

T0070907

PPAR ligand

Potent and selective PPAR γ antagonist. T0070907 blocks PPAR γ function in both cell-based reporter gene and adipocyte differentiation assays. It inhibits growth and survival of hepatocellular carcinoma cells and squamous cell carcinoma cell lines by interfering with cell adhesion "anoikis".

Citations: 7

[View Online »](#)

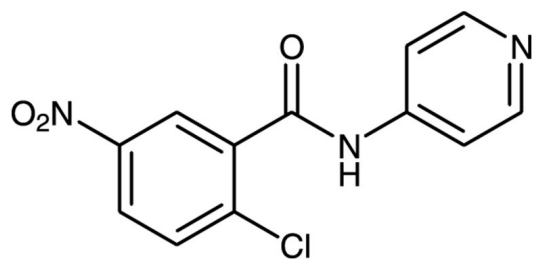
Ordering Information

[Order Online »](#)

BML-GR212-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at room temperature for up to 1 year. Store solutions at -20°C for up to 3 months.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	2-Chloro-5-nitro-N-4-pyridinylbenzamide
Appearance	Off-white solid.
CAS	313516-66-4
Couple Target	PPAR
Couple Type	Ligand
Formula	$C_{12}H_8ClN_3O_3$
MW	277.7
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (>25 mg/ml).

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-fr@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