

PRODUCT INFORMATION



Histone H3K27Me1 (23-34) (trifluoroacetate salt)

Item No. 28141

| | | |
|---------------------|--|---|
| Formal Name: | L-lysyl-L-alanyl-L-alanyl-L-arginyl-N ⁶ -methyl-L-lysyl-L-seryl-L-alanyl-L-prolyl-L-alanyl-L-threonylglycylglycine, trifluoroacetate salt | |
| Synonyms: | Histone H3 (23-34) (Lys ²⁷ me1), KAAR-K(Me1)-SAPATGG, [Lys(Me1)27]-Histone H3 (23-34) | H-Lys-Ala-Ala-Arg-Lys(Me1)-Ser-Ala-Pro-Ala-Thr-Gly-Gly-OH |
| MF: | C ₄₇ H ₈₅ N ₁₇ O ₁₅ • XCF ₃ COOH | • XCF ₃ COOH |
| FW: | 1,128.3 | |
| 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 H3K27Me1 (23-34) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the histone H3K27Me1 (23-34) (trifluoroacetate salt) in water. The solubility of histone H3K27Me1 (23-34) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Histone H3K27Me1 (23-34) is a peptide fragment of histone H3 that corresponds to amino acid residues 24-35 of the human histone H3.1 and H3.2 sequences. Monomethylation of histone H3 at lysine 27 is associated with actively transcribed genes and positively correlates with H3K36 trimethylation.¹ Levels of H3K27Me1 are increased in tumor tissue isolated from patients with metastatic hormone-naïve and castration-resistant prostate cancer.² Histone H3K27Me1 (23-34) has been used in epitope mapping of the lupus-derived monoclonal antibody BT164.³

References

1. Ferrari, K.J., Scelfo, A., Jammula, S., *et al.* Polycomb-dependent H3K27me1 and H3K27me2 regulate active transcription and enhancer fidelity. *Mol. Cell* **53**(1), 49-62 (2014).
2. Ellinger, J., Kahl, P., von der Gathen, J., *et al.* Global histone H3K27 methylation levels are different in localized and metastatic prostate cancer. *Cancer Invest.* **30**(2), 92-97 (2012).
3. van Bavel, C.C., Dieker, J., Kroeze, Y., *et al.* Apoptosis-induced histone H3 methylation is targeted by autoantibodies in systemic lupus erythematosus. *Ann. Rheum. Dis.* **70**(1), 201-207 (2011).

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, 10/11/2022

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