

4-

Hydroxyphenylretinamide

Retinoic acid receptor ligand

An analog of all *trans* retinoic acid that shows antiproliferative activity on cultured human breast cancer cells. Highly selective activator of retinoid receptors. Induces apoptosis in malignant hemopoietic cell lines.

Citations: 6

[View Online »](#)

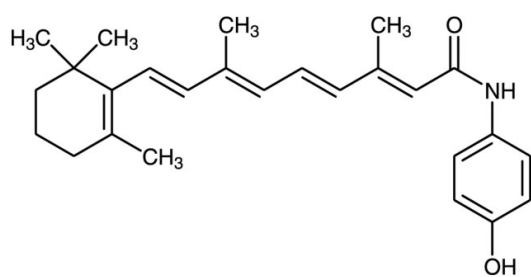
Ordering Information

[Order Online »](#)

BML-GR103-0005	5mg
----------------	-----

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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Fenretinide

Appearance Orange solid.

CAS 65646-68-6

Couple Target Retinoic acid receptor

Couple Type Ligand

Formula $C_{26}H_{33}NO_2$

MW 391.5

Purity ≥97% (TLC)

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

Source Synthetic.

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-
be@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)