



Product Data Sheet

Product Name: Histatin-5

Catalog Number: AS-61001 (1 mg) **Lot Number:** See label on vial

Sequence: H-Asp-Ser-His-Ala-Lys-Arg-His-His-Gly-Tyr-Lys-Arg-Lys-Phe-His-Glu-Lys-His-His-Ser-His-Arg-Gly-Tyr-OH (3-letter code)
DSHAKRHHGYKRKFHEKHHSHRGY (1-letter code)

Molecular Weight: 3037.3

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Histatin-5 peptide is freely soluble in H₂O.

Storage: Histatin-5 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: Histatin 5 is a human basic salivary anti-microbial peptide with strong fungicidal properties. Ref: Helmerhorst, E. et al. *J. Biol. Chem.* **274**, 7286 (1999).

For Research Use Only