

# PRODUCT INFORMATION



## Histone H3K27Ac (21-44)-GK-biotin amide (trifluoroacetate salt)

Item No. 27764

**Formal Name:** L-alanyl-L-threonyl-L-lysyl-L-alanyl-L-alanyl-L-arginyl-N<sup>6</sup>-acetyl-L-lysyl-L-seryl-L-alanyl-L-prolyl-L-seryl-L-threonylglycylglycyl-L-valyl-L-lysyl-L-lysyl-L-prolyl-L-histidyl-L-arginyl-L-tyrosyl-L-arginyl-L-prolylglycylglycyl-L-lysine-biotin amide, trifluoroacetate salt

**Synonym:** ATKAAR-K(Ac)-SAPSTGGVKKPHRYRPGG-K(Biotin)-NH<sub>2</sub>

**MF:** C<sub>129</sub>H<sub>217</sub>N<sub>45</sub>O<sub>34</sub>S • XCF<sub>3</sub>COOH

**FW:** 2,974.5

**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 H3K27Ac (21-44)-GK-biotin amide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the histone H3K27Ac (21-44)-GK-biotin amide (trifluoroacetate salt) in water. The solubility of histone H3K27Ac (21-44)-GK-biotin amide (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Histone H3K27Ac (21-44)-GK-biotin amide is a peptide fragment of histone H3 that corresponds to amino acid residues 22-45 of the human histone H3.3 sequence. It is acetylated at lysine 27 and biotinylated via a C-terminal GK linker. Acetylation of histone H3 at lysine 27 is associated with active gene transcription.<sup>1</sup> H3K27 acetylation and H3K4 monomethylation are enriched near active enhancers and H3K27 acetylation and H3K4 trimethylation together mark transcription start sites of actively transcribed genes.

### Reference

1. Kimura, H. Histone modifications for human epigenome analysis. *J. Hum. Genet.* **58(7)**, 439-445 (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, 12/02/2022

#### 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