



## Product Data Sheet

---

<b>Product Name:</b>	Histone H3 (73-83)
<b>Catalog Number:</b>	AS-65437-1 (1 mg); AS-65437-5 (5 mg) Lot Number: See label on vial
<b>Sequence:</b>	H-Glu-Ile-Ala-Gln-Asp-Phe-Lys-Thr-Asp-Leu-Arg-OH (3-letter code)  EIAQDFKTDLR (1-letter code)
<b>Molecular Weight:</b>	1336.5
<b>Peptide Purity:</b>	Peak Area by HPLC $\geq$ 95%
<b>Storage:</b>	This peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at $-20^{\circ}\text{C}$ or lower. Reconstituted peptide can be aliquoted and stored at $-20^{\circ}\text{C}$ or lower.
<b>Description:</b>	This peptide is based on amino acids 73 to 83 of histone H3. Lysine 79 of Histone H3 (H3K79) has been identified as a residue for acetylation or methylation.
<b>Related Products:</b>	[Lys(Ac) <sup>79</sup> ]-Histone H3(73-83), H3K79(Ac), Acetylated, Cat# AS-65438
<b>References:</b>	Bheda, P. et al. <i>PNAS</i> <b>109</b> , E916 (2012).

*For Research Use Only*