

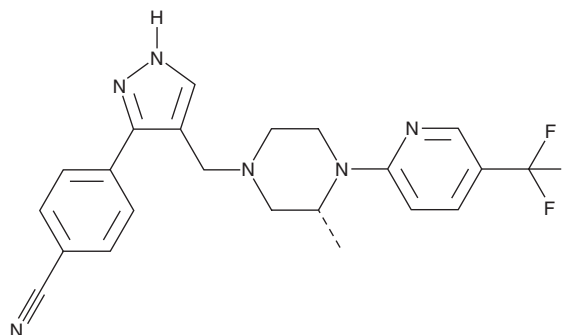
PRODUCT INFORMATION



GNF-362

Item No. 35866

CAS Registry No.: 1003019-41-7
Formal Name: 4-[4-[[[(3R)-3-methyl-4-[5-(trifluoromethyl)-2-pyridinyl]-1-piperazinyl]methyl]-1H-pyrazol-3-yl]-benzonitrile
MF: C₂₂H₂₁F₃N₆
FW: 426.4
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

GNF-362 is supplied as a solid. A stock solution may be made by dissolving the GNF-362 in the solvent of choice, which should be purged with an inert gas. GNF-362 is soluble in DMSO, ethanol, and methanol.

Description

GNF-362 is an inositol triphosphate 3 kinase (ITPK) inhibitor (IC₅₀s = 20, 9, and 19 nM for ITPKA, ITPKB, and ITPKC, respectively).¹ It induces calcium influx in isolated mouse splenocytes (EC₅₀ = 12 nM). *In vivo*, GNF-362 (10 and 25 mg/kg) reduces the number of thymic CD4⁺ T cells in mice. It also reduces inflammatory cell infiltration, joint erosion, and proteoglycan loss, in a rat model of antigen-induced arthritis when administered at a dose of 20 mg/kg.

Reference

1. Miller, A.T., Dahlberg, C., Sandberg, M.L., *et al.* Inhibition of the inositol kinase Itpkb augments calcium signaling in lymphocytes and reveals a novel strategy to treat autoimmune disease. *PLoS One* **10(6)**, e0131071 (2015).

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, 02/14/2024

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