

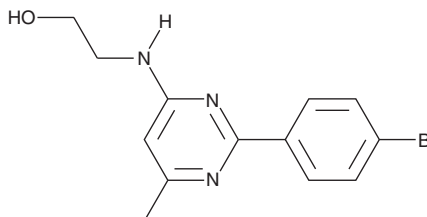
PRODUCT INFORMATION



AS-1269574

Item No. 10626

CAS Registry No.: 330981-72-1
Formal Name: 2-[[2-(4-bromophenyl)-6-methyl-4-pyrimidinyl]amino]-ethanol
MF: C₁₃H₁₄BrN₃O
FW: 308.2
Purity: ≥98%
UV/Vis.: λ_{max}: 253 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

AS-1269574 is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of AS-1269574 in these solvents is approximately 20 and 30 mg/ml, respectively.

AS-1269574 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of AS-1269574 should be diluted with the aqueous buffer of choice. AS-1269574 has a solubility of 0.1 mg/ml in a 1:10 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

GPR119, a glucose-dependent insulinotropic receptor, is a Gs protein-coupled receptor that is expressed in pancreas β-cells and intestinal L-cells.¹ Activation of GPR119 by the endogenous ligands lyso-phosphatidylcholine and oleoyl ethanolamide increases intracellular cAMP levels and promotes glucose-stimulated insulin secretion.^{1,2} AS-1269574 is an agonist of GPR119 that is effective both in isolated cells and *in vivo*.³ It increases cAMP levels in HEK293 cells transfected with human GPR119 (EC₅₀ = 2.5 μM) and glucose-stimulated insulin secretion in mice (100 mg/kg).³

References

- Overton, H.A., Babbs, A.J., Doel, S.M., *et al.* Deorphanization of a G protein-coupled receptor for oleoylethanolamide and its use in the discovery of small-molecule hypophagic agents. *Cell Metab.* **3**(3), 167-175 (2006).
- Soga, T., Ohishi, T., Matsui, T., *et al.* Lysophosphatidylcholine enhances glucose-dependent insulin secretion via an orphan G-protein-coupled receptor. *Biochem. Biophys. Res. Commun.* **326**(4), 744-751 (2005).
- Yoshida, S., Ohishi, T., Matsui, T., *et al.* Identification of a novel GPR119 agonist, AS1269574, with *in vitro* and *in vivo* glucose-stimulated insulin secretion. *Biochem. Biophys. Res. Commun.* **400**(3), 437-441 (2010).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/27/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM