

PRODUCT INFORMATION



Histone H3K36Me2 (21-44)-GK-biotin (trifluoroacetate salt)

Item No. 27531

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-N ⁶ ,N ⁶ -dimethyl-L-lysyl-L-lysyl-L-prolyl-L-histidyl-L-arginyl-L-tyrosyl-L-arginyl-L-prolylglycylglycyl-L-lysine-biotin, trifluoroacetate salt	
Synonyms:	ATKAARKSAPATGGV-K(Me2)-KPHRYRPG-GK(Biotin), [Lys(Me2)36]-Histone H3 (21-44)-GK(Biotin), Histone H3 (21-44) (Lys ³⁶ me2)-biotin	H—Ala—Thr—Lys—Ala—Ala—Arg—Lys—Ser—Ala—Pro—Ala—Thr—Gly—Gly—Val—Lys(Me2)—Lys—Pro—His—Arg—Tyr—Arg—Pro—Gly—Gly—Lys(Biotin)—OH
MF:	C ₁₂₉ H ₂₁₈ N ₄₄ O ₃₃ S • XCF ₃ COOH	• XCF ₃ COOH
FW:	2,945.5	
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.

Laboratory Procedures

Histone H3K36Me2 (21-44)-GK-biotin (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the histone H3K36Me2 (21-44)-GK-biotin (trifluoroacetate salt) in water. The solubility of histone H3K36Me2 (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 H3K36Me2 (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 is dimethylated at lysine 36 and biotinylated via a C-terminal GK linker. Dimethylation of H3K36 is enriched at coding regions of genes and correlates with the initiation, but not maintenance, of active transcription.^{1,2}

References

- Li, B., Jackson, J., Simon, M.D., *et al.* Histone H3 lysine 36 dimethylation (H3K36me2) is sufficient to recruit the Rpd3s histone deacetylase complex and to repress spurious transcription. *J. Biol. Chem.* **284**(12), 7970-7976 (2009).
- Rao, B., Shibata, Y., Strahl, B.D., *et al.* Dimethylation of histone H3 at lysine 36 demarcates regulatory and nonregulatory chromatin genome-wide. *Mol. Cell. Biol.* **25**(21), 9447-9459 (2005).

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