

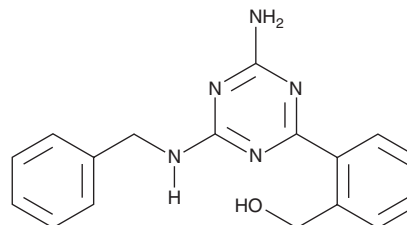
PRODUCT INFORMATION



Ogerin

Item No. 34228

CAS Registry No.: 1309198-71-7
Formal Name: 2-[4-amino-6-[(phenylmethyl)amino]-1,3,5-triazin-2-yl]-benzenemethanol
Synonym: ZINC67740571
MF: C₁₇H₁₇N₅O
FW: 307.3
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Ogerin is supplied as a solid. A stock solution may be made by dissolving the ogerin in the solvent of choice, which should be purged with an inert gas. Ogerin is soluble in the organic solvent DMSO at a concentration of approximately 50 mg/ml.

Description

Ogerin is a positive allosteric modulator of G protein-coupled receptor 68 (GPR68).¹ It increases proton-induced cAMP production in HEK293T cells expressing the human receptor in a reporter assay. Ogerin inhibits proton-mediated calcium release in the same cells when used at a concentration of 10 μM. It impairs contextual, but not cued, fear conditioning in wild-type, but not *Gpr68* knockout, mice when administered at a dose of 10 mg/kg.

Reference

1. Huang, X.-P., Karpiak, J., Kroeze, W.K., *et al.* Allosteric ligands for the pharmacologically dark receptors GPR68 and GPR65. *Nature* **527(7579)**, 477-483 (2015).

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, 12/08/2022

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