

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 »](#)

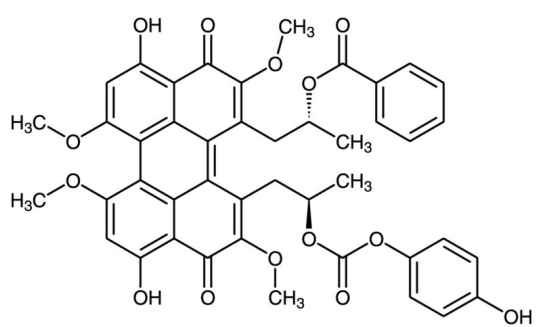
Ordering Information

[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. |

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-fr@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