

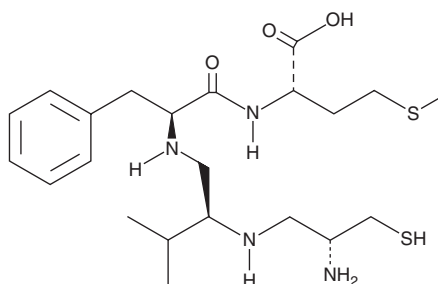
PRODUCT INFORMATION



FTase Inhibitor I

Item No. 22755

CAS Registry No.: 149759-96-6
Formal Name: N-[(2S)-2-[[[(2R)-2-amino-3-mercaptopropyl]amino]-3-methylbutyl]-L-phenylalanyl-L-methionine
Synonym: Farnesyltransferase Inhibitor I
MF: C₂₂H₃₈N₄O₃S₂
FW: 470.7
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Special Conditions: Hygroscopic



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

Laboratory Procedures

FTase inhibitor I is supplied as a solid. A stock solution may be made by dissolving the FTase inhibitor I in water. We do not recommend storing the aqueous solution for more than one day.

Description

FTase inhibitor I is a potent inhibitor of farnesyltransferase (FTase; IC₅₀ = 21 nM) that is greater than 30-fold selective for FTase over geranylgeranyl transferase (GGTase; IC₅₀ = 790 nM).¹ It prevents farnesylation of Ras and inhibits proliferation of cells transformed by Ras farnesylation.^{1,2} In a dog model of subarachnoid hemorrhage, it reduces GTP-Ras in spastic basilar arteries, decreases angiographic vasospasm, and improves clinical scores.³

References

1. Garcia, A.M., Rowell, C., Ackermann, K., *et al.* Peptidomimetic inhibitors of Ras farnesylation and function in whole cells. *J. Biol. Chem.* **268(25)**, 18415-18418 (1993).
2. Cox, A.D., Garcia, A.M., Westwick, J.K., *et al.* The CAAX peptidomimetic compound B581 specifically blocks farnesylation, but not geranylgeranylated or myristylated, oncogenic ras signaling and transformation. *J. Biol. Chem.* **269(30)**, 19203-19206 (1994).
3. Yamaguchi, M., Zhou, C., Nanda, A., *et al.* Ras protein contributes to cerebral vasospasm in a canine double-hemorrhage model. *Stroke* **35(7)**, 1750-1755 (2004).

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, 11/28/2022

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