

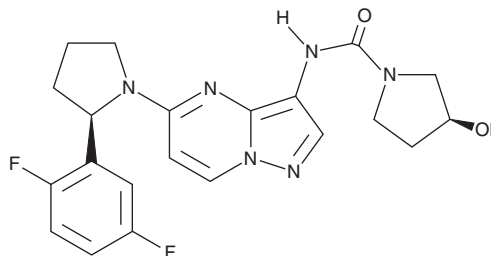
PRODUCT INFORMATION



LOXO-101

Item No. 27056

CAS Registry No.: 1223403-58-4
Formal Name: (3S)-N-[5-[(2R)-2-(2,5-difluorophenyl)-1-pyrrolidinyl]pyrazolo[1,5-a]pyrimidin-3-yl]-3-hydroxy-1-pyrrolidinecarboxamide
Synonyms: ARRY-470, Larotrectinib
MF: C₂₁H₂₂F₂N₆O₂
FW: 428.4
Purity: ≥98%
UV/Vis.: λ_{max}: 264 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

LOXO-101 is supplied as a crystalline solid. A stock solution may be made by dissolving the LOXO-101 in the solvent of choice, which should be purged with an inert gas. LOXO-101 is soluble in the organic solvent DMSO at a concentration of approximately 4 mg/ml.

Description

LOXO-101 is an inhibitor of the tropomyosin-related kinases TrkA, TrkB, and TrkC (IC₅₀s = 2-20 nM).¹ It is selective for TrkA, -B, and -C over a panel of 226 kinases at 1 μM. LOXO-101 inhibits the growth of CUTO-3.29, KM12, and MO-91 patient-derived cancer cell lines (IC₅₀s = <100, <10, and <10 nM, respectively). *In vivo*, LOXO-101 (60 and 200 mg/kg) reduces tumor growth in a KM12 mouse xenograft model.

Reference

1. Doebele, R.C., Davis, L.E., Vaishnavi, A., *et al.* An oncogenic NTRK fusion in a patient with soft-tissue sarcoma with response to the tropomyosin-related kinase inhibitor LOXO-101. *Cancer Discov.* **5**(10), 1049-1057 (2015).

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, 12/07/2022

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