

9-cis-Retinoic acid

Retinoid X receptor ligand

A high affinity ligand for the retinoid X receptor (RXR). Destabilizes VDR-RXR heterodimer-DNA complex.

Citations: 7

[View Online »](#)

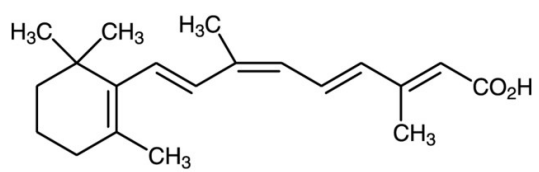
Ordering Information

[Order Online »](#)

BML-GR101-0005	5mg
BML-GR101-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Store tightly sealed in dark. Protect from light.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Alitretinoin

Appearance Light yellow to orange crystalline powder.

CAS 5300-03-8

Couple Target Retinoid X receptor

Couple Type Ligand

Formula $C_{20}H_{28}O_2$

MW 300.4

Purity ≥98% (UHPLC)

Solubility Soluble in DMSO (25mg/ml) or 100% ethanol (25mg/ml).

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-
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)