

Carnosic acid

PPAR γ activator. Antioxidant.

Naturally occurring phenolic compound with antioxidant properties. Inhibits lipid peroxidation induced by NADH or NADPH oxidation. Peroxisome proliferator-activated receptor γ (PPAR γ) activator. Anti-inflammatory. Antimicrobial.

Citations: 15

[View Online »](#)

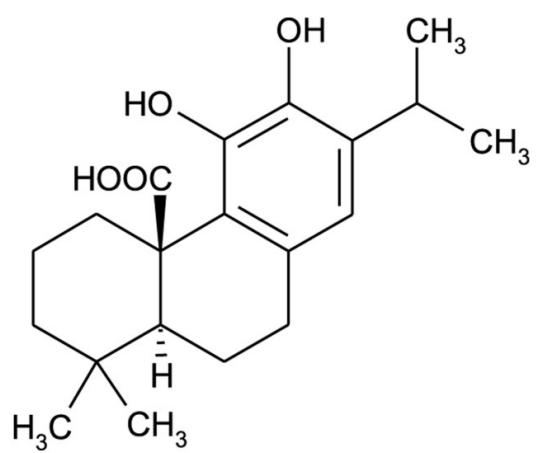
Ordering Information

[Order Online »](#)

ALX-270-264-M010	10mg
ALX-270-264-M050	50mg

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	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Yellow powder.
CAS	3650-09-7
Couple Target	PPAR
Couple Type	Activator
Formula	$C_{20}H_{28}O_4$
MW	332.4
Purity	≥95% (HPLC)
Source	Isolated from <i>Rosmarinus officinalis</i> .

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)