

# Calcitriol

Active form of vitamin D, regulates dietary Ca<sup>2+</sup> absorption; VDR agonist.

The active hormonal form of vitamin D<sub>3</sub>. Induces cell differentiation and prevents proliferation of cancer cells. Along with dietary calcium, reduces the risk of colon cancer.

## Ordering Information

[Order Online »](#)

LKT-C0145-5050	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- $\alpha$ , 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](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)