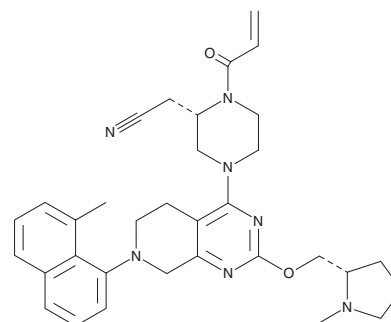


PRODUCT INFORMATION



MRTX-1257
Item No. 33527

CAS Registry No.: 2206736-04-9
Formal Name: (2S)-1-(1-oxo-2-propen-1-yl)-4-[5,6,7,8-tetrahydro-7-(8-methyl-1-naphthalenyl)-2-[[[(2S)-1-methyl-2-pyrrolidinyl]methoxy]pyrido[3,4-d]pyrimidin-4-yl]-2-piperazineacetonitrile
MF: C₃₃H₃₉N₇O₂
FW: 565.7
Purity: ≥98%
UV/Vis.: λ_{max}: 226 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

MRTX-1257 is supplied as a crystalline solid. A stock solution may be made by dissolving the MRTX-1257 in the solvent of choice, which should be purged with an inert gas. MRTX-1257 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of MRTX-1257 in these solvents is approximately 10 mg/ml in ethanol and approximately 30 mg/ml in DMSO and DMF.

Description

MRTX-1257 is a covalent inhibitor of K-Ras^{G12C}.¹ It inhibits proliferation of NCI H358 cells when used at concentrations ranging from 0.001 to 1 μM.

Reference

1. Kettle, J.G. and Cassar, D.J. Covalent inhibitors of the GTPase KRAS^{G12C}: A review of the patent literature. *Expert. Opin. Ther. Pat.* **30(2)**, 103-120 (2020).

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, 01/03/2023

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