

Product Data Sheet

Product Name: LL-37, Antimicrobial Peptide, human

Catalog Number: AS-61302 (1.0 mg) Lot Number: See label on vial

Sequence: H-Leu-Leu-Gly-Asp-Phe-Phe-Arg-Lys-Ser-Lys-Glu-Lys-Ile-Gly-Lys-Glu-

Phe-Lys-Arg-Ile-Val-Gln-Arg-Ile-Lys-Asp-Phe-Leu-Arg-Asn-Leu-Val-Pro-

Arg-Thr-Glu-Ser-OH (3-letter code)

LLGDFFRKSKEKIGKEFKRIVQRIKDFLRNLVPRTES (1-letter code)

Molecular Weight: 4494.3

% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: LL-37 Antimicrobial peptide is freely soluble in H₂O.

Storage: LL-37 Antimicrobial 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: LL - 37, Antimicrobial Peptide, human belongs to the cathelicidin family. It is the first amphipathic alpha-helical peptide isolated from human. It plays an important role in the first line of defense against local infection and systemic invasion of pathogens at sites of inflammation and wounds. Cytotoxic to both bacterial and normal eukaryotic cells, LL-37 is significantly resistant to proteolytic degradation in solution. Ref: Dürr, UHN. et al. *Biochim. Biophys. Acta* 1758, 1408 (2006); Neville, F. et al. *Biophys. J.* 90, 1275 (2006); Oren, Z. et al. *Biochem. J.* 341, 501(1999).

Published Citations:

Bourbon, C. et al. *Analy. Biochem.* **128**, 1077 (2008).

For Research Use Only