

Methoprene acid

Retinoid X receptor ligand

Methoprene acid is an analog of the insect growth regulator, juvenile hormone. It binds directly to RXR and is a transcriptional activator in both insect and mammalian cells. It is RXR specific and does not activate the RAR pathway.

Citations: 1

[View Online »](#)

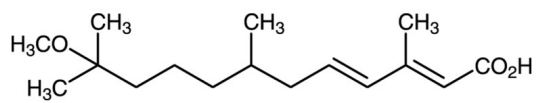
Ordering Information

[Order Online »](#)

BML-GR106-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. 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	Pale yellow oil.
CAS	53092-52-7
Couple Target	Retinoid X receptor
Couple Type	Ligand
Formula	$C_{16}H_{28}O_3$
MW	268.4
Purity	≥98% (TLC)
Solubility	Soluble in DMSO or methanol.

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-be@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