

Product Data Sheet

Product Name: Hexa His

Catalog Number: AS-24420 (1 mg) Lot Number: See label on vial

AS-24421 (5 mg)

Sequence: H-His-His-His-His-His-His-OH (3-letter code)

HHHHHH (1-letter code)

Molecular Weight: 840.9

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Hexa His is freely soluble in water.

Storage: Hexa His 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.

For Research Use Only