

PRODUCT INFORMATION



Histatin 8 (trifluoroacetate salt)

Item No. 36891

Formal Name: L-lysyl-L-phenylalanyl-L-histidyl-L- α -glutamyl-L-lysyl-L-histidyl-L-histidyl-L-seryl-L-histidyl-L-arginylglycyl-L-tyrosine, trifluoroacetate salt

Synonym: Hst-8

Peptide Sequence: KFHEKHHSHRGY-OH

MF: C₇₀H₉₉N₂₅O₁₇ • XCF₃COOH

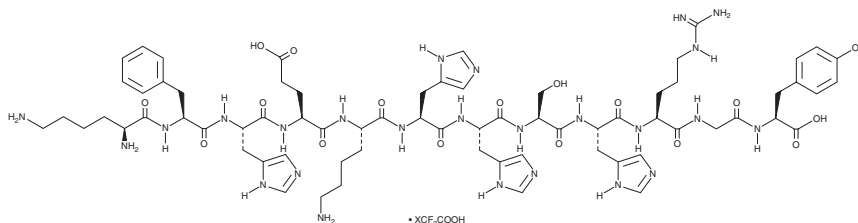
FW: 1,562.7

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

Histatin 8 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the histatin 8 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Histatin 8 (trifluoroacetate salt) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of histatin 8 (trifluoroacetate salt) in these solvents is approximately 10 mg/ml. Histatin 8 (trifluoroacetate salt) is slightly soluble in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of histatin 8 (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of histatin 8 (trifluoroacetate salt) in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Histatin 8 is a salivary antimicrobial peptide and proteolytic fragment of histatin 3 and histatin 5 (Item No. 36721).¹ It is active against *A. baumannii* strains, including strains that are susceptible or resistant to colistin (Item No. 17584), when used at a concentration of 32 μ g/ml. Histatin 8 is also active against *C. albicans* and *C. tropicalis* (MICs = 1 and 5 mg/ml, respectively).²

References

1. Neshani, A., Sedighian, H., Mirhosseini, S.A., *et al.* Antimicrobial peptides as a promising treatment option against *Acinetobacter baumannii* infections. *Microb. Pathog.* **146**, 104238 (2020).
2. Matheson, S.M.G., Cheeptham, N., and Huttunen-Hennelly, H.E.K. Investigating the effects of hydrophobicity and charge on the therapeutic ability of the antimicrobial histatin 8 peptide for potential use in oral applications. *Int. J. Biol.* **5(2)**, 85-91 (2013).

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, 03/25/2026

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