

# 27-Hydroxycholesterol

**Cholesterol metabolite linked to breast cancer, neurodegenerative disease, and metabolism**

27-Hydroxycholesterol (27-HC) is a metabolite resulting from the action of sterol 27-Hydroxylase on cholesterol. It is an endogenous selective estrogen receptor modulator and ligand for the liver X receptor (LXR). 27-Hydroxycholesterol plays a role in Alzheimer's disease, oxidative stress, metabolism, cell-signaling, and heart disease.

**Citations:** 10

[View Online »](#)

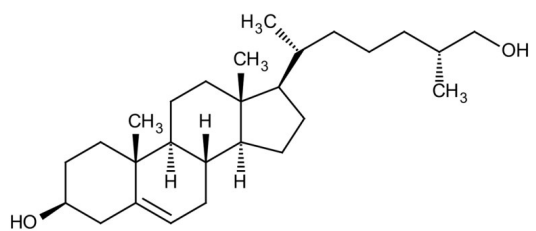
## Ordering Information

[Order Online »](#)

ENZ-CHM102-0001	1mg
ENZ-CHM102-0005	5mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable, as supplied, for 2 years. Solutions in DMSO or ethanol may be stored at -20°C for up to 2 months.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	5,25R- $\alpha$ -Cholest $\alpha$ -en- $\alpha$ 3 $\beta$ , $\alpha$ 26- $\alpha$ -diol, (25R)-26-Hydroxycholesterol
Appearance	White solid.
CAS	20380-11-4
Formula	C <sub>27</sub> H <sub>46</sub> O <sub>2</sub>
MW	402.7
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (up to 10 mg/ml with warming) or in ethanol (up to 10 mg/ml). Insoluble in water.

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)