

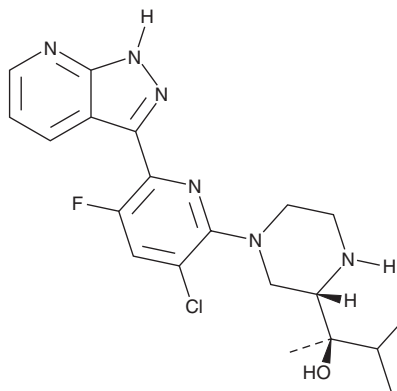
PRODUCT INFORMATION



VTX-27

Item No. 36938

CAS Registry No.: 1321924-70-2
Formal Name: (αR,2S)-4-[3-chloro-5-fluoro-6-(1H-pyrazolo[3,4-b]pyridin-3-yl)-2-pyridinyl]-α-methyl-α-(1-methylethyl)-2-piperazinemethanol
MF: C₂₀H₂₄ClFN₆O
FW: 418.9
Purity: ≥95%
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

VTX-27 is supplied as a solid. A stock solution may be made by dissolving the VTX-27 in the solvent of choice, which should be purged with an inert gas. VTX-27 is soluble in DMSO.

Description

VTX-27 is a PKCθ inhibitor ($K_i = 0.08$ nM).¹ It is selective for PKCθ over PKCα, -βI, -βII, -γ, -δ, -ε, -η, -μ, and -ζ with K_i values ranging from 1 to >5,000 nM as well as a panel of seven kinases from the Src, spleen tyrosine kinase (Syk), Tec, and MAP kinase families (K_i s = >1,000 nM for all). VTX-27 inhibits CD28-induced IL-2 release in isolated human peripheral blood mononuclear cells (PBMCs; $IC_{50} = 11$ nM). Oral administration of VTX-27 (12.5, 25, and 50 mg/kg) prevents staphylococcal enterotoxin B-induced increases in IL-2 plasma levels in mice.

Reference

1. Jimenez, J.-M., Boyall, D., Brenchley, G., *et al.* Design and optimization of selective protein kinase C θ (PKCθ) inhibitors for the treatment of autoimmune diseases. *J. Med. Chem.* **56**(5), 1799-1810 (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, 12/02/2022

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