

PRODUCT INFORMATION



Histone H3K36Me3 (26-46)-K-biotin (trifluoroacetate salt)

Item No. 28144

Formal Name: L-arginyl-L-lysyl-L-alanyl-L-alanyl-L-prolyl-L-alanyl-L-threonylglycylglycyl-L-valyl-L-6-(trimethylammonio)-L-norleucyl-L-lysyl-L-prolyl-L-histidyl-L-arginyl-L-tyrosyl-L-arginyl-L-prolylglycyl-L-threonyl-L-valyl-L-lysine-biotin, trifluoroacetate salt

Synonyms: Histone H3 (26-46) (Lys³⁶me₃), [Lys(Me₃)₃₆]-Histone H3 (26-46)-K(Biotin), RKAAPATGGV-K(Me₃)-KPHRYRPGTV-K(biotin)



MF: C₁₁₈H₁₉₉N₃₉O₂₈S • XCF₃COOH

FW: 2,644.2

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 H3K36Me3 (26-46)-K-biotin (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the histone H3K36Me3 (26-46)-K-biotin (trifluoroacetate salt) in water. The solubility of histone H3K36Me3 (26-46)-K-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 (26-46)-K-biotin is a peptide fragment of histone H3 that contains a serine-to-alanine substitution at position 29 and corresponds to amino acid residues 27-47 of the human histone H3.1 and H3.2 sequences. It is trimethylated at lysine 36 and biotinylated via a C-terminal lysine linker. Histone H3K36Me3 (26-46)-K-biotin corresponds directly to amino acid residues 27-47 of the yeast H3.1 sequence in *S. octosporus* and *S. pombe* and of the plant pathogenic fungus histone sequence in *P. striiformis*.

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, 12/08/2022

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