

PRODUCT INFORMATION

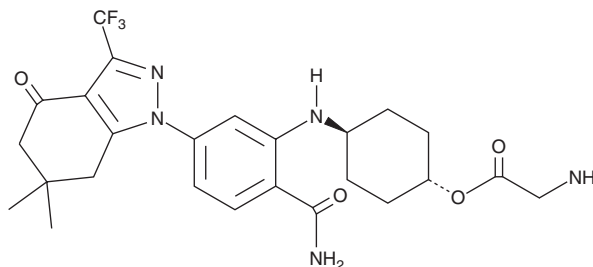


PF-04929113

Item No. 18270

CAS Registry No.: 908115-27-5
Formal Name: glycine, *trans*-4-[[2-(aminocarbonyl)-5-[4,5,6,7-tetrahydro-6,6-dimethyl-4-oxo-3-(trifluoromethyl)-1H-indazol-1-yl]phenyl]amino]cyclohexyl ester

Synonym: SNX-5422
MF: C₂₅H₃₀F₃N₅O₄
FW: 521.5
Purity: ≥98%
UV/Vis.: λ_{max}: 250, 359 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

PF-04929113 is supplied as a crystalline solid. A stock solution may be made by dissolving the PF-04929113 in the solvent of choice, which should be purged with an inert gas. PF-04929113 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of PF-04929113 in these solvents is approximately 1 and 0.25 mg/ml, respectively.

PF-04929113 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PF-04929113 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. PF-04929113 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

PF-04929113 is a prodrug for the Hsp90 inhibitor, PF-04928473 (SNX-2112), which binds both Hsp90 α and Hsp90 β with an IC₅₀ value of 30 nM.^{1,2} The prodrug, PF-04929113, is rapidly absorbed and converted into the active inhibitor after oral administration.² The active inhibitor causes degradation of Hsp90 client proteins, including HER2, and reduces phosphorylation of downstream kinases, including Akt and ERK1/2, leading to apoptosis in cancer cells.^{1,3} Oral administration of the prodrug, PF-04929113, reduces tumor growth and prolongs survival in mouse models of multiple myeloma and prostate cancer.^{3,4}

References

- Chandarlapaty, S., Sawai, A., Ye, Q., *et al.* SNX2112, a synthetic heat shock protein 90 inhibitor, has potent antitumor activity against HER kinase-dependent cancers. *Clin. Cancer Res.* **14**(1), 240-248 (2008).
- Jain, L., Gardner, E.R., Venitz, J., *et al.* Determination of PF-04928473 in human plasma using liquid chromatography with tandem mass spectrometry. *J. Chromatogr. B Analyt. Technol. Biomed. Life Sci.* **878**(30), 3187-3192 (2010).
- Okawa, Y., Hideshima, T., Steed, P., *et al.* SNX-2112, a selective Hsp90 inhibitor, potently inhibits tumor cell growth, angiogenesis, and osteoclastogenesis in multiple myeloma and other hematologic tumors by abrogating signaling via Akt and ERK. *Blood.* **113**(4), 846-855 (2009).
- Lamoureux, F., Thomas, C., Yin, M.-J., *et al.* A novel HSP90 inhibitor delays castrate-resistant prostate cancer without altering serum PSA levels and inhibits osteoclastogenesis. *Clin. Cancer Res.* **17**(8), 2301-2313 (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