

PRODUCT INFORMATION

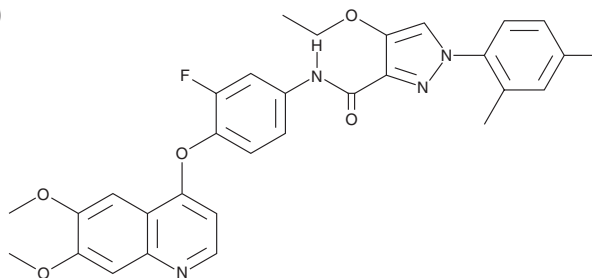


LDC-1267

Item No. 29459

CAS Registry No.: 1361030-48-9
Formal Name: N-[4-[(6,7-dimethoxy-4-quinolinyl)oxy]-3-fluorophenyl]-4-ethoxy-1-(4-fluoro-2-methylphenyl)-1H-pyrazole-3-carboxamide

MF: C₃₀H₂₆F₂N₄O₅
FW: 560.6
Purity: ≥98%
UV/Vis.: λ_{max}: 236, 294, 322 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

LDC-1267 is supplied as a crystalline solid. A stock solution may be made by dissolving the LDC-1267 in the solvent of choice, which should be purged with an inert gas. LDC-1267 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of LDC-1267 in these solvents is approximately 1 and 3 mg/ml, respectively. LDC-1267 is also slightly soluble in ethanol.

LDC-1267 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, LDC-1267 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. LDC-1267 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

LDC-1267 is a TAM family kinase inhibitor (IC₅₀s = 5, 8, and 29 nM for Mer, Tyro3, and Axl, respectively).¹ It is selective for TAM family kinases over a panel of 456 kinases at 1 μM, but does inhibit Met, Aurora B, LCK, and Src (IC₅₀s = 35, 36, 51, and 338 nM, respectively). It inhibits suppression of proliferation and IFN-γ production induced by growth arrest-specific protein 6 (GAS6) in C-type lectin-like receptor NKG2D-activated natural killer (NK) cells. Adoptive transfer of NK cells pre-incubated with LDC-1267 decreases tumor growth but not metastasis in a B16/F10 murine melanoma model. LDC-1267 (20 mg/kg, i.p.) reduces the number of metastases in B16/F10 murine melanoma and 4T1 murine mammary carcinoma models in an NK cell-dependent manner.

Reference

1. Paolino, M., Choidas, A., Wallner, S., *et al.* The E3 ligase Cbl-b and TAM receptors regulate cancer metastasis via natural killer cells. *Nature* **507**(7493), 508-512 (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, 10/20/2022

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