

GW9508

High potency synthetic agonist at GPR40

First high potency synthetic agonist at GPR40.

GPR40 was formerly an orphan GPCR whose endogenous ligands have been shown to be free fatty acids and has therefore been designated the free fatty acid receptor 1 (FFAR1). It is abundantly expressed on insulin-expressing beta cells and has been shown to mediate the majority of the effects of free fatty acids on insulin secretion. GW9508 enhanced insulin secretion in INS-1E cells and activates ATP-sensitive potassium channels in rat pancreatic beta cells (40µM). It is a useful tool for exploring the biology of GPR40 mediated signaling and physiology.

Citations: 5

[View Online »](#)

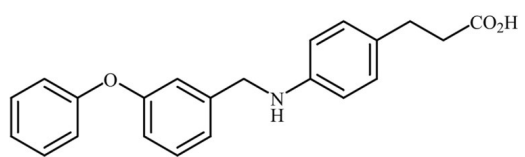
Ordering Information

[Order Online »](#)

BML-GP102-0010	10mg
BML-GP102-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for 1 year as supplied. Solutions can be stored 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	4-[[[(3-Phenoxyphenyl)methyl]amino]benzene propanoic acid
Appearance	Off-white solid.
CAS	885101-89-3
Formula	$C_{22}H_{21}NO_3$
MW	347.4
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (15mg/ml).

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-
be@enzolifesciences.com](mailto:info-be@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)