

PRODUCT INFORMATION



Lonafarnib

Item No. 11746

CAS Registry No.: 193275-84-2
Formal Name: 4-[2-[4-[(11R)-3,10-dibromo-8-chloro-6,11-dihydro-5H-benzo[5,6]cyclohepta[1,2-b]pyridin-11-yl]-1-piperidinyl]-2-oxoethyl]-1-piperidinecarboxamide

Synonyms: Sarasar, SCH 66336

MF: C₂₇H₃₁Br₂ClN₄O₂

FW: 638.8

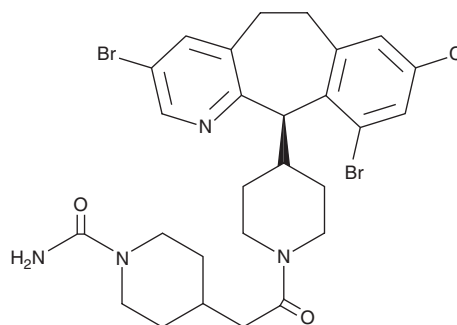
Purity: ≥98%

UV/Vis.: λ_{max}: 278 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

Lonafarnib is supplied as a crystalline solid. A stock solution may be made by dissolving the lonafarnib in the solvent of choice, which should be purged with an inert gas. Lonafarnib is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of lonafarnib in ethanol and DMF is approximately 14 mg/ml and approximately 3 mg/ml in DMSO.

Lonafarnib is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, lonafarnib should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Lonafarnib has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Farnesyltransferase drives the post-translational farnesylation of a number of target proteins, including Ras, a component of the MAPK and Akt signaling pathways involved in cell proliferation, and Ras homolog enriched in brain (Rheb), which stimulates mTOR activity.¹ By arresting growth, farnesyltransferase inhibitors have great potential as antitumor agents. Lonafarnib is a farnesyltransferase inhibitor that blocks the post-translational lipid modification of oncogenic Ras isoforms H-Ras, N-Ras, and K-Ras (IC₅₀s = 1.9, 2.8, and 5.2 nM, respectively) as well as Rheb (IC₅₀ = 10-100 nM).^{2,3} It inhibits H-Ras signaling in whole cells (IC₅₀ = 10 nM) and blocks the transformed growth properties of fibroblasts (IC₅₀ = 75 nM) and human tumor cell lines (IC₅₀ = 400 nM) expressing activated K-Ras proteins. In the nude mouse, lonafarnib demonstrates potent dose-dependent oral activity in an array of human tumor xenograft models including tumors originating from colon, lung, pancreas, prostate, and urinary bladder.² Combination treatments with various cytotoxic agents (cyclophosphamide, 5-fluorouracil, and vincristine) or MAPK or Akt pathway inhibitors can enhance the *in vivo* efficacy of lonafarnib.^{2,4}

References

1. Appels, N.M.G.M., Beijnen, J.H., and Schellens, J.H.M. *Oncologist* **10**(8), 565-578 (2005).
2. Liu, M., Bryant, M.S., Chen, J., et al. *Cancer Res.* **58**(21), 4947-4956 (1998).
3. Basso, A.D., Mirza, A., Liu, G., et al. *J. Biol. Chem.* **280**(35), 31101-31108 (2005).
4. Niessner, H., Beck, D., Sinnberg, T., et al. *J. Invest. Dermatol.* **131**(2), 468-479 (2011).

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, 12/15/2022

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