

# PRODUCT INFORMATION



## Histatin 1 (trifluoroacetate salt)

Item No. 45085

**Formal Name:** L- $\alpha$ -aspartyl-O-phosphono-L-seryl-L-histidyl-L-glutamyl-L-lysyl-L-arginyl-L-histidyl-L-histidylglycyl-L-tyrosyl-L-arginyl-L-arginyl-L-lysyl-L-phenylalanyl-L-histidyl-L- $\alpha$ -glutamyl-L-lysyl-L-histidyl-L-histidyl-L-seryl-L-histidyl-L-arginyl-L- $\alpha$ -glutamyl-L-phenylalanyl-L-prolyl-L-phenylalanyl-L-tyrosylglycyl-L- $\alpha$ -aspartyl-L-tyrosylglycyl-L-seryl-L-asparaginyl-L-tyrosyl-L-leucyl-L-tyrosyl-L-aspartyl-L-asparagine Hst-1

**Synonym:** Hst-1  
**Peptide Sequence:** H-D-pS-HEKRHHGYRRKFHEKHHSHPFPYGDYGSNYLYDN-OH

**MF:** C<sub>217</sub>H<sub>298</sub>N<sub>69</sub>O<sub>64</sub>P • XCF<sub>3</sub>COOH

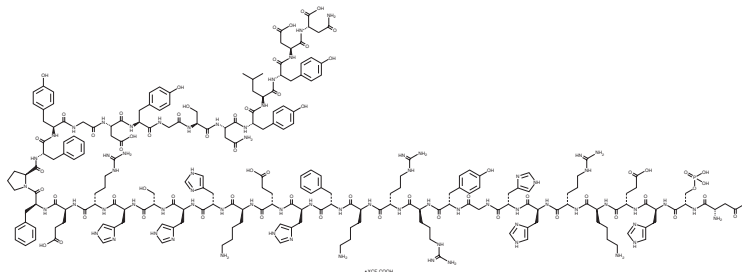
**FW:** 4,928.2

**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 1 (trifluoroacetate salt) is supplied as a solid. Aqueous solutions of histatin 1 (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. Histatin 1 (trifluoroacetate salt) is soluble (≥10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

Histatin 1 is an endogenous peptide secreted by the salivary glands.<sup>1,2</sup> It accelerates wound closure in a scratch assay using TR146 epithelial-like cells and increases adhesion, migration, and the formation of vascular networks in human umbilical vein endothelial cells (HUVECs) and EA.hy926 endothelial cells. Histatin 1 increases angiogenesis in a chick chorioallantoic membrane (CAM) assay.<sup>2</sup> It is active against *C. albicans* (LC<sub>50</sub> = 6  $\mu$ M).<sup>1</sup>

### References

1. Oudhoff, M.J., Bolscher, J.G., Nazmi, K., *et al.* Histatins are the major wound-closure stimulating factors in human saliva as identified in a cell culture assay. *FASEB J.* **22**(11), 3805-3812 (2008).
2. Torres, P., Díaz, J., Arce, M., *et al.* The salivary peptide Histatin 1 promotes endothelial cell adhesion, migration, and angiogenesis. *FASEB J.* **31**(11), 4946-4958 (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, 06/01/2026

#### 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](http://WWW.CAYMANCHEM.COM)