

PRODUCT INFORMATION



Histone H3K36Me3 (21-44)-GK-biotin (trifluoroacetate salt)

Item No. 27532

Formal Name: N²-((S)-2-((S)-2-(2-(2-((2S,3R)-2-((S)-2-((S)-1-(L-alanyl-L-threonyl-L-lysyl-L-alanyl-L-alanyl-L-arginyl-L-lysyl-L-seryl-L-alanyl)pyrrolidine-2-carboxamido)propanamido)-3-hydroxybutanamido)acetamido)acetamido)-3-methylbutanamido)-6-(trimethylammonio)hexanoyl)-L-lysyl-L-prolyl-L-histidyl-L-arginyl-L-tyrosyl-L-arginyl-L-prolylglycylglycyl-N⁶-(5-((3aS,4S,6aR)-2-oxohexahydro-1H-thieno[3,4-d]imidazol-4-yl)pentanoyl)-L-lysinate, trifluoroacetate salt

Synonyms: ATKAARKSAPATGGV-K(Me3)-KPHRYRPG-GK(Biotin), Histone H3 (21-44) (Lys³⁶me3), [Lys(Me3)₃₆]-Histone H3 (21-44)-GK(Biotin)

MF: C₁₃₀H₂₂₀N₄₄O₃₃S • XCF₃COOH

FW: 2,959.5

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.

H—Ala—Thr—Lys—Ala—Ala—Arg—Lys—Ser—Ala—Pro—

Ala—Thr—Gly—Gly—Val—Lys(Me3)—Lys—Pro—His—Arg—

Tyr—Arg—Pro—Gly—Gly—Lys(Biotin)—OH

• XCF₃COOH

Laboratory Procedures

Histone H3K36Me3 (21-44)-GK-biotin (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the histone H3K36Me3 (21-44)-GK-biotin (trifluoroacetate salt) in water. The solubility of histone H3K36Me3 (21-44)-GK-biotin (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Histone H3K36Me3 (21-44) is a peptide fragment of histone H3 that corresponds to amino acid residues 22-45 of the human histone H3.1 and 3.2 sequences. It is trimethylated at lysine 36 and biotinylated via a C-terminal GK linker. Trimethylation of H3K36 is increased in the G₁ and early S phases of the cell cycle where it binds to and recruits the mismatch recognition protein MutSa in preparation for DNA replication and mismatch repair.¹ H3K36 trimethylation is found in greater amounts on exons compared with introns in *C. elegans*, mouse, and human genome-wide maps of histone H3 tail methylations.²

References

1. Li, F., Mao, G., Tong, D., *et al.* The histone mark H3K36me3 regulates human DNA mismatch repair through its interaction with MutSa. *Cell* **153**(3), 590-600 (2013).
2. Kolasinska-Zwierz, P., Down, T., Latorre, I., *et al.* Differential chromatin marking of introns and expressed exons by H3K36me3. *Nat. Genet.* **41**(3), 376-381 (2009).

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