

PRODUCT INFORMATION

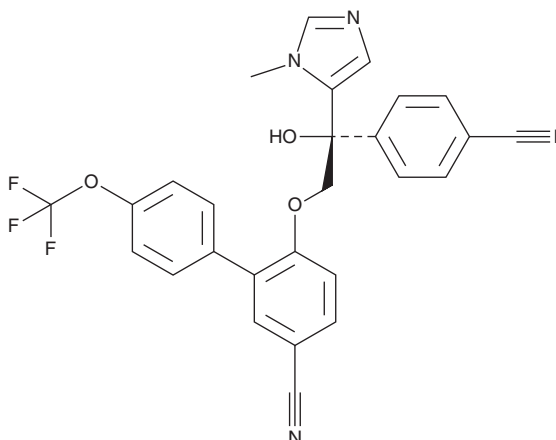


ABT-100

Item No. 41355

CAS Registry No.: 450839-40-4
Formal Name: 6-[(2S)-2-(4-cyanophenyl)-2-hydroxy-2-(1-methyl-1H-imidazol-5-yl)ethoxy]-4'-(trifluoromethoxy)-[1,1'-biphenyl]-3-carbonitrile

Synonym: A-367074
MF: C₂₇H₁₉F₃N₄O₃
FW: 504.5
Purity: ≥98%
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

ABT-100 is supplied as a solid. A stock solution may be made by dissolving the ABT-100 in the solvent of choice, which should be purged with an inert gas. ABT-100 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

ABT-100 is an inhibitor of farnesyltransferase (FTase).¹ It inhibits Ras farnesylation but not Rap1A geranylgeranylation in HepG2 and Huh7 cells when used at a concentration of 20 μM. ABT-100 (2-20 μM) inhibits the proliferation of the same cells. *In vivo*, ABT-100 (39 mg/kg per day) increases body weight and percent survival, improves grip strength, and decreases the number of rib fractures in a *Zmpste24*^{-/-} mouse model of progeria.²

References

1. Carloni, V., Vizzutti, F., and Pantaleo, P. Farnesyltransferase inhibitor, ABT-100, is a potent liver cancer chemopreventive agent. *Clin. Cancer Res.* **11**(11), 4266-4274 (2005).
2. Fong, L.G., Frost, D., Meta, M., et al. A protein farnesyltransferase inhibitor ameliorates disease in a mouse model of progeria. *Science* **311**, 1621-1623 (2006).

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, 07/26/2024

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