

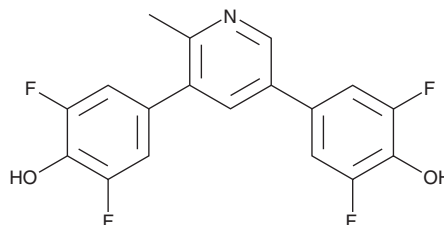
PRODUCT INFORMATION



VRK-IN-1

Item No. 44649

CAS Registry No.: 2378855-09-3
Formal Name: 4,4'-(2-methylpyridine-3,5-diyl)bis(2,6-difluorophenol)
Synonyms: Vaccinia-related Kinase 1-IN-1, Vaccinia-related Kinase-1 Inhibitor 1, VRK Inhibitor 1
MF: C₁₈H₁₁F₄NO₂
FW: 349.3
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

VRK-IN-1 is supplied as a solid. A stock solution may be made by dissolving the VRK-IN-1 in water. VRK-IN-1 is slightly soluble (0.1-1 mg/ml) in water. We do not recommend storing the aqueous solution for more than one day.

Description

VRK-IN-1 is an inhibitor of vaccinia-related kinase 1 (VRK1; IC₅₀ = 150 nM).¹ It is selective for VRK1 over a panel of 48 additional kinases at 1 μM.

Reference

1. Serafim, R.A.M., de Souza Gama, F.H., Dutra, L.A., *et al.* Development of pyridine-based inhibitors for the human vaccinia-related kinases 1 and 2. *ACS Med. Chem. Lett.* **10(9)**, 1266-1271 (2019).

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, 12/10/2025

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