

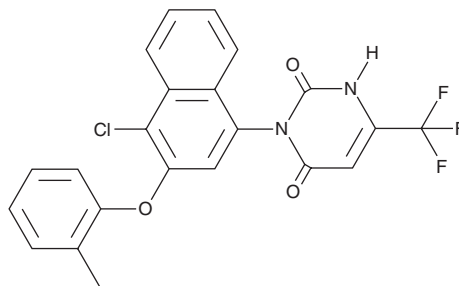
PRODUCT INFORMATION



BAY-069

Item No. 41149

CAS Registry No.: 2639638-66-5
Formal Name: 3R-[4-chloro-3-(2-methylphenoxy)-1-naphthalenyl]-6-(trifluoromethyl)-2,4(1H,3H)-pyrimidinedione
MF: C₂₂H₁₄ClF₃N₂O₃
FW: 446.8
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

BAY-069 is supplied as a solid. A stock solution may be made by dissolving the BAY-069 in the solvent of choice, which should be purged with an inert gas. BAY-069 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

BAY-069 is an inhibitor of branched-chain amino acid transaminase 1 (BCAT1) and BCAT2 (IC₅₀s = 31 and 153 nM, respectively).¹ It is selective for BCAT1 and BCAT2 over a panel of 77 enzymes, receptors, and transporters at 10 μM. BAY-069 increases the levels of branched-chain amino acids (BCAAs) in U87MG brain and MDA-MB-231 breast cancer cells (IC₅₀s = 358 and 874 nM, respectively) but not in the presence of human serum albumin (HSA) or BSA (IC₅₀s = >50 μM for both).

Reference

1. Günther, J., Hillig, R.C., Zimmermann, K., *et al.* BAY-069, a novel (trifluoromethyl)pyrimidinedione-based BCAT1/2 inhibitor and chemical probe. *J. Med. Chem.* **65**(21), 14366-14390 (2024).

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, 06/26/2024

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