

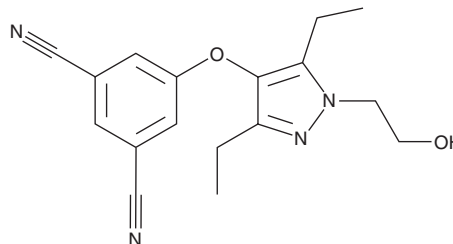
PRODUCT INFORMATION



Lersivirine

Item No. 29497

CAS Registry No.: 473921-12-9
Formal Name: 5-[[3,5-diethyl-1-(2-hydroxyethyl)-1H-pyrazol-4-yl]oxy]-1,3-benzenedicarbonitrile
Synonym: UK-453,061
MF: C₁₇H₁₈N₄O₂
FW: 310.4
Purity: ≥98%
UV/Vis.: λ_{max}: 213 nm
Supplied as: A crystalline 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

Lersivirine is supplied as a crystalline solid. A stock solution may be made by dissolving the lersivirine in the solvent of choice, which should be purged with an inert gas. Lersivirine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of lersivirine in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Description

Lersivirine is a non-nucleoside reverse transcriptase inhibitor (NNRTI; mean IC₅₀ = 118 nM for the wild-type HIV-1 enzyme).¹ It inhibits a panel of 14 single-point mutant HIV-1 reverse transcriptases with mean IC₅₀ values within 10-fold of the wild-type enzyme. Lersivirine inhibits HIV-1 virus replication in MT-2 cells with EC₅₀ values of 5.04, 13.1, and 34.9 nM at multiplicities of infection (MOIs) of 0.005, 0.05, and 0.5, respectively, and exhibits synergistic or additive effects when used in combination with a variety of antiretroviral agents.

Reference

1. Corbau, R., Mori, J., Phillips, C., *et al.* Lersivirine, a nonnucleoside reverse transcriptase inhibitor with activity against drug-resistant human immunodeficiency virus type 1. *Antimicrob. Agents Chemother.* **54**(10), 4451-4463 (2010).

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, 11/03/2022

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