

Apigenin

Antioxidant and chemopreventive

Apigenin is an antioxidant flavonoid. Has chemopreventive and antitumor properties. Induces apoptosis. Inhibits the proliferation of malignant tumor cells by G2/M arrest. Is a MAP-Kinase (MAPK/ERK) inhibitor. Inhibits hypoxia-inducible factor-1 (HIF-1) and vascular endothelial growth factor (VEGF) expression. Induces gastric relaxation in mouse stomach, mainly by suppressing Ca^{2+} influx through voltage-dependent calcium channels. Apigenin is a selective inhibitor of CK2. In WEHI 231 cells inhibition of CK2 by apigenin decreased the rate of I κ B α turnover and nuclear levels of NF- κ B (20 μ M). It blocks proliferation of Wnt-1 transfected cells and abrogates phosphorylation of β -catenin. In a rat model of glomerulonephritis, inhibition of CK2 by apigenin ameliorates renal dysfunction.

Citations: 16

[View Online »](#)

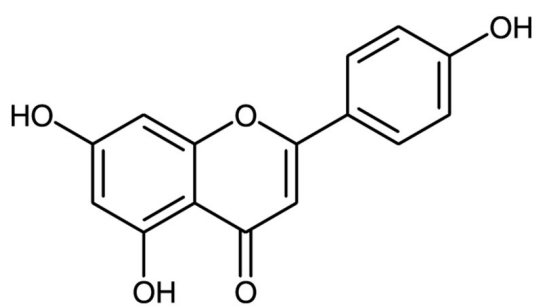
Ordering Information

[Order Online »](#)

BML-EI345-0100	100mg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance light yellow to slightly brown powder

CAS 520-36-5

Couple Target Casein kinase, HIF, MAPK

Couple Type Inhibitor

Formula $C_{15}H_{10}O_5$

MW 270.2

Purity ≥97% (HPLC)

Solubility Soluble in diluted potassium hydroxyde or DMSO, slightly soluble in 100% ethanol (hot). 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-
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)