



Product Data Sheet

Product Name:	[Lys(Me ₁) ²⁷ , Lys(Me ₂) ³⁶]-Histone H3 (21-44)-GK(Biotin), H3K27(Me1)K36(Me2), biotin-labeled	
Catalog Number:	65447-1 (1 mg); 65447-5 (5 mg)	Lot Number: See label on vial
Sequence:	H-Ala-Thr-Lys-Ala-Ala-Arg-Lys(Me1)-Ser-Ala-Pro-Ala-Thr-Gly-Gly-Val-Lys(Me2)-Lys-Pro-His-Arg-Tyr-Arg-Pro-Gly-Gly-Lys(Biotin)-OH ATKAAR-K(Me1)-SAPATGGV-K(Me2)-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 monomethylated at Lys27 and dimethylated at 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.	
Related Products:	Histone H3 (21-44)-GK(Biotin), biotin-labeled, Cat# 64641	
References:	Jung, HR. et al. <i>Mol Cell Proteomics</i> 9 , 838 (2010).	

For Research Use Only