

25-Hydroxyvitamin D₂

High purity vitamin D₂ intermediate

25-Hydroxyvitamin D₂ is an intermediate produced from vitamin D₂ by cytochrome P450 isoforms in serum. This primary metabolite of vitamin D₂ can be used to indirectly measure the concentration of vitamin D via HPLC and UV.

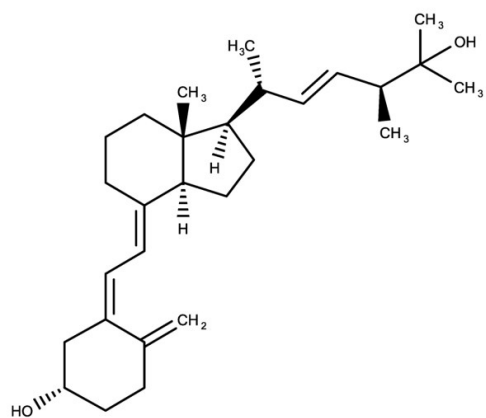
Ordering Information

[Order Online »](#)

BML-DM101-0001	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 and oxygen.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Ercalcidiol, 25-Hydroxyergocalciferol, 25-Hydroxycalciferol
Appearance	White crystalline powder.
CAS	21343-40-8
Couple Target	Vitamin D receptor
Couple Type	Ligand
Formula	$C_{28}H_{44}O_2$
MW	412.65
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO, ethanol or chloroform.

Last modified: March 19, 2025



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)