



Product Data Sheet

Product Name:	[Lys(Me1) ^{27/36}]-Histone H3 (21-44)-GK(Biotin), H3K27(Me1)K36(Me1), biotin-labeled
Catalog Number:	AS-65446-025 (0.2)
Sequence:	H-Ala-Thr-Lys-Ala-Ala-Arg-Lys -Ser-Ala-Pro-Ala-Thr-Gly-Gly-Val-Lys - Lys-Pro-His-Arg-Tyr-Arg-Pro-Gly-Gly-Lys(Biotin)-OH (3-letter code) ATKAAR-K(Me1)-SAPATGGV-K(Me1)-KPHRYRPG-GK(Biotin) (1-letter code)
Molecular Weight:	2946.5
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 amino acid residues 21 to 44. It is mono-methylated at Lys27 and 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. <i>Mol Cell Proteomics</i> 9 , 838 (2010).

For Research Use Only