

24(S),25-Epoxycholesterol

LXR ligand

Endogenous agonist for the nuclear receptor LXR.

Citations: 3

[View Online »](#)

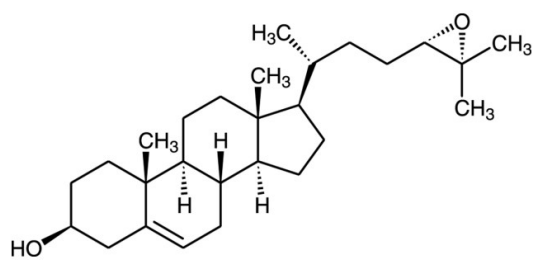
Ordering Information

[Order Online »](#)

BML-GR231-0001	1mg
BML-GR231-0005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Product should be used immediately following dissolution.
Handling	Store with desiccant. Product packaged under inert gas.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White solid.
CAS	77058-74-3
Couple Target	LXR
Couple Type	Ligand
Formula	$C_{27}H_{44}O_2$
Identity	Determined by NMR.
MW	400.6
MeltingPoint	156-159°C
Purity	≥98% (TLC)
Solubility	Soluble in ethanol (10mg/ml) or DMSO (10mg/ml – requires warming to ~40°C which has been shown to lead to significant decomposition of the product).

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@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)