

Wogonin

Anti-inflammatory and antioxidant

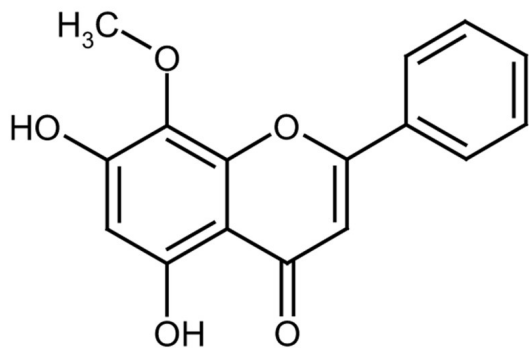
Cell permeable and orally available flavonoid. Induces apoptosis in tumor cells but not in healthy cells by increasing peroxide production. Displays anti-inflammatory properties. Suppresses the release of nitric oxide (NO) by inducible nitric oxide synthase (iNOS; NOS II), PGE₂ by cyclooxygenase-2 (COX-2), proinflammatory cytokines, MCP-1 gene expression and NF-κB activation. Antioxidant.

Citations: 15

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	5,7-Dihydroxy-8-methoxyflavone
Appearance	White to yellow solid.
CAS	632-85-9
Couple Target	COX, NOS
Couple Type	Inhibitor
Formula	$C_{16}H_{12}O_5$
Identity	Determined by 1H-NMR.
MW	284.3
Purity	≥98% (HPLC)
RTECS	LK8331000
Solubility	Soluble in DMSO (10mg/ml).

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-
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)