

## **Product Data Sheet**

Product Name: LL17-32

Catalog Number: AS-65452-1 (1 mg); AS-65452-5 (5 mg) Lot Number: See label on vial

Sequence: H-Phe-Lys-Arg-Ile-Val-Gln-Arg-Ile-Lys-Asp-Phe-Leu-Arg-Asn-Leu-Val-

ОН

**FKRIVQRIKDFLRNLV** 

Molecular Weight: 2045.5

Peptide Purity: Peak Area by HPLC ≥95%

Storage: This 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: 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.

**Related Products:** 

LL-37, Antimicrobial Peptide, human, Cat# AS-61302

References: 1. Dürr, UHN. et al. Biochim Biophys Acta 1758, 1408 (2006).

2. To, KK. et al. Peptides 40,13 (2013).

For Research Use Only