

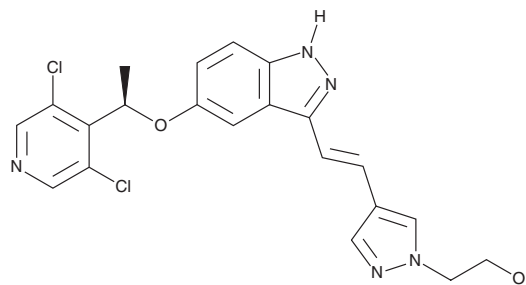
PRODUCT INFORMATION



LY2874455

Item No. 20843

CAS Registry No.: 1254473-64-7
Formal Name: 4-[(1E)-2-[5-[(1R)-1-(3,5-dichloro-4-pyridinyl)ethoxy]-1H-indazol-3-yl]ethenyl]-1H-pyrazole-1-ethanol
MF: C₂₁H₁₉Cl₂N₅O₂
FW: 444.3
Purity: ≥98%
UV/Vis.: λ_{max}: 222, 279, 333 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

LY2874455 is supplied as a crystalline solid. A stock solution may be made by dissolving the LY2874455 in the solvent of choice, which should be purged with an inert gas. LY2874455 is soluble in the organic solvent DMSO at a concentration of approximately 4 mg/ml.

Description

LY2874455 is a pan-inhibitor of fibroblast growth factor receptor (FGFR) tyrosine kinases with IC₅₀ values of 2.8, 2.6, 6.4, 6, and 7 nM for FGFR1, FGFR2, FGFR3, FGFR4, and VEGFR2, respectively.¹ It has been shown to inhibit proliferation of KMS-11, OPM-2, SNU-16, and KATO-III cancer cells *in vitro* and in tumor xenografts in mice.¹

Reference

1. Zhao, G., Li, W.-y., Chen, D., *et al.* A novel, selective inhibitor of fibroblast growth factor receptors that shows a potent broad spectrum of antitumor activity in several tumor xenograft models. *Mol. Cancer Ther.* **10(11)**, 2200-2210 (2011).

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, 12/01/2022

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