

PRODUCT INFORMATION



Histone H4 Peptide Substrate (15-24)

Item No. 10380

Supplied As: A lyophilized peptide
Synonym: Histone H4K20 Peptide, SET8 Methyltransferase Acceptor Peptide
Peptide Sequence: Human histone H4 amino acids 15-24: AKRHRKVL RD-NH₂
FW: 1,280
Storage: -20°C
Stability: ≥1 year
Solubility: 20 mg/ml with water or aqueous buffers

Description

Histone H4 undergoes many modifications including acetylation, methylation, and phosphorylation, each important for regulation of gene transcription.^{1,2} This histone H4 peptide contains a lysine at position 20 which is a substrate or acceptor peptide for the lysine methyltransferases KMT5A (Set8) and KMT5B (SUV4-20H1).^{3,4}

References

1. Kouzarides, T. Chromatin modifications and their function. *Cell* **128**, 693-705 (2007).
2. Pennisi, E. Opening the way to gene activity. *Science* **275**, 155-157 (1997).
3. Yin, Y., Liu, C., Tsai, S.N., *et al.* SET8 recognizes the sequence RHRK20VLRDN within the N terminus of histone H4 and mono-methylates lysine 20. *J. Biol. Chem.* **280**(34), 30025-30031 (2005).
4. Couture, J.-F., Collazo, E., Brunzelle, J.S., *et al.* Structural and functional analysis of SET8, a histone H4 Lys-20 methyltransferase. *Genes Dev.* **19**, 1455-1465 (2005).

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, 10/23/2019

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