

Baicalein

Lipoxygenase inhibitor

Inhibits cellular 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 (IC_{50} =0.64 μM for 12-LO and 1.6 μM for 15-LO). Inhibits leukotriene biosynthesis (at higher concentrations) and release of lysosomal enzymes. Inhibits adjuvant-induced arthritis. Also inhibits prolyl oligopeptidase (POP/PREP)(IC_{50} =36 μM).

Citations: 21

[View Online »](#)

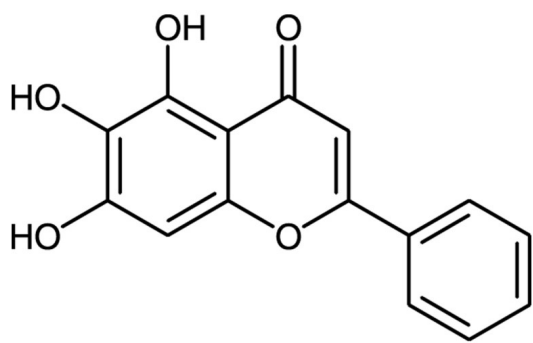
Ordering Information

[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.

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-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-uk@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com