

Alfacalcidol

Vitamin D₃ precursor

Converts to active calcitriol (1 α ,25-dihydroxyvitamin D₃; Prod. No. BML-DM200) *in vivo*. Inhibits bone resorption.

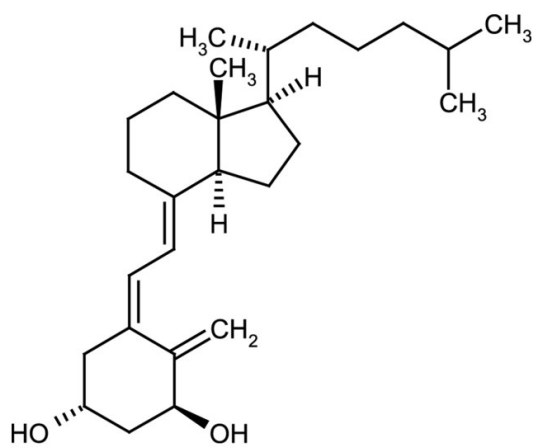
Ordering Information

[Order Online »](#)

ALX-460-036-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Keep under inert gas.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Hydroxyvitamin D3, 1-Hydroxycholecalciferol, (1 α ,3 β ,5Z,7E)-9,10-Secocholesta-5,7,10(19)-triene-1,3-diol
Appearance	White to off-white crystals.
CAS	41294-56-8
Formula	C ₂₇ H ₄₄ O ₂
MW	400.6
Purity	≥97% (Assay)
RTECS	VS2851000
Solubility	Soluble in 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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)