

# PRODUCT INFORMATION



**XL647**

Item No. 27049

**CAS Registry No.:** 781613-23-8

**Formal Name:** N-(3,4-dichloro-2-fluorophenyl)-6-methoxy-7-[[[(3 $\alpha$ ,5 $\beta$ ,6 $\alpha$ )-octahydro-2-methylcyclopenta[c]pyrrol-5-yl]methoxy]-4-quinazolinamine

**Synonyms:** EXEL-7647, Tesevatinib

**MF:** C<sub>24</sub>H<sub>25</sub>Cl<sub>2</sub>FN<sub>4</sub>O<sub>2</sub>

**FW:** 491.4

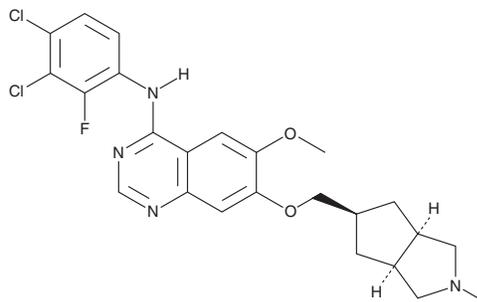
**Purity:**  $\geq$ 98%

**UV/Vis.:**  $\lambda_{\text{max}}$ : 233, 250, 331 nm

**Supplied as:** A solid

**Storage:** -20°C

**Stability:**  $\geq$ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

## Laboratory Procedures

XL647 is supplied as a solid. A stock solution may be made by dissolving the XL647 in the solvent of choice, which should be purged with an inert gas. XL647 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of XL647 in these solvents is approximately 10 mg/ml.

## Description

XL647 is a multi-kinase inhibitor ( $IC_{50}$ s = 0.3, 16, 1.5, 8.7, and 1.4 nM for EGFR, ErbB2, KDR, FLT4, and EphB4, respectively).<sup>1</sup> It is selective for these kinases over a panel of 10 tyrosine kinases and 55 serine/threonine kinases at 10  $\mu$ M. XL647 inhibits growth of A431 cells expressing wild-type EGFR and H1975 non-small cell lung cancer (NSCLC) cells expressing both the activating mutant EGFR<sup>L858R</sup> and the drug resistance-associated mutant EGFR<sup>T790M</sup> ( $IC_{50}$ s = 13 and 920 nM, respectively). *In vivo*, XL647 (10, 30, and 100 mg/kg) inhibits tumor growth and EGFR phosphorylation in an H1975 mouse xenograft model in a dose-dependent manner.

## Reference

1. Gendreau, S.B., Ventura, R., Keast, P., *et al.* Inhibition of the T790M gatekeeper mutant of the epidermal growth factor receptor by EXEL-7647. *Clin. Cancer Res.* **13**(12), 3713-3723 (2007).

### 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, 12/19/2022

## 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