

# Baicalein

## Lipoxygenase inhibitor

Inhibits cellular  $\text{Ca}^{2+}$  uptake and calcium mobilization. Inhibitor of protein tyrosine kinase in leukemia (CEM) cells. Induces cell cycle arrest and apoptosis. Anti-inflammatory compound. Has anti-thrombotic, anti-proliferative and anti-mitogenic effects. Originally reported as a selective inhibitor of 12-lipoxygenase, baicalein has since been shown to inhibit both 12-LO and 15-LO ( $\text{IC}_{50}$ =0.64 $\mu\text{M}$  for 12-LO and 1.6 $\mu\text{M}$  for 15-LO). Inhibits leukotriene biosynthesis (at higher concentrations) and release of lysosomal enzymes. Inhibits adjuvant-induced arthritis. Also inhibits prolyl oligopeptidase (POP/PREP) ( $\text{IC}_{50}$ =36 $\mu\text{M}$ ).

Citations: 21

[View Online »](#)

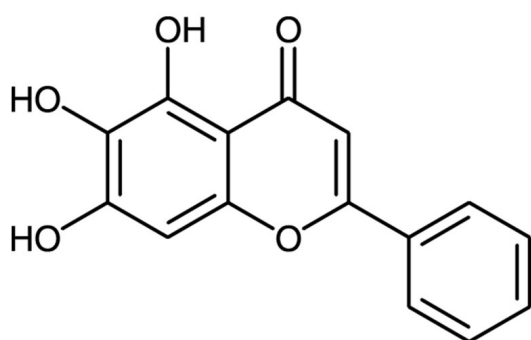
## Ordering Information

[Order Online »](#)

BML-EI106-0050	50mg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Subject to oxidation.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	5,6,7-Trihydroxyflavone
Appearance	Yellow to yellow-green crystalline powder.
CAS	491-67-8
Couple Target	Endopeptidase, Lipoxygenase
Couple Type	Inhibitor
Formula	$C_{15}H_{10}O_5$
MI	14: 942
MW	270.2
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO, acetone, 100% ethanol or methanol; almost insoluble in water.



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)