

PRODUCT INFORMATION



Histone H3 (1-35) (human, mouse, rat, bovine) (trifluoroacetate salt) Item No. 27524

Formal Name: L-alanyl-L-arginyl-L-threonyl-L-lysyl-L-glutaminy-
L-threonyl-L-alanyl-L-arginyl-L-lysyl-L-seryl-L-
threonylglycylglycyl-L-lysyl-L-alanyl-L-prolyl-L-arginyl-
L-lysyl-L-glutaminy-L-leucyl-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-valine,
trifluoroacetate salt



Synonym: ARTKQTARKSTGGKAPRKQLATKAARKSAPATGGV

MF: $\text{C}_{149}\text{H}_{269}\text{N}_{55}\text{O}_{45} \bullet \text{XCF}_3\text{COOH}$

FW: 3,551.1

Purity: $\geq 95\%$

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 H3 (1-35) (human, mouse, rat, bovine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the histone H3 (1-35) (human, mouse, rat, bovine) (trifluoroacetate salt) in water. The solubility of histone H3 (1-35) (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 H3 (1-35) is an N-terminal peptide fragment of histone H3 that corresponds to amino acid residues 2-36 of the human histone H3.1 and 3.2 sequences. Histone H3 (1-35) contains lysine residues at positions 4, 9, and 27 that are subject to methylation and acetylation and a serine residue at position 10 that is subject to phosphorylation, all of which have a role in the regulation of gene expression.¹

Reference

1. Bhaumik, S.R., Smith, E., and Shilatifard, A. Covalent modifications of histones during development and disease pathogenesis. *Nat. Struct. Mol. Biol.* **14(11)**, 1008-1016 (2007).

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, 12/01/2022

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