

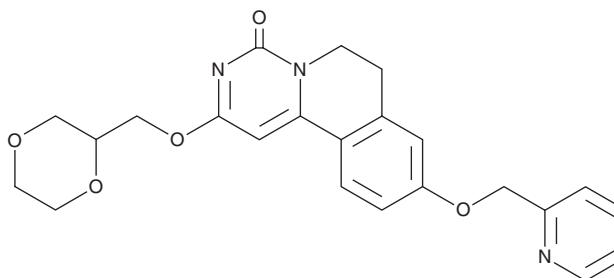
PRODUCT INFORMATION



GPR84 Antagonist 8

Item No. 43485

CAS Registry No.: 1445846-30-9
Formal Name: 2-(1,4-dioxan-2-ylmethoxy)-6,7-dihydro-9-(2-pyridinylmethoxy)-4H-pyrimido[6,1-a]isoquinolin-4-one
Synonyms: EN 818470, G Protein-coupled Receptor 84 Antagonist 8
MF: C₂₃H₂₃N₃O₅
FW: 421.5
Purity: ≥98%
Supplied as: A crystalline 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

GPR84 antagonist 8 is supplied as a crystalline solid. A stock solution may be made by dissolving the GPR84 antagonist 8 in the solvent of choice, which should be purged with an inert gas. GPR84 antagonist 8 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

GPR84 antagonist 8 is an antagonist of G protein-coupled receptor 84 (GPR84).¹ It inhibits calcium mobilization induced by 6-OAU (Item No. 17687) in HEK293 cells expressing human GPR84 and Gα₁₆ (IC₅₀ = 716 nM). GPR84 antagonist 8 (10 μM) inhibits 6-OAU-enhanced phagocytosis of *E. coli* by primary mouse bone marrow-derived macrophages (BMDMs).²

References

1. Chen, L.-H., Zhang, Q., Xiao, Y.-F., *et al.* Phosphodiesterases as GPR84 antagonists for the treatment of ulcerative colitis. *J. Med. Chem.* **65**(5), 3991-4006 (2022).
2. Recio, C., Lucy, D., Purvis, G.S.D., *et al.* Activation of the immune-metabolic receptor GPR84 enhances inflammation and phagocytosis in macrophages. *Front. Immunol.* **9**, 1419 (2018).

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, 07/14/2025

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