

Retinoic acid, all trans

Retinoic acid receptor ligand

Retinoic acid is the active metabolite of vitamin A. It is a potent modulator of growth and differentiation exerting its effects by binding to nuclear retinoic acid receptors (RARs) which directly regulate gene expression. All trans retinoic acid binds to RAR α , β and γ subtypes with K_d values of 0.2, 0.4 and 0.2 nM respectively. RARs function as ligand-dependent transcription factors. Retinoic Acid also promotes the differentiation of mouse embryonic stem cells into adipocytes, neurons and glia in vitro. It guides development of the posterior part of the embryo.

Citations: 13

[View Online »](#)

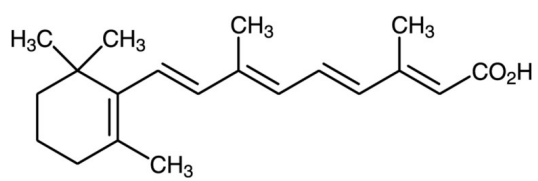
Ordering Information

[Order Online »](#)

BML-GR100-0500	500mg
BML-GR100-5000	5g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Yellow solid.

CAS 302-79-4

Couple Target Retinoic acid receptor

Couple Type Ligand

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

MW 300.4

Purity ≥95%

Solubility Soluble in DMSO (25 mg/ml) or ethanol (10 mg/ml).



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)