



## Product Data Sheet

---

|                          |                                                                                                                                                                                                                                                                       |                               |
|--------------------------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|-------------------------------|
| <b>Product Name:</b>     | LL17-32                                                                                                                                                                                                                                                               |                               |
| <b>Catalog Number:</b>   | AS-65452-1 (1 mg); AS-65452-5 (5 mg)                                                                                                                                                                                                                                  | Lot Number: See label on vial |
| <b>Sequence:</b>         | H-Phe-Lys-Arg-Ile-Val-Gln-Arg-Ile-Lys-Asp-Phe-Leu-Arg-Asn-Leu-Val-OH<br><br>FKRIVQRIKDFLRNLV                                                                                                                                                                          |                               |
| <b>Molecular Weight:</b> | 2045.5                                                                                                                                                                                                                                                                |                               |
| <b>Peptide Purity:</b>   | Peak Area by HPLC $\geq$ 95%                                                                                                                                                                                                                                          |                               |
| <b>Storage:</b>          | This peptide is shipped at ambient temperature. Upon receipt, store lyophilized peptide at $-20^{\circ}\text{C}$ or lower. Reconstituted peptide can be aliquoted and stored at $-20^{\circ}\text{C}$ or lower.                                                       |                               |
| <b>Description:</b>      | This peptide is an active segment of LL-37, a peptide derived from the C-terminal domain of human cathelicidin antimicrobial peptide. It has been reported that LL17-32 peptide exhibits reversal effect on ABCG2-mediated multidrug resistance in cancer cell lines. |                               |
| <b>Related Products:</b> | LL-37, Antimicrobial Peptide, human, Cat# AS-61302                                                                                                                                                                                                                    |                               |
| <b>References:</b>       | 1. Dürr, UHN. et al. <i>Biochim Biophys Acta</i> <b>1758</b> , 1408 (2006).<br>2. To, KK. et al. <i>Peptides</i> <b>40</b> ,13 (2013).                                                                                                                                |                               |

*For Research Use Only*