

PRODUCT INFORMATION



PFI-653

Item No. 40955

CAS Registry No.: 2173134-00-2

Formal Name: 8-oxa-2-azaspiro[4.5]dec-2-yl[2-[(2-pyrazinylmethyl)amino]-5-pyrimidinyl]-methanone

Synonyms: Vanin 1-IN-1, Vanin 1 Inhibitor 1, VNN1-IN-1, VNN1 Inhibitor 1

MF: $C_{18}H_{22}N_6O_2$

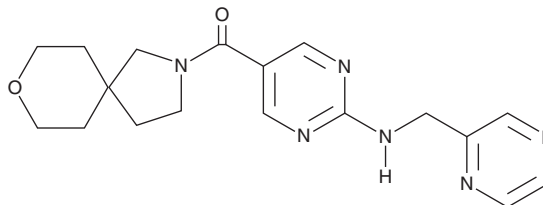
FW: 354.4

Purity: $\geq 98\%$

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

PFI-653 is supplied as a solid. A stock solution may be made by dissolving the PFI-653 in the solvent of choice, which should be purged with an inert gas. PFI-653 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

PFI-653 is an inhibitor of the pantetheinase vanin-1 ($IC_{50} = 6.85$ nM for the human enzyme).¹ *In vivo*, PFI-653 (50 mg/kg) prevents reductions in body weight, as well as reduces disease severity and serum vanin-1 levels in a mouse model of colitis induced by dextran sulfate sodium (DSS; Item No. 23250). See the Structural Genomics Consortium (SGC) website for more information.

Reference

1. Casimiro-Garcia, A., Allais, C., Brennan, A., *et al.* Discovery of a series of pyrimidine carboxamides as inhibitors of vanin-1. *J. Med. Chem.* **65**(1), 757-784 (2022).

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, 05/30/2024

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