

# Myricetin

## Chemopreventive agent

Antioxidant flavonoid. Has antitumor and chemopreventive properties. Anti-inflammatory. Inhibits NF- $\kappa$ B activation.  
PIM1 kinase inhibitor  $IC_{50}=0.78\mu\text{M}$ .

Citations: 10

[View Online »](#)

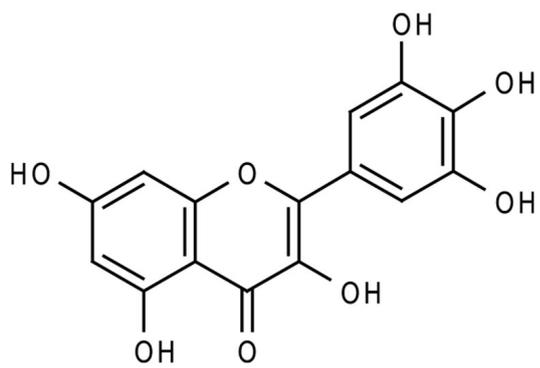
## Ordering Information

[Order Online »](#)

ALX-385-012-M010	10mg
------------------	------

## 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	-20°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	3,3',4',5,5',7-Hexahydroxyflavone
Appearance	Light yellow-green crystals.
CAS	529-44-2
Couple Target	NF-kappaB
Couple Type	Inhibitor
Formula	$C_{15}H_{10}O_8$
MI	14: 6332
MW	318.2
Purity	≥90% (HPLC)
RTECS	LK8646000
Solubility	Soluble in alcohol; slightly soluble in water (warm); insoluble in acetic acid.

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)