

PRODUCT INFORMATION



Lestaurtinib

Item No. 12094

CAS Registry No.: 111358-88-4
Formal Name: 2,3,9S,10S,11,12R-hexahydro-10-hydroxy-10-(hydroxymethyl)-9-methyl-9,12-epoxy-1H-diindolo[1,2,3-fg:3',2',1'-kl]pyrrolo[3,4-i][1,6]benzodiazocin-1-one

Synonyms: A 154475.0, CEP-701, KT 5555, NSC 621867, SP 924

MF: C₂₆H₂₁N₃O₄

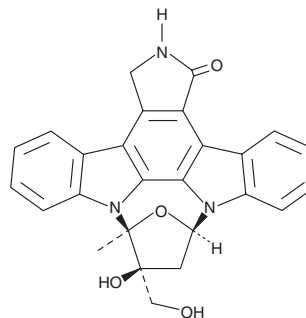
FW: 439.5

Purity: ≥98%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Lestaurtinib is supplied as a crystalline solid. A stock solution may be made by dissolving the lestaurtinib in the solvent of choice. Lestaurtinib is soluble in organic solvents such as DMSO and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of lestaurtinib in these solvents is approximately 3 and 5 mg/ml, respectively.

Lestaurtinib is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, lestaurtinib should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Lestaurtinib has a solubility of approximately 0.05 mg/ml in a 1:20 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Janus-associated kinases (JAKs) are cytoplasmic tyrosine kinases that are required for activating the signaling of certain cytokines and growth factor receptors.^{1,2} A JAK2 gene fusion mutation, JAK2^{V617F}, that causes unchecked activation of various growth factors and cytokines, has been linked to myeloproliferative disorders.³ Lestaurtinib is a staurosporine analog that potently inhibits JAK2 kinase (IC₅₀ = 1 nM) and downstream targets STAT5 (IC₅₀ = 10-30 nM) and STAT3 in a human erythroleukemic cell line expressing the JAK2^{V617F} mutation.^{3,4} It has been used to reduce tumor growth in a xenograft model of pancreatic ductal adenocarcinoma, to inhibit cell growth in a trk-B overexpressing neuroblastoma xenograft model, and to block proliferation and induce apoptosis in Hodgkin lymphoma cell lines.⁵⁻⁷ Lestaurtinib has since been reported to potently inhibit the epigenetic kinase PRK1 *in vitro* (IC₅₀ = 8.6 nM), inducing hypophosphorylation of histone H3 threonine 11 and blocking androgen-dependent gene expression in prostate cancer cells at a concentration of 5 μM.⁸

References

1. Leonard, W.J. and O'Shea, J.J. *Annu. Rev. Immunol.* **16**, 293-322 (1998).
2. Yamaoka, K., Saharinen, P., Pesu, M., et al. *Genome Biol.* **5**(12), 1-6 (2004).
3. Verstovsek, S. *Hematology Am. Soc. Hematol. Educ. Program* **2009**(1), 636-642 (2009).
4. Hexner, E.O., Serdikoff, C., Jan, M., et al. *Blood* **111**(12), 5663-5671 (2008).
5. Miknyoczki, S.J., Chang, H., Klein-Szanto, A., et al. *Clin. Cancer Res.* **5**, 2205-2212 (1999).
6. Iyer, R., Evans, A.E., Qi, X., et al. *Clin. Cancer Res.* **16**(5), 1478-1485 (2010).
7. Diaz, T., Navarro, A., Ferrer, G., et al. *PLoS One* **6**(4), (2011).
8. Köhler, J., Erlenkamp, G., Eberlin, A., et al. *PLoS One* **7**(4), (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, 05/02/2023

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