

GW1929

PPAR γ agonist

Potent (1'000-fold) PPAR γ agonist which does not contain a thiazolidinedione moiety. The glucose-lowering effect of GW1929 in rats is 100-fold more potent than that of troglitazone (Prod. No. BML-GR210).

Citations: 5

[View Online »](#)

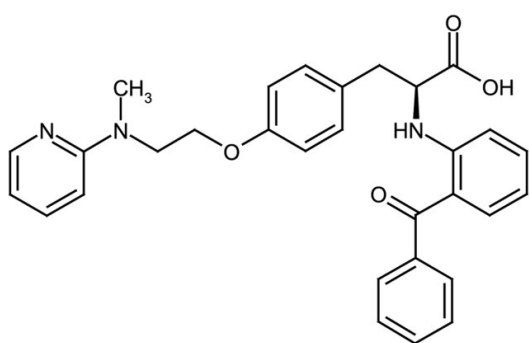
Ordering Information

[Order Online »](#)

ALX-420-029-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-(2-Benzoylphenyl)-L-tyrosine PPAR γ Agonist, (2S)-((2-Benzoylphenyl)amino-3-[4-[2-(methylpyridin-2-ylamino)ethoxy]phenyl]propionic acid
Appearance	Off-white to yellow crystalline solid.
CAS	196808-24-9
Couple Target	PPAR
Couple Type	Activator, Ligand
Formula	C ₃₀ H ₂₉ N ₃ O ₄
MW	495.6
Purity	≥97% (HPLC)
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-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