

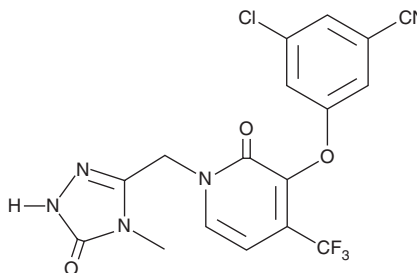
PRODUCT INFORMATION



Doravirine

Item No. 26405

CAS Registry No.: 1338225-97-0
Formal Name: 3-chloro-5-[[1-[(4,5-dihydro-4-methyl-5-oxo-1H-1,2,4-triazol-3-yl)methyl]-1,2-dihydro-2-oxo-4-(trifluoromethyl)-3-pyridinyl]oxy]-benzonitrile
Synonym: MK-1439
MF: C₁₇H₁₁ClF₃N₅O₃
FW: 425.8
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

Doravirine is supplied as a solid. A stock solution may be made by dissolving the doravirine in the solvent of choice, which should be purged with an inert gas. Doravirine is soluble in the organic solvent DMSO. Doravirine is slightly soluble in methanol.

Description

Doravirine is a non-nucleoside reverse transcriptase inhibitor (NNRTI) that inhibits HIV reverse transcriptase (IC₅₀ = 11 nM).¹ It inhibits the spread of wild-type HIV and K103N, Y181C, and K103N/Y181C mutant HIV strains in cells cultured with 50% normal human serum (IC₉₅s = 19, 42, 25, and 54 nM, respectively). Formulations containing doravirine have been used in the treatment of HIV-1 infection.

Reference

1. Côté, B., Burch, J.D., Asante-Appiah, E., et al. Discovery of MK-1439, an orally bioavailable non-nucleoside reverse transcriptase inhibitor potent against a wide range of resistant mutant HIV viruses. *Bioorg. Med. Chem. Lett.* **24(3)**, 917-922 (2014).

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, 12/07/2022

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