

PRODUCT INFORMATION



CEP-28122

Item No. 18406

CAS Registry No.: 1022958-60-6
Formal Name: (1S,2S,3R,4R)-3-[[5-chloro-2-[[[(7S)-6,7,8,9-tetrahydro-1-methoxy-7-(4-morpholinyl)-5H-benzocyclohepten-2-yl]amino]-4-pyrimidinyl]amino]-bicyclo[2.2.1]hept-5-ene-2-carboxamide

MF: C₂₈H₃₅ClN₆O₃

FW: 539.1

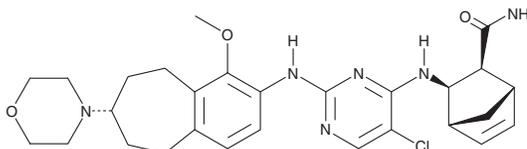
Purity: ≥98%

UV/Vis.: λ_{max}: 274 nm

Supplied as: A crystalline 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

CEP-28122 is supplied as a crystalline solid. A stock solution may be made by dissolving the CEP-28122 in the solvent of choice, which should be purged with an inert gas. CEP-28122 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of CEP-28122 in these solvents is approximately 30 and 12 mg/ml, respectively.

CEP-28122 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CEP-28122 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CEP-28122 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

CEP-28122 is a potent inhibitor of anaplastic lymphoma kinase (ALK; IC₅₀ = 1.9 nM) that less potently inhibits other tyrosine and serine/threonine kinases.¹ It blocks the phosphorylation of ALK and other ALK substrates in cells and induces concentration-dependent growth inhibition or cytotoxicity of ALK-positive cancer cells.¹ CEP-28122 is orally active, inhibiting ALK tyrosine phosphorylation in tumor xenografts in mice for more than 12 hours following single oral dosing at 30 mg/kg.¹

Reference

1. Chang, M., Quail, M.R., Gingrich, D.E., *et al.* CEP-28122, a highly potent and selective orally active inhibitor of anaplastic lymphoma kinase with antitumor activity in experimental models of human cancers. *Mol. Cancer Ther.* **11**(3), 670-679 (2012).

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, 11/18/2022

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