

25-Hydroxyvitamin D3

Precursor to 1 α ,25-dihydroxyvitamin D3

25-Hydroxyvitamin D3 is an Intermediate in the biosynthesis of 1 α ,25-dihydroxyvitamin D₃ and a major circulating metabolite of vitamin D₃. Plasma or serum levels of 25-hydroxyvitamin D3 can be determined by ELISA (Prod. No. ADI-900-215) and used as indicators of the vitamin D status of an individual.

Citations: 17

[View Online »](#)

Ordering Information

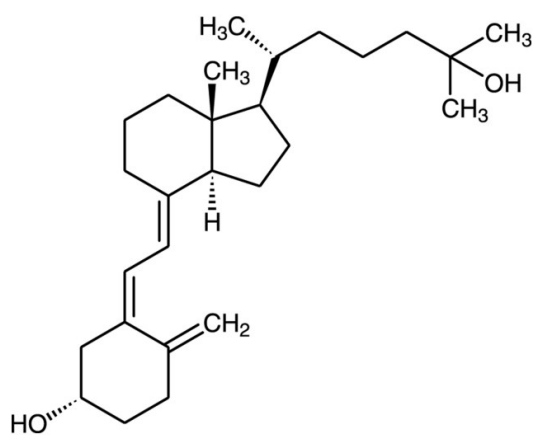
[Order Online »](#)

BML-DM100-0001	1mg
BML-DM100-0005	5mg

Manuals, SDS & CofA

[View Online »](#)

- Ultra-high purity
- Key indicator of vitamin D levels



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light and oxygen.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Calcifediol, 25-Hydroxycholecalciferol
Appearance	White to off-white crystalline solid.
CAS	19356-17-3
Couple Target	Vitamin D receptor
Couple Type	Ligand
Formula	$C_{27}H_{44}O_2$
Identity	Identity determined by MS and NMR.
MW	400.6
Purity	≥98% (HPLC)
Solubility	Soluble in 100% ethanol.
Source	Synthetic.

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)