

Phloretin

PKC inhibitor

Flavonoid. Protein kinase C (PKC) inhibitor. PGF_{2a} receptor antagonist. In astrocytes, these receptors are linked to phospholipase C, thus affecting phosphoinositide hydrolysis and intracellular Ca^{2+} levels. Inhibits *myo*-inositol uptake and 5'-iodothyronine deiodinase. Prevents TNF- α stimulated upregulation of VCAM-1, ICAM-1 and E-selectin. Enhances adiponectin expression.

Citations: 14

[View Online »](#)

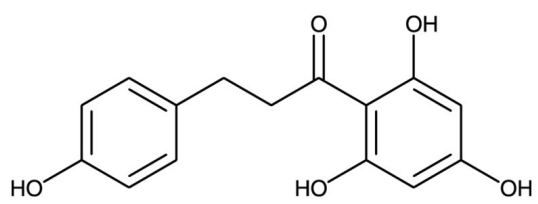
Ordering Information

[Order Online »](#)

BML-EI154-0200	200mg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Product may air oxidize in solution. Make solutions fresh daily.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Off-white to beige solid.
CAS	60-82-2
Couple Target	PKC, Prostaglandin receptor
Couple Type	Inhibitor
Formula	$C_{15}H_{14}O_5$
Identity	Determined by 1H -NMR and MS.
MI	14: 7326
MW	274.3
Purity	≥98% (TLC)
Solubility	Soluble in acetone, DMSO, 100% ethanol or methanol.

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)