

PRODUCT INFORMATION

Histone H2A (1-20) (human, mouse, rat, bovine) (trifluoroacetate salt)
Item No. 27440

Formal Name: L-serylglycyl-L-arginylglycyl-L-lysyl-L-glutaminyglycylglycyl-L-lysyl-L-alanyl-L-arginyl-L-alanyl-L-lysyl-L-alanyl-L-lysyl-L-threonyl-L-arginyl-L-seryl-L-seryl-L-arginine, trifluoroacetate salt

Synonyms: H-Ser-Gly-Arg-Gly-Lys-Gln-Gly-Gly-Lys-Ala-Arg-Ala-Lys-Ala-Lys-Thr-Arg-Ser-Ser-Arg-OH, SGRGKQGKARAKAKTRSSR

H-Ser-Gly-Arg-Gly-Lys-Gln-Gly-Gly-Lys-Ala-Arg-Ala-Lys-Ala-Lys-Thr-Arg-Ser-Ser-Arg-OH
• XCF₃COOH

MF: C₈₃H₁₅₅N₃₇O₂₆ • XCF₃COOH
FW: 2,087.4
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 H2A (1-20) (human, mouse, rat, bovine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the histone H2A (1-20) (human, mouse, rat, bovine) (trifluoroacetate salt) in water. The solubility of histone H2A (1-20) (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 H2A (1-20) is an N-terminal peptide fragment of histone H2A that corresponds to amino acid residues 2-21 of the human histone H2A sequence. Methylation of the arginine at position 3 of histone H2A by protein arginine methyltransferase 1 (PRMT1), PRMT5, or PRMT6 is associated with both transcriptional activation and repression.¹ It is associated with DNA damage repair when methylated by PRMT7. The arginine residue at position 3 of histone H2A is also subject to citrullination by protein arginine deiminase 4 (PAD4).

Reference

1. Fuhrmann, J. and Thompson, P.R. Protein arginine methylation and citrullination in epigenetic regulation. *ACS Chem. Biol.* **11**(3), 654-668 (2016).

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, 11/30/2022

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