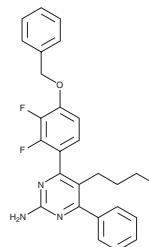


PRODUCT INFORMATION

5-Butyl-4-[2,3-difluoro-4-(phenylmethoxy)phenyl]-6-phenyl-2-pyrimidinamine

Item No. 31338

CAS Registry No.: 2253745-78-5
MF: C₂₇H₂₅F₂N₃O
FW: 445.5
Purity: ≥98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

5-Butyl-4-[2,3-difluoro-4-(phenylmethoxy)phenyl]-6-phenyl-2-pyrimidinamine is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol and chloroform purged with an inert gas can be used. The solubility of 5-butyl-4-[2,3-difluoro-4-(phenylmethoxy)phenyl]-6-phenyl-2-pyrimidinamine in these solvents is approximately 10 mg/ml.

Description

5-Butyl-4-[2,3-difluoro-4-(phenylmethoxy)phenyl]-6-phenyl-2-pyrimidinamine is an inhibitor of microsomal prostaglandin E synthase-1 (mPGES-1; IC₅₀s = 66 and 2,730 nM for the human and mouse enzymes, respectively).¹ It is selective for mPGES-1 over COX-1 but does inhibit COX-2 at 20 μM. It also inhibits LPS-induced prostaglandin E₂ (PGE₂; Item No. 14010) production in mouse peritoneal cells (IC₅₀ = 12 nM). *In vivo*, 5-butyl-4-[2,3-difluoro-4-(phenylmethoxy)phenyl]-6-phenyl-2-pyrimidinamine (25 mg/kg) reduces carrageenan-induced paw edema in rats.

Reference

1. Kalčić, F., Kolman, V., Ajani, H., *et al.* Polysubstituted pyrimidines as mPGES-1 inhibitors: Discovery of potent inhibitors of PGE₂ production with strong anti-inflammatory effects in carrageenan-induced rat paw edema. *ChemMedChem* **15**(15), 1398-1407 (2020).

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, 08/18/2020

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