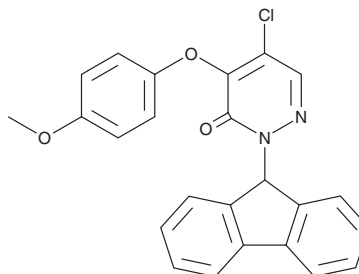


PRODUCT INFORMATION



CYM 50769
Item No. 34484

CAS Registry No.: 1421365-63-0
Formal Name: 5-chloro-2-(9H-fluoren-9-yl)-4-(4-methoxyphenoxy)-3(2H)-pyridazinone
MF: C₂₄H₁₇ClN₂O₃
FW: 416.9
Purity: ≥98%
UV/Vis.: λ_{max}: 269 nm
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

CYM 50769 is supplied as a solid. A stock solution may be made by dissolving the CYM 50769 in the solvent of choice, which should be purged with an inert gas. CYM 50769 is soluble in the organic solvent DMSO.

Description

CYM 50769 is a nonpeptide antagonist of neuropeptides B and W receptor 1 (NPBWR1).¹ It inhibits neuropeptide W-induced activation of NPBWR1 (IC₅₀ = 0.12 μM). CYM 50769 is selective for NPBWR1 over a panel of 30 G protein-coupled receptors (GPCRs), enzymes, transporters, and ion channels at 30 μM.

Reference

- Guerrero, M., Urbano, M., Schaeffer, M.-T., *et al.* SAR analysis of novel non-peptidic NPBWR1 (GPR7) antagonists. *Bioorg. Med. Chem. Lett.* **23**(3), 614-619 (2013).

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, 11/04/2022

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