



Product Data Sheet

Product Name: Biotin - LC - LL - 37

Catalog Number: AS-63693 (0.5mg) **Lot Number:** See label on vial

Sequence: Biotin-LC-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)
Biotin-LC-LLGDFFRKSKEKIGKEFKRIVQRIKDFLRNLPRTES (1-letter code)

Molecular Weight: 4832.8

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Biotin - LC - LL - 37 is freely soluble in water.

Storage: Biotin - LC - LL - 37 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.

Published Citations:

Into, T. et al. *Cell. Immunol.* **264**, 104 (2010)

Related Products:

Name	Cat #	Size
LL-37, Antimicrobial Peptide, human LLGDFFRKSKEKIGKEFKRIVQRIKDFLRNLPRTES	AS-61302	1 mg
Cys-LC-LL-37 C-LC-LLGDFFRKSKEKIGKEFKRIVQRIKDFLRNLPRTES	AS-63692	1 mg
5-FAM-LC-LL-37 5-FAM-LC-LLGDFFRKSKEKIGKEFKRIVQRIKDFLRNLPRTES	AS-63694	0.1 mg

For Research Use Only