

PRODUCT INFORMATION

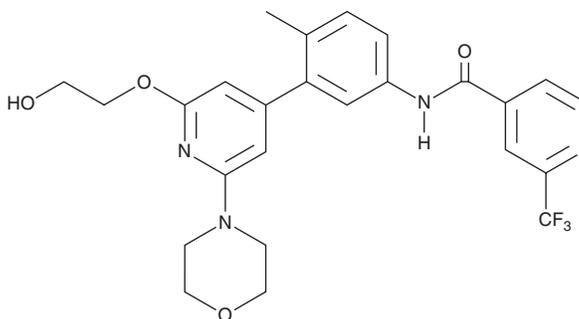


LXH254

Item No. 28231

CAS Registry No.: 1800398-38-2
Formal Name: N-[3-[2-(2-hydroxyethoxy)-6-(4-morpholinyl)-4-pyridinyl]-4-methylphenyl]-2-(trifluoromethyl)-4-pyridinecarboxamide

MF: C₂₅H₂₅F₃N₄O₄
FW: 502.5
Purity: ≥98%
UV/Vis.: λ_{max}: 250, 305 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

LXH254 is supplied as a crystalline solid. A stock solution may be made by dissolving the LXH254 in the solvent of choice, which should be purged with an inert gas. LXH254 is soluble in the organic solvent DMSO.

Description

LXH254 is a pan-RAF inhibitor that selectively inhibits B-RAF and C-RAF over a panel of 456 human kinases.¹ It inhibits MAPK signaling in tumor cells expressing mutant B-RAF^{V600}. LXH254 induces tumor regression in various mouse xenograft models.

Reference

1. Stuart, D.D., Shao, W., Mishina, Y., *et al.* Abstract DDT01-04: Pharmacological profile and anti-tumor properties of LXH254, a highly selective RAF kinase inhibitor. *Cancer Res.* **78(Suppl. 13)** (2018).

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, 11/11/2022

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