

PRODUCT INFORMATION



TPX-0022

Item No. 36118

CAS Registry No.: 2271119-26-5

Formal Name: (7S)-3-amino-14-ethyl-11-fluoro-4,5,6,7,13,14-hexahydro-7-methyl-4-oxo-1,15-etheno-1H-pyrazolo[4,3-f][1,4,8,10]benzoxatriazacyclotridecine-12-carbonitrile
CSF1R-IN-2, CSF1R Inhibitor 2, Elzovantinib

Synonyms:

MF: C₂₀H₂₀FN₇O₂

FW: 409.4

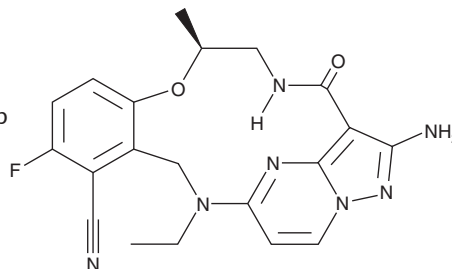
Purity: ≥98%

UV/Vis.: λ_{max}: 215, 271 nm

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-0022 is supplied as a solid. A stock solution may be made by dissolving the TPX-0022 in the solvent of choice, which should be purged with an inert gas. TPX-0022 is soluble in the organic solvent ethanol at a concentration of approximately 5 mg/ml.

Description

TPX-0022 is an inhibitor of Src, MET, and CSF-1 receptor tyrosine kinase (FMS) kinases (IC₅₀s = 0.12, 0.14, and 0.76 nM, respectively).¹ It inhibits the proliferation of MKN45 and SNU-5 human gastric cancer cells (IC₅₀s = 0.2 and 0.17 nM, respectively), as well as Ba/F3 pro-B cells expressing TEL-CSF1R (IC₅₀ = 19.3 nM). TPX-0022 (100 nM) inhibits cell migration of HCC827 human lung cancer cells. It inhibits tumor growth in an MKN45 mouse xenograft model when administered at doses of 10 and 30 mg/kg. TPX-0022 (10 mg/kg) also inhibits phosphorylation of MET at Tyr¹²³⁴ and Tyr¹³⁴⁹ in the same model.

Reference

1. Cui, J.J., Rogers, E.W., Ung, J., *et al.* Macrocyclic compounds and uses thereof. *TP Therapeutics, Inc.* WO2019/023417A1 (2019).

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, 10/26/2022

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