

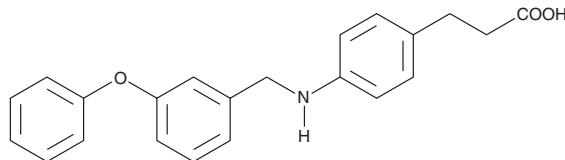
Product Information



GW 9508

Catalog No. 10008907

CAS Registry No.: 885101-89-3
Formal Name: 4-[[[(3-phenoxyphenyl)methyl]amino]-benzenepropanoic acid
MF: C₂₂H₂₁NO₃
FW: 347.4
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that GW 9508 be stored as supplied at -20°C. It will be stable for at least one year.

GW 9508 is supplied as a crystalline solid. A stock solution may be made by dissolving the GW 9508 in an organic solvent purged with an inert gas. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of GW 9508 in these solvents is approximately 15 mg/ml.

GW 9508 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GW 9508 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. GW 9508 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

GPR40 and GPR120 are G protein-coupled receptors (GPCRs) that are activated by medium and long-chain fatty acids. In addition, there is evidence of a link between GPR40 and the ability of fatty acids to acutely potentiate insulin secretion. GW 9508 is a small-molecule agonist of GPR40 and GPR120. Stimulation of intracellular Ca²⁺ mobilization by GW9508 in human embryonic kidney (HEK) 293 cells expressing GPR40 and GPR120 was observed with EC₅₀ values of 47 nM and 2.2 μM, respectively.¹ GW 9508 dose dependently potentiates glucose-stimulated insulin secretion and the potassium chloride-mediated increase in insulin secretion in MIN6 cells.¹

Reference

1. Briscoe, C.P., Peat, A.J., McKeown, S.C., *et al.* Pharmacological regulation of insulin secretion in MIN6 cells through the fatty acid receptor GPR40: Identification of agonist and antagonist small molecules. *Br. J. Pharmacol.* **148**, 619-628 (2006)

Related Product

GPR40 Polyclonal Antibody - Cat. No. 10007205

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 09/09/2009