

# PRODUCT INFORMATION



## FTI-277 (hydrochloride)

Item No. 10011667

**CAS Registry No.:** 180977-34-8  
**Formal Name:** N-[[5-[[[(2R)-2-amino-3-mercaptopropyl]amino][1,1'-biphenyl]-2-yl]carbonyl]-L-methionine, methyl ester, hydrochloride

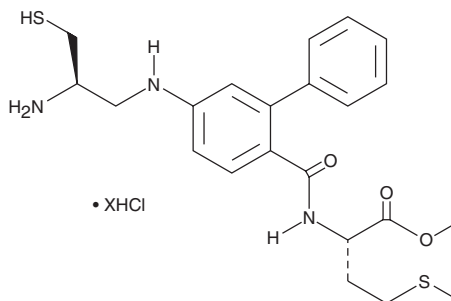
**Synonyms:** Farnesyltransferase Inhibitor 277, FTase Inhibitor 277

**MF:** C<sub>22</sub>H<sub>29</sub>N<sub>3</sub>O<sub>3</sub>S<sub>2</sub> • XHCl  
**FW:** 447.6

**UV/Vis.:** λ<sub>max</sub>: 240, 273 nm

**Storage:** -20°C

**Stability:** ≥4 years



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

### Description

FTI-277 is an inhibitor of farnesyltransferase (FTase; IC<sub>50</sub> = 0.1 μM).<sup>1</sup> It is selective for FTase over geranylgeranyl transferase (GGTase; IC<sub>50</sub> = >10 μM). FTI-277 (10 μM) inhibits IL-1β-induced production of nitric oxide (NO) in primary rat pulmonary artery smooth muscle cells.<sup>2</sup> It decreases the proliferation of A549 lung cancer cells when used at a concentration of 10 μM and induces cell cycle arrest at the G<sub>2</sub>/M phase in the same cells at 20 μM.<sup>3</sup> Intracisternal administration of FTI-277 increases the diameter of basilar arteries, improves appetite, and increases activity in a dog model of subarachnoid hemorrhage-induced vasospasm.<sup>4</sup>

### 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. Finder, J.D., Litz, J.L., Blaskovich, M.A., *et al.* Inhibition of protein geranylgeranylation causes a superinduction of nitric-oxide synthase-2 by interleukin-1β in vascular smooth muscle cells. *J. Biol. Chem.* **272**, 13484-13488 (1997).
3. Miquel, K., Pradines, A., Sun, J., *et al.* GGTI-298 induces G<sub>0</sub>-G<sub>1</sub> block and apoptosis whereas FTI-277 causes G<sub>2</sub>-M enrichment in A549 cells. *Cancer Res.* **57**, 1846-1850 (1997).
4. 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, 05/29/2025

#### 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