

PRODUCT INFORMATION



Shepherdin (trifluoroacetate salt)

Item No. 38327

Formal Name: L-lysyl-L-histidyl-L-seryl-L-serylglycyl-L-cysteinyl-L-alanyl-L-phenylalanyl-L-leucine, trifluoroacetate salt

Synonyms: Lys-His-Ser-Ser-Gly-Cys-Ala-Phe-Leu-OH, Survivin (79-87)

Peptide Sequence: KHSSGCAFL-OH

MF: $C_{41}H_{64}N_{12}O_{12}S \cdot XCF_3COOH$

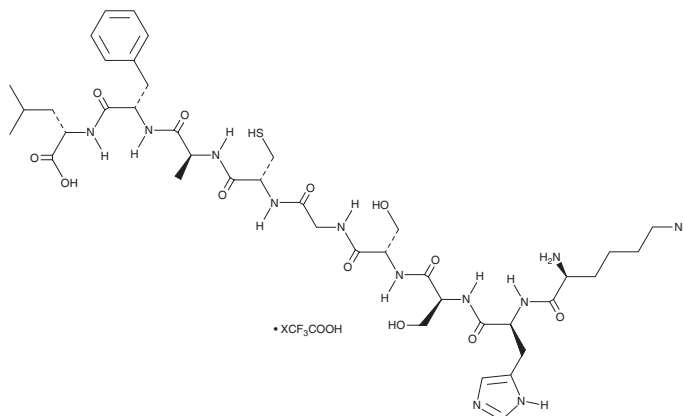
FW: 949.1

Purity: $\geq 90\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

Shepherdin (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the shepherdin (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Shepherdin (trifluoroacetate salt) is soluble in the organic solvent methanol.

Description

Shepherdin is a synthetic peptide that corresponds to amino acids 79-87 of human survivin and an inhibitor of the protein-protein interaction between survivin and heat shock protein 90 (Hsp90).¹ It binds to the N-terminal domain of Hsp90 in a concentration-dependent manner and inhibits the interaction between recombinant human survivin and Hsp90 in isolated human reticulocyte extracts when used at a concentration of 150 μM . Peptide conjugates of shepherdin with the cell-permeable peptides penetratin (Item No. 36878) or HIV Tat sequence have anticancer activity *in vitro* and *in vivo*.

Reference

1. Plescia, J., Salz, W., Xia, F., *et al.* Rational design of shepherdin, a novel anticancer agent. *Cancer Cell* **7**(5), 457-468 (2005).

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, 09/28/2023

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