

Product Data Sheet

 $\label{eq:Lys} $[Lys(Me_2)^{27},\ Lys(Me_1)^{36}]$-Histone H3 (21-44)-GK(Biotin), $H3K27(Me2)K36(Me1),$ biotin-labeled$ **Product Name:**

Catalog Number: AS-65449-1 (1 mg) Lot Number: See label on vial

Sequence: H-Ala-Thr-Lys-Ala-Arg-Lys(Me2)-Ser-Ala-Pro-Ala-Thr-Gly-Gly-Val-

Lys(Me1)-Lys-Pro-His-Arg-Tyr-Arg-Pro-Gly-Gly-Lys(Biotin)-OH

ATKAAR-K(Me2)-SAPATGGV-K(Me1)-KPHRYRPGG-K(Biotin)

Molecular Weight: 2959.5

Peptide Purity: Peak Area by HPLC ≥95%

Storage: This peptide is shipped at ambient temperature. Upon receipt, store

lyophilized peptide at -20°C or lower. Reconstituted peptide can be

aliquoted and stored at -20°C or lower.

Description: This peptide is Histone 3, with amino acid residues 21 to 44. It is

dimethylated at Lys27 and monomethylated Lys36, with a C-terminal G linker followed by a biotinylated lysine. Histone methylation plays an important role in the regulation of chromatin structure and function.

References:

Jung, HR. et al. Mol Cell Proteomics 9, 838 (2010).

For Research Use Only