

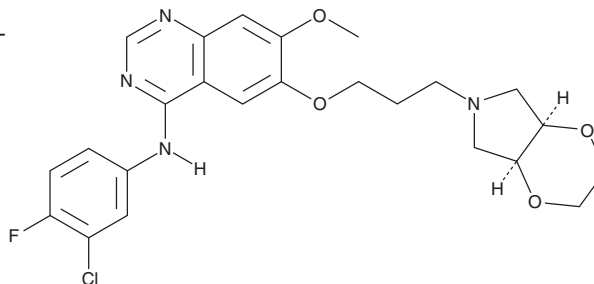
# PRODUCT INFORMATION



## Larotinib

Item No. 44444

**CAS Registry No.:** 1438072-11-7  
**Formal Name:** *rel*-N-(3-chloro-4-fluorophenyl)-6-[3-[(4aR,7aS)-hexahydro-6H-1,4-dioxino[2,3-c]pyrrol-6-yl]propoxy]-7-methoxy-4-quinazolinamine  
**Synonym:** NX-125  
**MF:** C<sub>24</sub>H<sub>26</sub>ClFN<sub>4</sub>O<sub>4</sub>  
**FW:** 488.9  
**Purity:** ≥98%  
**Supplied as:** A 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

Larotinib is supplied as a solid. A stock solution may be made by dissolving the larotinib in the solvent of choice, which should be purged with an inert gas. Larotinib is sparingly soluble (1-10 mg/ml) in DMSO.

### Description

Larotinib is a multi-kinase inhibitor.<sup>1</sup> It inhibits wild-type EGFR (IC<sub>50</sub> = 0.6 nM), mutant EGFRs, including EGFR<sup>L858R</sup> and EGFR<sup>L861Q</sup> (IC<sub>50</sub>s = 0.563 and 0.423 nM, respectively), HER2 (IC<sub>50</sub> = 253 nM), and HER4 (IC<sub>50</sub> = 84 nM). It also inhibits receptor-interacting serine/threonine kinase 2 (RIPK2), IL-1 receptor-associated kinase 1 (IRAK1), B-lymphoid kinase (Blk), and Bruton's tyrosine kinase (BTK; IC<sub>50</sub>s = 26.6, 167, 102, and 196 nM, respectively).

### Reference

1. Liu, J., Zhang, H., Zhu, X., *et al.* Phase I trial to evaluate the tolerance, pharmacokinetics and efficacy of the broad-spectrum ErbB family inhibitor larotinib mesylate in patients with advanced solid tumors. *Front. Pharmacol.* **12**, 636324 (2021).

#### 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, 12/19/2025

#### 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