

(+)-Catechin . monohydrate

Free radical scavenger

Antioxidant flavonoid. Free radical scavenger. Has chemopreventive and antitumor properties.

Citations: 10

[View Online »](#)

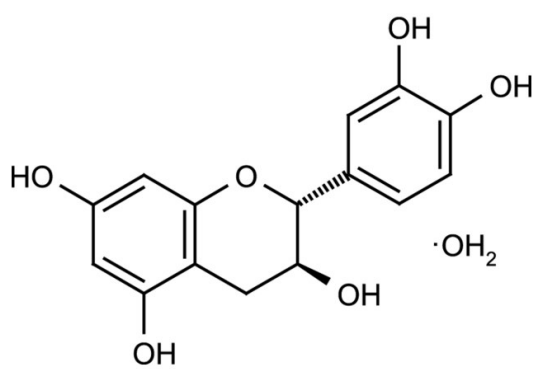
Ordering Information

[Order Online »](#)

ALX-385-017-G001	1g
------------------	----

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 (+)-3,3',4',5,7-Flavanpentol . H₂O

Appearance Off-white to yellow with a gray to tan cast.

CAS 225937-10-0

Formula $C_{15}H_{14}O_6 \cdot H_2O$

MI **14:** 1902

MW 290.3 (anhydrous basis)

Purity ≥98% (HPLC)

Solubility Soluble in 100% ethanol (50mg/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-
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)