

Calcitriol

Active form of vitamin D, regulates dietary Ca²⁺ absorption; VDR agonist.

The active hormonal form of vitamin D3. Induces cell differentiation and prevents proliferation of cancer cells. Along with dietary calcium, reduces the risk of colon cancer.

Ordering Information

[Order Online »](#)

LKT-C0145-M001	1mg
LKT-C0145-C050	50µg
LKT-C0145-5050	5x50µg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Air, light and moisture sensitive.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1- α , 25-Dihydroxyvitamin D3

CAS 32222-06-3

Formula $C_{27}H_{44}O_3$

MI 13: 1643

MW 416.64

Purity $\geq 97\%$

Solubility Slightly soluble in ethanol, ethyl acetate. Soluble in DMSO.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com