

# PRODUCT INFORMATION

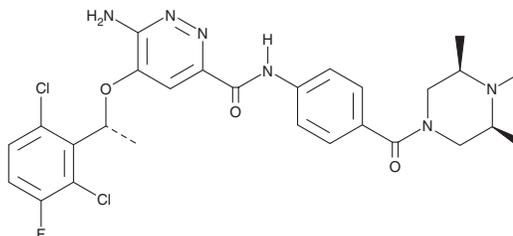


## X-396

Item No. 44938

**CAS Registry No.:** 1370651-20-9  
**Formal Name:** 6-amino-5-[(1R)-1-(2,6-dichloro-3-fluorophenyl)ethoxy]-N-[4-[[[(3R,5S)-3,5-dimethyl-1-piperazinyl]carbonyl]phenyl]-3-pyridinecarboxamide

**Synonym:** Ensartinib  
**MF:** C<sub>26</sub>H<sub>27</sub>Cl<sub>2</sub>FN<sub>6</sub>O<sub>3</sub>  
**FW:** 561.4  
**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

X-396 is supplied as a solid. A stock solution may be made by dissolving the X-396 in the solvent of choice, which should be purged with an inert gas. X-396 is slightly soluble (0.1-1 mg/ml) in DMSO.

### Description

X-396 is an inhibitor of anaplastic lymphoma kinase (ALK) and MET (IC<sub>50</sub>s = <0.4 and 0.74 nM, respectively).<sup>1</sup> It selectively inhibits the growth of cancer cell lines harboring gain-of-function ALK-fusion proteins (IC<sub>50</sub>s = 9-106 nM) over those with cell growth driven by other mutant kinases (IC<sub>50</sub>s = 757->3,000 nM). X-396 (25 mg/kg) reduces tumor growth in a lung cancer mouse xenograft model using NCI H3122 cells, which express a gain-of-function EML4-ALK fusion protein.

### Reference

1. Lovly, C.M., Heuckmann, J.M., de Stanchina, E., *et al.* Insights into ALK-driven cancers revealed through development of novel ALK tyrosine kinase inhibitors. *Cancer Res.* **71(14)**, 4920-4931 (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, 01/22/2026

#### 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