

PRODUCT INFORMATION

Histone H3 (5-23) (human, mouse, rat, porcine, bovine) (trifluoroacetate salt)

Item No. 28048

Synonym: QTARKSTGGKAPRKQLATK
MF: $C_{85}H_{155}N_{31}O_{26} \cdot XCF_3COOH$ H—Gln—Thr—Ala—Arg—Lys—Ser—Thr—Gly—Gly—Lys—
FW: 2,027.4 Ala—Pro—Arg—Lys—Gln—Leu—Ala—Thr—Lys—OH
Purity: $\geq 95\%$ • XCF_3COOH
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years

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

Laboratory Procedures

Histone H3 (5-23) (human, mouse, rat, porcine, bovine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the histone H3 (5-23) (human, mouse, rat, porcine, bovine) (trifluoroacetate salt) in water. The solubility of histone H3 (5-23) (human, mouse, rat, porcine, 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 (5-23) is a peptide fragment of histone H3 that corresponds to amino acid residues 6-24 of the human histone H3.3 sequence. Histone H3 contains a lysine residue at position 14 that is subject to acetylation and has a role in transcriptional activation and regulation of gene expression.^{1,2}

References

1. Holbert, M.A., Sikorski, T., Carten, J., *et al.* The human monocytic leukemia zinc finger histone acetyltransferase domain contains DNA-binding activity implicated in chromatin targeting. *J. Biol. Chem.* **282(50)**, 36603-36613 (2007).
2. Karmodiya, K., Krebs, A.R., Oulad-Abdelghani, M., *et al.* H3K9 and H3K14 acetylation co-occur at many gene regulatory elements, while H3K14ac marks a subset of inactive inducible promoters in mouse embryonic stem cells. *BMC Genomics* **13:424** (2012).

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