

# Phloretin

## PKC inhibitor

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

Citations: 14

[View Online »](#)

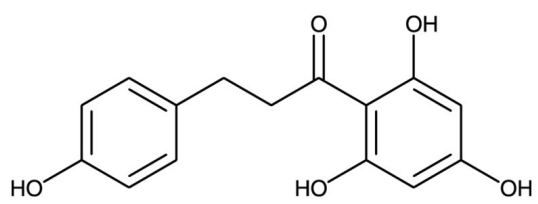
## Ordering Information

[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	<b>14:</b> 7326
MW	274.3
Purity	≥98% (TLC)
Solubility	Soluble in acetone, DMSO, 100% ethanol or methanol.



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)