

# PRODUCT INFORMATION



## Histone H3K27Me3 (23-34) (human, mouse, rat, bovine) (trifluoroacetate salt) Item No. 28143

<b>Formal Name:</b>	L-lysyl-L-alanyl-L-alanyl-L-arginyl-L-6-(trimethylammonio)-L-norleucyl-L-seryl-L-alanyl-L-prolyl-L-alanyl-L-threonylglycylglycine, trifluoroacetate salt	
<b>Synonyms:</b>	[Lys(Me3)27]-Histone H3 (23-34), Histone H3 (23-34) (Lys <sup>27</sup> me3), KAAR-K(Me3)-SAPATGG	H—Lys—Ala—Ala—Arg—Lys(Me3)—Ser—Ala—Pro—Ala—Thr—Gly—Gly—OH • XCF <sub>3</sub> COOH
<b>MF:</b>	C <sub>49</sub> H <sub>89</sub> N <sub>17</sub> O <sub>15</sub> • XCF <sub>3</sub> COOH	
<b>FW:</b>	1,156.3	
<b>Purity:</b>	≥95%	
<b>Supplied as:</b>	A solid	
<b>Storage:</b>	-20°C	
<b>Stability:</b>	≥4 years	

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

### Laboratory Procedures

Histone H3K27Me3 (23-34) (human, mouse, rat, bovine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the histone H3K27Me3 (23-34) (human, mouse, rat, bovine) (trifluoroacetate salt) in water. The solubility of histone H3K27Me3 (23-34) (human, mouse, rat, 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 H3K27Me3 (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. Trimethylation of histone H3 at lysine 27 is associated with gene silencing.<sup>1</sup> It is involved in tumor progression through its regulation by enhancer of zeste homolog 2 (EZH2) and transcriptional repression of tumor suppressor genes.<sup>2,3</sup> Levels of H3K27Me3 are reduced in 293 T-REX cells containing EED<sup>R236T</sup> and SUZ12<sup>G610V</sup> mutations and in lymphoblastoid cells isolated from patients with Weaver syndrome, a rare overgrowth disorder characterized by EZH2, EED, or SUZ12 mutations, cancer susceptibility, and various distinctive physical features.<sup>4</sup> Histone H3K27Me3 (23-34) has been used in AlphaLISA<sup>®</sup> assays to verify antibody specificity.<sup>5</sup>

### References

1. Becker, J.S., Nicetto, D., and Zaret, K.S. H3K9me3-dependent heterochromatin: Barrier to cell fate changes. *Trends Genet.* **32(1)**, 29-41 (2016).
2. Wu, Z., Lee, S.T., Qiao, Y., et al. Polycomb protein EZH2 regulates cancer cell fate decision in response to DNA damage. *Cell Death Differ.* **18(11)**, 1771-1779 (2011).
3. Gan, L., Yang, Y., Li, Q., et al. Epigenetic regulation of cancer progression by EZH2: From biological insights to therapeutic potential. *Biomark. Res.* **6**, 10 (2018).
4. Imagawa, E., Higashimoto, K., Sakai, Y., et al. Mutations in genes encoding polycomb repressive complex 2 subunits cause Weaver syndrome. *Hum. Mutat.* **38(6)**, 637-648 (2017).
5. Qian, J., Lu, L., Wu, J., et al. Development of multiple cell-based assays for the detection of histone H3 Lys27 trimethylation (H3K27me3). *Assay Drug Dev. Technol.* **11(7)**, 449-456 (2013).

#### 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, 12/05/2022

#### 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