

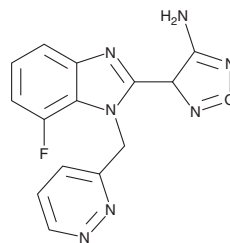
PRODUCT INFORMATION



CVN293

Item No. 40983

CAS Registry No.: 2815296-08-1
Formal Name: 4-[7-fluoro-1-(3-pyridazinylmethyl)-1H-benzimidazol-2-yl]-1,2,5-oxadiazol-3-amine
MF: C₁₄H₁₀FN₇O
FW: 311.3
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

CVN293 is supplied as a solid. A stock solution may be made by dissolving the CVN293 in the solvent of choice, which should be purged with an inert gas. CVN293 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

CVN293 is an inhibitor of the two-pore domain potassium channel K_{2p}13.1/THIK1 (IC₅₀ = 49 nM).¹ It is selective for K_{2p}13.1/THIK1 over K_{2p}6.1/TWIK2 and K_{2p}2.1/TREK-1 channels (IC₅₀s = >30,000 nM for both). CVN293 inhibits NOD-like receptor protein 3 (NLRP3) inflammasome-mediated IL-1β production in LPS-stimulated primary neonatal murine microglia (IC₅₀ = 24 nM).

Reference

1. Bürli, R.W., Doyle, K.J., Dickson, L., *et al.* Discovery of CVN293, a brain permeable KCNK13 (THIK-1) inhibitor suitable for clinical assessment. *ACS Med. Chem. Lett.* **15**(5), 646-652 (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, 12/10/2024

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