

# PRODUCT INFORMATION

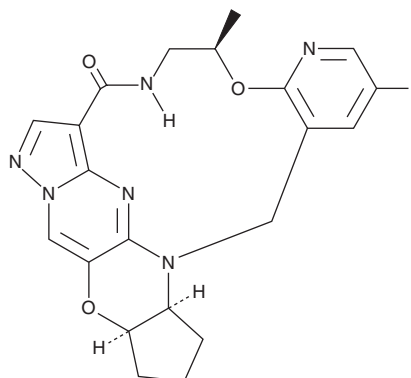


**TPX-0046**

Item No. 41990

**CAS Registry No.:** 2359650-19-2  
**Formal Name:** (3aS,11R,20aR)-7-fluoro-2,3,3a,12,13,20a-hexahydro-11-methyl-17,19-metheno-1H,5H-cyclopent[5,6][1,4]oxazino[3,4-i]pyrazolo[4,3-f]pyrido[3,2-l][1,4,8,10]oxatriazacyclotridecin-14(11H)-one

**Synonym:** Enbezotinib  
**MF:** C<sub>21</sub>H<sub>21</sub>FN<sub>6</sub>O<sub>3</sub>  
**FW:** 424.4  
**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

TPX-0046 is supplied as a solid. A stock solution may be made by dissolving the TPX-0046 in the solvent of choice, which should be purged with an inert gas. TPX-0046 slightly soluble (0.1-1 mg/ml) in DMSO.

## Description

TPX-0046 is a RET inhibitor ( $IC_{50}$  = 0.26 nM for the wild-type enzyme).<sup>1</sup> It is selective for RET over a panel of 217 additional kinases but does inhibit 39 tyrosine kinases by greater than 50% at 26 nM. TPX-0046 inhibits the phosphorylation of wild-type RET ( $IC_{50}$  = 21.9 nM) and 15 RET mutants ( $IC_{50}$ s = 2.69-108 nM) in Ba/F3 cells.

## Reference

1. Miyazaki, I., Odintsov, I., Ishida, K., *et al.* Vepafestinib is a pharmacologically advanced RET-selective inhibitor with high CNS penetration and inhibitory activity against RET solvent front mutations. *Nat. Cancer* **4**(9), 1345-1361 (2023).

### 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, 11/11/2024

## 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