

Quercetin . dihydrate

Antioxidant flavonoid

Antioxidant flavonoid. Inhibitor of mitochondrial ATPase, cAMP- and cGMP-phosphodiesterases. Inhibitor of protein tyrosine kinases and protein kinase C (PKC). Induces apoptosis. Blocks cells at the G0/G1 interface. Activator of human deacetylase SIRT1. Reversible inhibitor of fatty acid synthase (FAS). Inhibits the production of the inflammatory mediators nitric oxide (NO), TNF- α and IL-12 in activated macrophages. Induces gastric relaxation in mouse stomach, mainly by suppressing Ca^{2+} influx through voltage-dependent calcium channels.

Citations: 28

[View Online »](#)

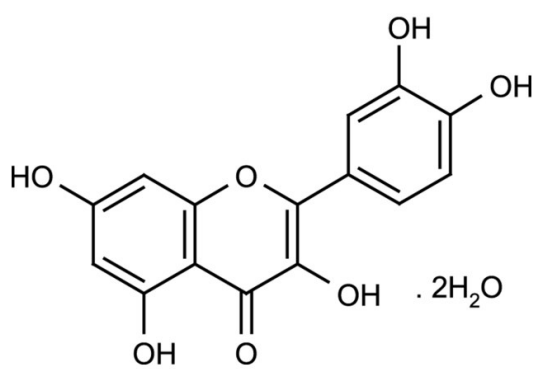
Ordering Information

[Order Online »](#)

ALX-385-001-G005	5g
------------------	----

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

Alternative Name 3,3',4',5,7-Pentahydroxyflavone . 2H₂O

Appearance Yellow powder with a green or brown cast.

CAS 6151-25-3

Couple Target ATPase, Fatty acid synthase, PDE, PKC, SIRT, SIRT1

Couple Type Activator, Inhibitor

Formula C₁₅H₁₀O₇ . 2H₂O

MI 14: 8034

MW 302.2 . 36.0

Purity ≥95% (HPLC)

RTECS LK8950000

Solubility Soluble in 100% ethanol or 1M sodium hydroxide; almost insoluble in water.

Source Isolated from *Sophora japonica* L.

Last modified: May 29, 2024



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)