

PRODUCT INFORMATION



Histatin 3 (trifluoroacetate salt)

Item No. 36983

Formal Name: L- α -aspartyl-L-seryl-L-histidyl-L-alanyl-L-lysyl-L-arginyl-L-histidyl-L-histidylglycyl-L-tyrosyl-L-lysyl-L-arginyl-L-lysyl-L-phenylalanyl-L-histidyl-L- α -glutamyl-L-lysyl-L-histidyl-L-histidyl-L-seryl-L-histidyl-L-arginylglycyl-L-tyrosyl-L-arginyl-L-seryl-L-asparaginyl-L-tyrosyl-L-leucyl-L-tyrosyl-L- α -aspartyl-L-asparagine, trifluoroacetate salt

Synonym: Hst-3

Peptide Sequence: DSHAKRHHGKYKRFHEKHSHRGRYSNYLYDN-OH

MF: C₁₇₈H₂₅₈N₆₄O₄₈ • XCF₃COOH

FW: 4,062.4

Purity: \geq 95%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years

H—Asp—Ser—His—Ala—Lys—Arg—His—His—Gly—Tyr—
Lys—Arg—Lys—Phe—His—Glu—Lys—His—His—Ser—
His—Arg—Gly—Tyr—Arg—Ser—Asn—Tyr—Leu—Tyr—
Asp—Asn—OH
• XCF₃COOH

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

Laboratory Procedures

Histatin 3 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the histatin 3 (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Histatin 3 is a salivary antimicrobial peptide.¹ It is active against *C. albicans* (LC₅₀ = 9.2 μ M). Histatin 3 also induces actin polymerization in a cell-free assay when used at a concentration of 12 μ M.²

References

1. Edgerton, M., Koshlukova, S.E., Lo, T.E., *et al.* Candidacidal activity of salivary histatins. *J. Biol. Chem.* **273**(32), 20438-20447 (1998).
2. Blotnick, E., Sol, A., Bachrach, G., *et al.* Interactions of histatin-3 and histatin-5 with actin. *BMC Biochem.* **18**(1), 3 (2017).

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, 10/10/2023

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