

PRODUCT INFORMATION

Histone H3K79Ac (73-83) (human, mouse, rat, porcine, bovine) (trifluoroacetate salt)

Item No. 27489

Formal Name:	L- α -glutamyl-L-isoleucyl-L-alanyl-L-glutaminyl-L- α -aspartyl-L-phenylalanyl-N ⁶ -acetyl-L-lysyl-L-threonyl-L- α -aspartyl-L-leucyl-L-arginine, trifluoroacetate salt
Synonyms:	EIAQDF-K(Ac)-TDLR, H-Glu-Ile-Ala-Gln-Asp-Phe-Lys(Ac)-Thr-Asp-Leu-Arg-OH, Histone H3 (73-83) (Lys ⁷⁹ ac), H3K79Ac, [Lys(Ac)79]-Histone H3 (73-83)
MF:	C ₆₀ H ₉₆ N ₁₆ O ₂₁ • XCF ₃ COOH
FW:	1,377.5
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 H3K79Ac (73-83) (human, mouse, rat, porcine, bovine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the histone H3K79Ac (73-83) (human, mouse, rat, porcine, bovine) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Histone H3K79Ac (73-83) (human, mouse, rat, porcine, bovine) (trifluoroacetate salt) is soluble in the organic solvent formic acid at a concentration of approximately 1 mg/ml.

Description

Histone H3K79Ac (73-83) is a peptide fragment of histone H3 that corresponds to amino acid residues 74-84 of the human histone H3 sequence. Acetylation of histone H3 at lysine 79 has been detected in humans and yeast and is associated with inactive chromatin.¹⁻³

References

1. Garcia, B.A., Hake, S.B., Diaz, R.L., *et al.* Organismal differences in post-translational modifications in histones H3 and H4. *J. Biol. Chem.* **282**(10), 7641-7655 (2007).
2. Bheda, P., Swatkoski, S., Fiedler, K.L., *et al.* Biotinylation of lysine method identifies acetylated histone H3 lysine 79 in *Saccharomyces cerevisiae* as a substrate for Sir2. *Proc. Natl. Acad. Sci. U.S.A.* **109**(16), E916-E925 (2012).
3. Gatta, R. and Mantovani, R. Single nucleosome ChIPs identify an extensive switch of acetyl marks on cell cycle promoters. *Cell Cycle* **9**(11), 2149-2159 (2010).

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