

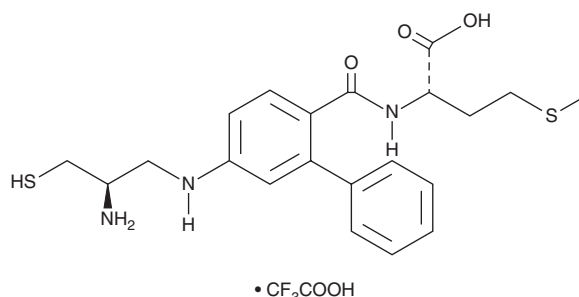
PRODUCT INFORMATION



FTI-276 (trifluoroacetate salt)

Item No. 10011661

CAS Registry No.: 1217471-51-6
Formal Name: N-[[5-[[[(2R)-2-amino-3-mercaptopropyl]amino][1,1'-biphenyl]-2-yl]carbonyl]-L-methionine, 2,2,2-trifluoroacetate
Synonym: Farnesyltransferase Inhibitor 276
MF: C₂₁H₂₇N₃O₃S₂ • CF₃COOH
FW: 547.6
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

FTI-276 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the FTI-276 (trifluoroacetate salt) in water. The solubility of FTI-276 (trifluoroacetate salt) in water is approximately 10 mM. We do not recommend storing the aqueous solution for more than one day.

Description

FTI-276 is a CAAX peptidomimetic inhibitor of farnesyltransferase (FTase; IC₅₀ = 500 pM).¹ It contains the C-terminal CAAX sequence, Cys-Val-Iso-Met (CVIM), of K-Ras4B, a form of Ras that is commonly mutated in human cancers, and is selective for FTase over geranylgeranyltransferase I (GGTase I; IC₅₀ = 50 nM). It inhibits the growth of NIH3T3 cells expressing oncogenic farnesylated H-Ras (RasF), but not geranylgeranylated H-Ras (RasGG) or Raf, when used at a concentration of 20 μM.² FTI-276 (50 mg/kg) also reduces tumor growth in a mouse xenograft model using Calu-1 cells that contain a K-Ras oncogenic mutation but not in a model using NCI H810 cells that do not contain a K-Ras mutation.

References

1. Lerner, E.C., Qian, Y., Blaskovich, M.A., *et al.* Ras CAAX peptidomimetic FTI-277 selectively blocks oncogenic Ras signaling by inducing cytoplasmic accumulation of inactive Ras-Raf complexes. *J. Biol. Chem.* **270**(45), 26802-26806 (1995).
2. Sun, J., Qian, Y., Hamilton, A.D., *et al.* Ras CAAX peptidomimetic FTI 276 selectively blocks tumor growth in nude mice of a human lung carcinoma with K-Ras mutation and p53 deletion. *Cancer Res.* **55**(19), 4243-4247 (1995).

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, 04/10/2023

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