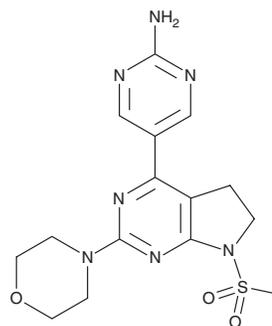


PRODUCT INFORMATION



CH5132799
Item No. 26410

CAS Registry No.: 1007207-67-1
Formal Name: 5-[6,7-dihydro-7-(methylsulfonyl)-2-(4-morpholinyl)-5H-pyrrolo[2,3-d]pyrimidin-4-yl]-2-pyrimidinamine
MF: C₁₅H₁₉N₇O₃S
FW: 377.4
Purity: ≥95%
UV/Vis.: λ_{max}: 276, 337 nm
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

CH5132799 is supplied as a crystalline solid. A stock solution may be made by dissolving the CH5132799 in the solvent of choice, which should be purged with an inert gas. CH5132799 is soluble in the organic solvent DMSO at a concentration of approximately 12 mg/ml.

Description

CH5132799 is an inhibitor of class I PI3K isoforms (IC₅₀s = 14, 120, 500, and 36 nM for PI3Kα, PI3Kβ, PI3Kδ, and PI3Kγ, respectively).¹ It is selective for these PI3K isoforms over PI3KC2α and PI3KC2β, which are class II PI3Ks, Vps34, a class III PI3K, and mTOR (IC₅₀s = 5,300, >10,000, >10,000, and 1,600 nM, respectively), as well as 26 additional protein kinases (IC₅₀s = >10,000 nM). CH5132799 inhibits proliferation of HCT116, KPL-4, T47D, and SKOV3 cancer cells expressing the activating mutation PI3Kα^{H1047R} (IC₅₀s = 200, 32, 56, and 120 nM, respectively). It also reduces tumor growth in a PC3 mouse xenograft model when administered at a dose of 25 mg/kg per day.

Reference

- Ohwada, J., Ebiike, H., Kawada, H., *et al.* Discovery and biological activity of a novel class I PI3K inhibitor, CH5132799. *Bioorg. Med. Chem. Lett.* **21(6)**, 1767-1772 (2011).

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, 01/03/2023

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