Baicalein

Lipoxygenase inhibitor

Inhibits cellular Ca²⁺ 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 μ M for 12-LO and 1.6 μ 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)(IC $_{50}$ =36 μ M).

Citations: 21

View Online »

Ordering Information

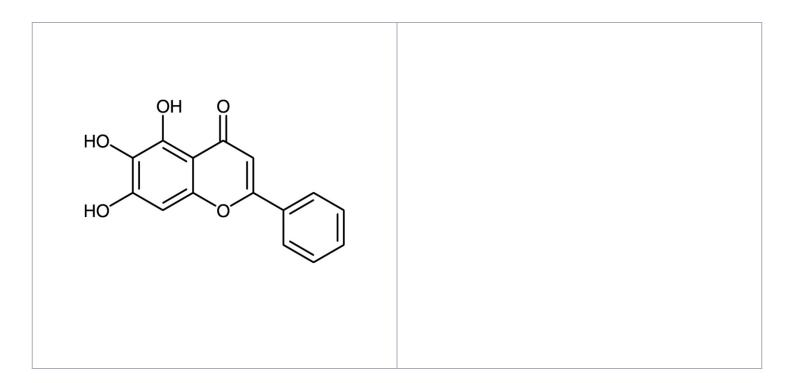
Order Online »

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)

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 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-

uk@enzolifesciences.com