

PRODUCT INFORMATION

Histone H3 (21-44) (human, mouse, rat, bovine) (trifluoroacetate salt)
Item No. 27760

Formal Name: L-alanyl-L-threonyl-L-lysyl-L-alanyl-L-alanyl-L-arginyl-L-lysyl-L-seryl-L-alanyl-L-prolyl-L-alanyl-L-threonylglycylglycyl-L-valyl-L-lysyl-L-lysyl-L-prolyl-L-histidyl-L-arginyl-L-tyrosyl-L-arginyl-L-prolyl-glycine, trifluoroacetate salt
Synonym: ATKAARKSAPATGGVKKPHRYRPG
MF: $C_{109}H_{185}N_{39}O_{29} \cdot XCF_3COOH$
FW: 2,505.9
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years

H—Ala—Thr—Lys—Ala—Ala—Arg—Lys—Ser—Ala—Pro—
Ala—Thr—Gly—Gly—Val—Lys—Lys—Pro—His—Arg—
Tyr—Arg—Pro—Gly—OH
• XCF₃COOH

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

Laboratory Procedures

Histone H3 (21-44) (human, mouse, rat, bovine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the histone H3 (21-44) (human, mouse, rat, bovine) (trifluoroacetate salt) in water. The solubility of histone H3 (21-44) (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 (21-44) is a peptide fragment of histone H3 that corresponds to amino acid residues 22-45 of the human histone H3.1 and H3.2 sequences. It has been used as a peptide substrate to quantify and compare the histone H3 lysine 27 methyltransferase activity and enzyme kinetics of recombinant polycomb repressive complex 2 (PRC2) complexes containing wild-type or mutant versions of the catalytic subunit EZH2.¹

Reference

1. Sneeringer, C.J., Scott, M.P., Kuntz, K.W., *et al.* Coordinated activities of wild-type plus mutant EZH2 drive tumor-associated hypertrimethylation of lysine 27 on histone H3 (H3K27) in human B-cell lymphomas. *Proc. Nat. Acad. Sci. USA* **107**(49), 20980-20985 (2010).

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, 12/02/2022

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