

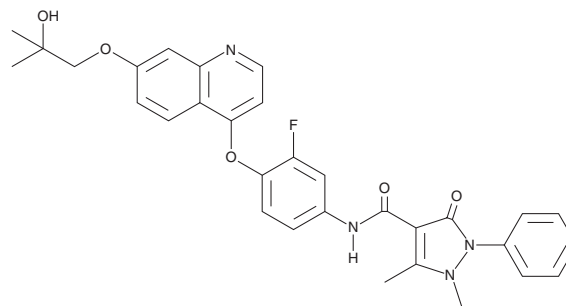
PRODUCT INFORMATION



Ningetinib

Item No. 25233

CAS Registry No.: 1394820-69-9
Formal Name: N-[3-fluoro-4-[[7-(2-hydroxy-2-methylpropoxy)-4-quinolinyl]oxy]phenyl]-2,3-dihydro-1,5-dimethyl-3-oxo-2-phenyl-1H-pyrazole-4-carboxamide
MF: C₃₁H₂₉FN₄O₅
FW: 556.6
Purity: ≥98%
UV/Vis.: λ_{max}: 234, 294 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

Ningetinib is supplied as a crystalline solid. A stock solution may be made by dissolving the ningetinib in the solvent of choice, which should be purged with an inert gas. Ningetinib is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of ningetinib in these solvents is approximately 2 and 15 mg/ml, respectively.

Ningetinib is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ningetinib should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Ningetinib has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Ningetinib is a multi-kinase inhibitor that inhibits c-Met, VEGFR2/KDR, and Axl (IC₅₀s = 19, 37, and 11 nM, respectively).¹ *In vivo*, ningetinib (20 mg/kg) inhibits tumor growth in MKN45, Caki-1, NCI-H441, Huh-7, U87MG, and MDA-MB-231 mouse xenograft models by 53-97%.

Reference

1. Xi, N. Substituted quinoline compounds and methods of use. *Sunshine Lake Pharma Co. Ltd. US20120219522A1*, (2012).

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, 11/02/2022

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