

GW-9662

PPAR ligand

Potent, selective and irreversible PPAR γ antagonist. Blocks the PPAR γ -induced differentiation of monocytes to osteoclasts by >90% at a dose of 0.1 μ M. Much more potent antagonist than BADGE (Prod. No. ALX-270-353). Can be used to antagonize the effects of PPAR γ agonists in a variety of cell types.

Citations: 14

[View Online »](#)

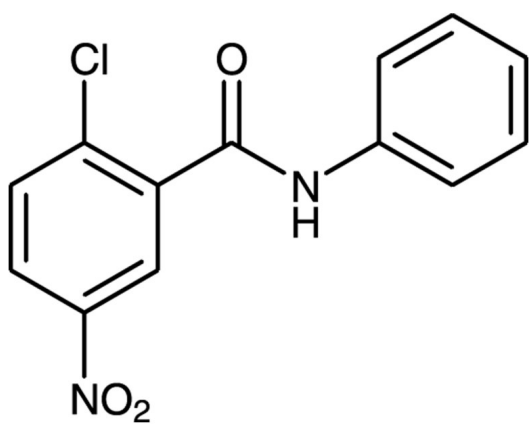
Ordering Information

[Order Online »](#)

BML-GR234-0010	10mg
BML-GR234-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at +20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-Chloro-5-nitrobenzanilide
Appearance	White to off-white solid.
CAS	22978-25-2
Couple Target	PPAR
Couple Type	Ligand
Formula	$C_{13}H_9N_2O_3Cl$
MW	276.7
MeltingPoint	155-157°C
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (50mg/ml), 100% ethanol (10mg/ml) or dimethyl formamide. Insoluble in water.
Technical Info / Product Notes	Note: Product is not sterile.

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)