

Isorhamnetin

Antitumor agent

Antiviral agent. Antioxidant. Antitumor compound. Apoptosis inducer.

Citations: 8

[View Online »](#)

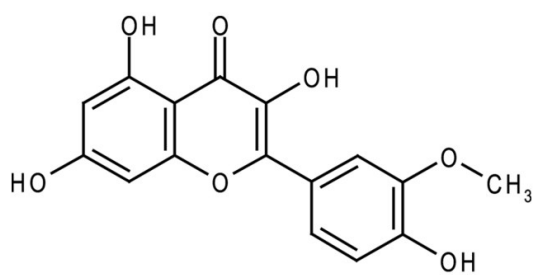
Ordering Information

[Order Online »](#)

ALX-385-024-M010	10mg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 3-Methylquercetin, 3'-Methoxy-3,4',5,7-tetrahydroxyflavone, 3,5,7-Trihydroxy-2-(4-hydroxy-3-methoxyphenyl)-4-benzopyrone

Appearance Yellow crystalline solid.

CAS 480-19-3

Formula $C_{16}H_{12}O_7$

Identity Determined by NMR and MS.

MW 316.3

Purity ≥98% (HPLC)

Solubility Soluble in acetone, methanol, ethylacetate or DMSO. Only sparingly soluble in water.

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)