inhibitor

Potent and selective inhibitor of the protein kinase C (PKC). Inhibition is via interaction with the regulatory DAG binding site and phorbol ester binding site. Inhibition of PKC has been found to be light dependent. At higher concentrations inhibits myosin light chain kinase, cAMP-dependent protein kinase, protein kinase G, pp60v-src protein tyrosine kinase and DAG kinase. It also inhibits Phospholipase D1 and D2. Induces apoptotic DNA fragmentation and cell death. Kills breast cancer cells.

Has antiviral potential. Inhibits cardiac L-type Ca^{2+} channels. Cell permeable. Inhibits cell proliferation of malignant glioma cells in light-treated conditions *in vitro*.

Citations: 18

[View Online »](#)

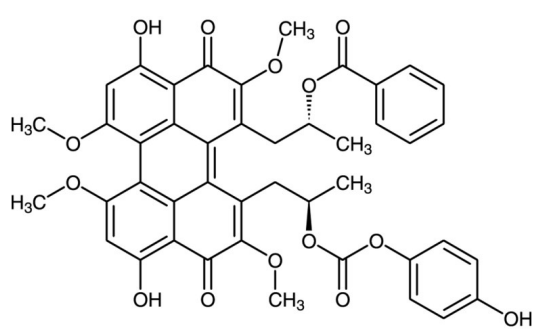
Ordering Information

[Order Online »](#)

BML-EI198-0100	100µg
----------------	-------

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 0-4°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	UCN-1028C
Appearance	Dark red solid.
CAS	121263-19-2
Couple Target	Calcium channel, Diacylglycerol kinase, MLCK, Phospholipase, PKA, PKC, PKG, PLD, Src kinase
Couple Type	Blocker, Inhibitor
Formula	$C_{44}H_{38}O_{14}$
MW	790.8
Purity	≥95% (HPLC)
Solubility	Soluble in ethanol, methanol, DMF or DMSO. Poor water solubility.



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)