

Captopril

Inhibitor of ACE and leukotriene hydrolase

First orally active angiotensin-converting enzyme (ACE) inhibitor. IC_{50} = 2.14 nM. Reversible and competitive inhibitor of LTA_4 hydrolase (IC_{50} = 11 μ M). Inhibits angiogenesis. Antihypotensive agent.

Citations: 3

[View Online »](#)

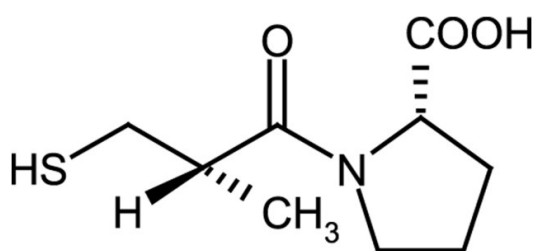
Ordering Information

[Order Online »](#)

ALX-270-212-G001	1g
ALX-270-212-G005	5g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (2S)-1-(3-Mercapto-2-methylpropionyl)-L-proline

Appearance White to off-white crystalline powder.

CAS 62571-86-2

Couple Target Angiotensin converting enzyme, Leukotriene hydrolase

Couple Type Inhibitor

Formula $C_9H_{15}NO_3S$

MI **14:** 1774

MW 217.3

Purity ≥98% (HPLC)

RTECS UY0550000

Solubility Soluble in water, chloroform, dichloromethane or 100% ethanol.

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)