

24(R),25-Dihydroxyvitamin D3

Major metabolite of vitamin D3

Major metabolite of vitamin D₃. Mediates calcium and phosphorous homeostasis. Inhibits 1 α ,25-dihydroxyvitamin D3 and testosterone potentiation of calcium channels in osteosarcoma cells.

Citations: 4

[View Online »](#)

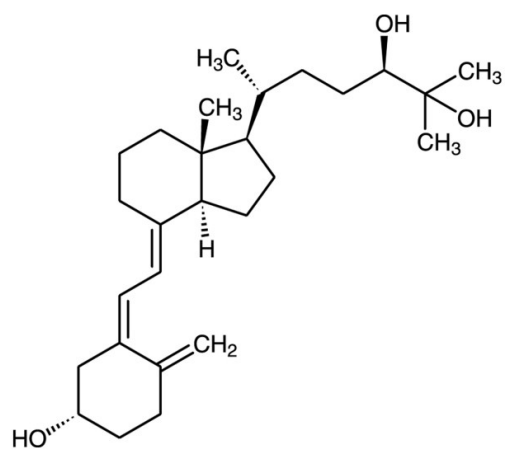
Ordering Information

[Order Online »](#)

BML-DM300-0050	50 μ g
----------------	------------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -80°C for up to 1 year.
Long Term Storage	-80°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

CAS	55721-11-4
Formula	$C_{27}H_{44}O_3$
Formulation	White solid dissolved in ethanol (50µg/50µl).
MW	416.6
Purity	≥98% (TLC)

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-
be@enzolifesciences.com](mailto:info-be@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)