



Product Data Sheet

Product Name:	[Lys(biotin)79]-Histone H3 (73-83), H3K79(biotin), Biotinylated
Catalog Number:	65440-025 (0.25 mg); 65440-1 (1 mg) Lot Number: See label on vial
Sequence:	H-Glu-Ile-Ala-Gln-Asp-Phe-Lys(Biotin)-Thr-Asp-Leu-Arg-OH (3-letter code) EIAQDF-K(biotin)-TDLR (1-letter code)
Molecular Weight:	1562.8
Peptide Purity:	Peak Area by HPLC $\geq 95\%$
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 H3 amino acid residues 73 to 83, biotinylated at Lysine 79.
Related Products:	Histone H3 (73-83), H3K79 peptide. Cat# 65437
References:	Bheda, P. et al. <i>PNAS</i> 109 , E916 (2012).

For Research Use Only