

Product Data Sheet

Product Name: [Lys(Ac)⁷⁹]-Histone H3 (69-89), H3K79(Ac), biotin-labeled

Catalog Number: AS-65441-025 (0.2

Lot Number: See label on vial

Sequence: H-Arg-Leu-Val-Arg-Glu-Ile-Ala-Gln-Asp-Phe-Lys(Ac)-Thr-Asp-Leu-Arg-

Phe-Gln-Ser-Ser-Ala-Val-Lys(biotin)-OH (3-letter code)

RLVREIAQDF-K(Ac)-TDLRFQSSAV-K(biotin) (1-letter code)

Molecular Weight: 2877.4

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 a synthetic peptide corresponding to amino acids 69 to 89

of histone H3. It is acetylated at Lys79 and biotinylated on the C-terminal Lys90. The acetylation of histone H3 at Lysine (K) 79 is observed in vivo

in S. cerevisiae.

References: Bheda, P. et al. PNAS 109, E916 (2012).

For Research Use Only