

24(S)- Hydroxycholesterol

LXR ligand

Endogenous agonist for the nuclear receptor LXR. Important metabolite of cholesterol in the brain. Displays neurotoxic effects on human neuroblastoma cells. Cerebrospinal fluid concentrations of 24(S)-hydroxycholesterol are increased in patients with Alzheimer's disease and mild cognitive impairment.

Citations: 9

[View Online »](#)

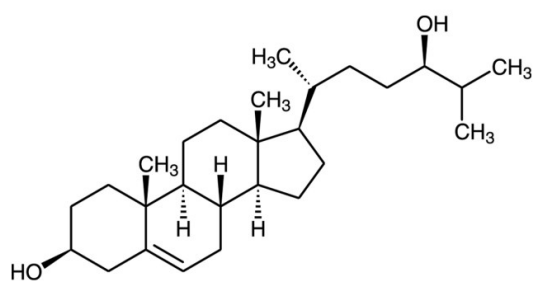
Ordering Information

[Order Online »](#)

BML-GR230-0001	1mg
BML-GR230-0005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store solutions at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	474-73-7
Couple Target	LXR
Couple Type	Ligand
Formula	$C_{27}H_{46}O_2$
MW	402.7
MeltingPoint	177-180°C
Purity	≥98% (TLC)
Solubility	Soluble in ethanol (10mg/ml) or DMSO (10mg/ml – requires warming to ~40°C).

Last modified: May 29, 2024



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-fr@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