

PRODUCT INFORMATION

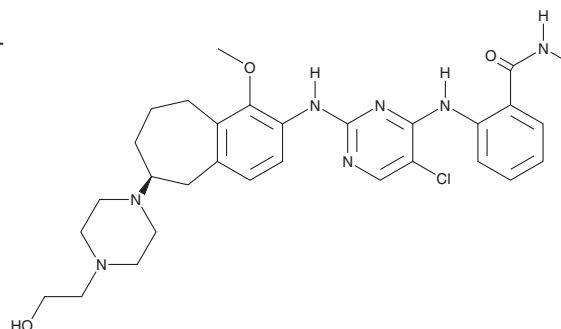


CEP-37440

Item No. 21278

CAS Registry No.: 1391712-60-9
Formal Name: 2-[[5-chloro-2-[[[(6S)-6,7,8,9-tetrahydro-6-[4-(2-hydroxyethyl)-1-piperazinyl]-1-methoxy-5H-benzocyclohepten-2-yl]amino]-4-pyrimidinyl]amino]-N-methyl-benzamide

MF: C₃₀H₃₈ClN₇O₃
FW: 580.1
Purity: ≥98%
UV/Vis.: λ_{max}: 279, 333 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-37440 is supplied as a crystalline solid. A stock solution may be made by dissolving the CEP-37440 in the solvent of choice, which should be purged with an inert gas. CEP-37440 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of CEP-37440 in these solvents is approximately 30 and 12 mg/ml, respectively.

CEP-37440 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, CEP-37440 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. CEP-37440 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-37440 is a dual inhibitor of focal adhesion kinase 1 (FAK1; IC₅₀s = 2.0 and 80 nM for *in vitro* enzyme and cellular inhibition, respectively) and anaplastic lymphoma kinase (ALK; IC₅₀s = 3.1 and 22 nM for *in vitro* enzyme and cellular inhibition, respectively).^{1,2} It blocks the proliferation of inflammatory and triple-negative breast cancer cell lines but not normal breast epithelial cells.² CEP-37440 reduces the growth and metastasis of breast cancer xenografts in mice when given by oral gavage.²

References

- Ott, G.R., Cheng, M., Learn, K.S., *et al.* Discovery of clinical candidate CEP-37440, a selective inhibitor of focal adhesion kinase (FAK) and anaplastic lymphoma kinase (ALK). *J. Med. Chem.* **59(16)**, 7478-7496 (2016).
- Salem, I., Alsalahi, M., Chervoneva, I., *et al.* The effects of CEP-37440, an inhibitor of focal adhesion kinase, *in vitro* and *in vivo* on inflammatory breast cancer cells. *Breast Cancer Res.* **18(1)**, 37 (2016).

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, 11/29/2022

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