

PRODUCT INFORMATION



PSI (trifluoroacetate salt)

Item No. 28382

Formal Name: *tert*-butyl (S)-4-((2S,3S)-2-(((benzyloxy) carbonyl)amino)-3-methylpentanamido)-5-(((S)-1-(((S)-4-methyl-1-oxopentan-2-yl)amino)-1-oxopropan-2-yl)amino)-5-oxopentanoate, trifluoroacetate salt

Synonyms: Proteasome Inhibitor I, Z-Ile-Glu(OtBu)-Ala-Leu-aldehyde, Z-Ile-Glu(OtBu)-Ala-Leu-CHO

MF: C₃₂H₅₀N₄O₈ • XCF₃COOH

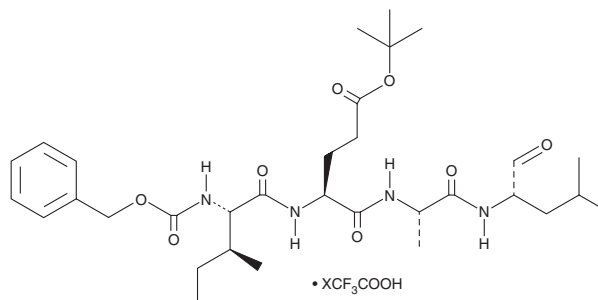
FW: 618.8

Purity: ≥95%

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

PSI (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the PSI (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. PSI (trifluoroacetate salt) is soluble in the organic solvent formic acid at a concentration of approximately 1 mg/ml.

Description

PSI is a synthetic peptide proteasome inhibitor.¹⁻³ It inhibits the acidic and neutral chymotrypsin-like activities of the 20S proteasome (IC₅₀s = 0.25 and 6.5 μM, respectively).³ PSI is cytotoxic to HL-60, KG-1a, RWLeu-4, AR230, and K562 human leukemia cells (IC₅₀s = 4.9, 17.5, 7.8, 7.3, and 10 nM, respectively).⁴ It selectively induces apoptosis in subconfluent bovine aortic endothelial cells (BAECs) and human umbilical vein endothelial cells (HUVECs; EC₅₀s = 24 and 7 nM, respectively) over confluent BAECs and HUVECs (EC₅₀s = 8,150 and 7,830 nM, respectively).⁵ PSI induces accumulation of ubiquitin-protein conjugates in HT4 mouse neuronal cells.⁶

References

- Collins, G.A. and Goldberg, A.L. The logic of the 26S proteasome. *Cell* **169**(5), 792-806 (2017).
- Orlowski, M. and Wilk, S. Catalytic activities of the 20 S proteasome, a multicatalytic proteinase complex. *Arch. Biochem. Biophys.* **383**(1), 1-16 (2000).
- Figueiredo-Pereira, M.E., Chen, W.-E., Yuan, H.-M., et al. A novel chymotrypsin-like component of the multicatalytic proteinase complex optimally active at acidic pH. *Arch. Biochem. Biophys.* **317**(1), 69-78 (1995).
- Soligo, D., Servida, F., Delia, D., et al. The apoptogenic response of human myeloid leukaemia cell lines and of normal and malignant haematopoietic progenitor cells to the proteasome inhibitor PSI. *Br. J. Haematol.* **113**(1), 126-135 (2001).
- Drexler, H.C., Risau, W., and Konerding, M.A. Inhibition of proteasome function induces programmed cell death in proliferating endothelial cells. *FASEB J.* **14**(1), 65-77 (2000).
- Figueiredo-Pereira, M.E., Berg, K.A., and Wilk, S. A new inhibitor of the chymotrypsin-like activity of the multicatalytic proteinase complex (20S proteasome) induces accumulation of ubiquitin-protein conjugates in a neuronal cell. *J. Neurochem.* **63**(4), 1578-1581 (1994).

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, 10/10/2022

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