

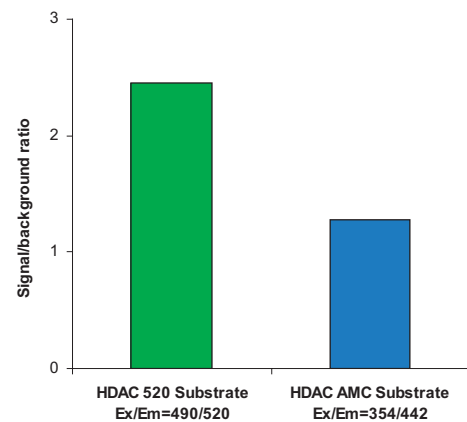
HDAC assays & Histone peptides

SensoLyte™ 520 HDAC Assay Kit (cat# 72084)

- Sensitive fluorimetric assay
- Convenient 2-step homogeneous cell based assay
- Industry's longest HDAC assay kit, Ex/Em=490 nm/520 nm
- Minimum autofluorescence from cell components and test compounds
- Cell permeant substrate

SensoLyte™ 440 HDAC Assay Kit (cat# 72083)

- Fluorimetric assay
- Convenient 2-step homogeneous cell based assay
- Ac-GAK(Ac)-AMC substrate, Ex/Em=354 nm/442 nm
- Cell permeant substrate



Related Products:

- Histone Deacetylase HDAC Substrate Ac-GAK(Ac)-AMC, Ex/Em=354 nm/442 nm, cat# 61855 (1 mg)
- HeLa Nuclear Extract, cat# 72085 (100 ul)

Histone Peptides

The core histones (H2A, H2B, H3 and H4) act as spools around which DNA winds. Covalent modification on the amino terminal of the core histones through de/acetylation, de/methylation and de/phosphorylation affect chromatin structure and exquisitely regulate gene expression [Jenuwein, T. and CD. Allis Science 293, 1074 (2001); Sterner, DE. et al. Microbiol. Mol. Biol. Rev. 64, 435 (2000); Lachner, M. et al. J. Cell Sci. 116, 2117 (2003)]. Histone hyperacetylation is correlated with increased transcription, whereas hypoacetylation correlates with transcriptional repression [Sterner, DE. et al. Microbiol. Mol. Biol. Rev. 64, 435 (2000)].

- Wide selection of histone peptides manufactured by AnaSpec
- H1 H2A, H3 and H4 histone peptide fragments
- Methylated and acetylated peptides



ISO 9001:2000 Company

Histone Peptides

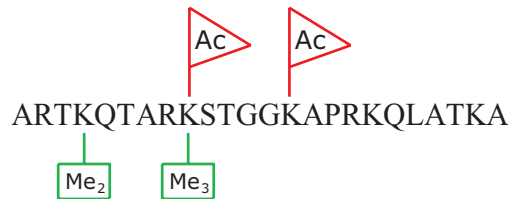
H1 Peptides

Product	Size	Cat #
Histone H1' (22-42), human, mouse STDHPKYSDMIVAAIQAEKNR	1 mg	63674
Histone H1 PKTPKKAKKL	1 mg	60026-1

H2A Peptides

Product	Size	Cat #
Histone H2A (1-20) STDHPKYSDMIVAAIQAEKNR	1 mg	63676
[Lys(Biotin) ⁹]-Histone H2A (1-12)-Cys SGRGKQGG-K(Biotin)-ARAC	1 mg	63675

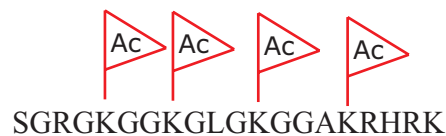
H3 Peptides



Product	Size	Cat #
Histone H3 (1-20), N-Terminal ARTKQTARKSTGGKAPRKQL	1 mg	62753
Histone H3 (1-21), N-Terminal ARTKQTARKSTGGKAPRKQLA	1 mg	61701
Histone H3 (1-21), Biotinylated ARTKQTARKSTGGKAPRKQLA-GG- K(BIOTIN)-NH ₂	1 mg	61702
Histone H3 (1-25), amide ARTKQTARKSTGGKAPRKQLATKAA-NH ₂	1 mg	61703
Histone H3 (15-39) APRKQLATKAARKSAPATGGVKKPH	1 mg	61705
Histone H3 (116-136), C116-136 KRVTIMPKDIQLARRIRGERA	1 mg	61704

Product	Size	Cat #
[Lys(Ac) ¹⁴]-Histone H3 (1-19), H3K14(Ac) ARTKQTARKSTGG-K(Ac)-APRKQL	1 mg	63682
[Lys(Ac) ⁹]-Histone H3 (1-20), H3K9(Ac) ARTKQTAR-K(Ac)-STGGKAPRKQL	1 mg	63680
[Lys(Ac) ¹⁴]-Histone H3 (9-19), H3K14(Ac) KSTGG-K(Ac)-APRKQ	1 mg	63681
[Lys(Me ₂) ⁴]-Histone H3 (1-21), H3K4(Me ₂) ART-K(Me ₂)-QTARKSTGGKAPRKQLA	1 mg	63677
[Lys(Me ₃) ⁹]-Histone H3 (1-24), H3K9(Me ₃) ARTKQTAR-K(Me ₃)-STGGKAPRKQLATKA	0.5 mg	63679

H4 Peptides



Product	Size	Cat #
Histone H4 (1-7), N-Terminal SGRGKGG	1 mg	62754
Histone H4 (1-20), PRMT7 Substrate, M1 SGRGKGGKGLGKGGAKRHRK-NH ₂	1 mg	62498
Histone H4 (1-21), p300/CBP Substrate SGRGKGGKGLGKGGAKRHRKV	1 mg	62499
[Lys(Ac) ⁵]-Histone H4 (1-20), H4K5(Ac) SGRG-K(Ac)-GGKGLGKGGAKRHRK	1 mg	63683
[Lys(Ac) ⁸]-Histone H4 (1-20), H4K8(Ac) SGRGKGG-K(Ac)-GLGKGGAKRHRK	1 mg	63684

Product	Size	Cat #
[Lys(Ac) ¹²]-Histone H4 (1-20), H4K12(Ac) SGRGKGGKGLG-K(Ac)-GGAKRHRK	1 mg	63685
[[Lys(Ac) ¹⁶]-Histone H4 (1-20), H4K16(Ac) SGRGKGGKGLGKGGK-Ac(Ac)-RHRK	1 mg	63687
[Lys(Ac) ¹⁶]-Histone H4 (14-21), Mca- labeled, H4K16(Ac) (14-21), pepH4 - AcK16 (Ex/Em = 325 nm/393 nm) Mca-GA-K(Ac)-RHRKV-NH ₂	1 mg	63686