

PRODUCT INFORMATION



Histone H3 (69-89) amide (human, mouse, rat, bovine) (trifluoroacetate salt)

Item No. 28146

Formal Name:	L-arginyl-L-leucyl-L-valyl-L-arginyl-L-glutamyl-L-isoleucyl-L-alanyl-L-glutamyl-L-aspartyl-L-phenylalanyl-L-lysyl-L-threonyl-L-aspartyl-L-leucyl-L-arginyl-L-phenylalanyl-L-glutamyl-L-seryl-L-seryl-L-alanyl-L-valinamide, trifluoroacetate salt	H — Arg — Leu — Val — Arg — Glu — Ile — Ala — Gln — Asp — Phe — Lys — Thr — Asp — Leu — Arg — Phe — Gln — Ser — Ser — Ala — Val — NH ₂
Synonym:	RLVREIAQDFKTDLRFSQSSAV-NH ₂	• XCF ₃ COOH
MF:	C ₁₀₉ H ₁₈₀ N ₃₄ O ₃₂ • XCF ₃ COOH	
FW:	2,478.8	
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

Histone H3 (69-89) amide (human, mouse, rat, bovine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the histone H3 (69-89) amide (human, mouse, rat, bovine) (trifluoroacetate salt) in water. The solubility of histone H3 (69-89) amide (human, mouse, rat, bovine) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Histone H3 (69-89) amide is a peptide fragment of histone H3 that corresponds to amino acid residues 70-90 of the human histone H3.1 and H3.2 sequences. Histone H3 (69-89) amide contains a lysine residue at position 79 that is subject to methylation and acetylation and has a role in regulation of gene expression.^{1,2}

References

1. Garcia, B.A., Hake, S.B., Diaz, R.L., *et al.* Organismal differences in post-translational modifications in histones H3 and H4. *J. Biol. Chem.* **282**(10), 7641-7655 (2007).
2. Barski, A., Cuddapah, S., Cui, K., *et al.* High-resolution profiling of histone methylations in the human genome. *Cell* **129**(4), 823-837 (2007).

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, 12/05/2022

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