

PRODUCT INFORMATION



Histone H3 (1-15) (Phospho-Ser¹⁰) (human, mouse, rat, porcine, bovine) (trifluoroacetate salt)

Item No. 27494

Formal Name:	N-L-alanyl-L-arginyl-L-threonyl-L-lysyl-L-glutamyl-L-threonyl-L-alanyl-L-arginyl-L-lysyl-O-phosphono-L-seryl-L-threonylglycylglycyl-L-lysyl-L-alanine, trifluoroacetate salt	H—Ala—Arg—Thr—Lys—Gln—Thr—Ala—Arg—Lys—pSer—
Synonyms:	ARTKQTARK-pS-TGGKA, [pSer10]-Histone H3 (1-15), 1-15 H3pS10, H3S10ph (1-15), H3 ₁₋₁₅ S10ph	Thr—Gly—Gly—Lys—Ala—OH • XCF ₃ COOH
MF:	C ₆₃ H ₁₁₈ N ₂₅ O ₂₄ P • XCF ₃ COOH	
FW:	1,640.8	
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 H3 (1-15) (phospho-Ser¹⁰) (human, mouse, rat, porcine, bovine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the histone H3 (1-15) (phospho-Ser¹⁰) (human, mouse, rat, porcine, bovine) (trifluoroacetate salt) in water. The solubility of histone H3 (1-15) (phospho-Ser¹⁰) (human, mouse, rat, porcine, 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 (1-15) (phospho-Ser¹⁰) is a peptide fragment that corresponds to amino acid residues 2-16 of the human histone H3 sequence. Phosphorylation of histone H3 at serine 10 is correlated with chromatin condensation during mitosis and with transcriptional activation of genes during interphase.^{1,2} Histone H3 (1-15) (phospho-Ser¹⁰) does not inhibit binding of the ADD domain of the chromatin-remodeling protein ATRX to histone H3 due to its positioning away from the core peptide binding sequence but does decrease binding of the HP1α chromodomain (CD) when histone H3 is also methylated at lysine 9.³

References

1. Sawicka, A. and Seiser, C. Histone H3 phosphorylation - A versatile chromatin modification for different occasions. *Biochimie* **94(11)**, 2193-2201 (2012).
2. Prigent, C. and Dimitrov, S. Phosphorylation of serine 10 in histone H3, what for? *J. Cell. Sci.* **116 (Pt 18)**, 3677-3685 (2003).
3. Noh, K.M., Maze, I., Zhao, D., *et al.* ATRX tolerates activity-dependent histone H3 methyl/phos switching to maintain repetitive element silencing in neurons. *Proc. Natl. Acad. Sci. USA* **112(22)**, 6820-6827 (2015).

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