

## **Product Data Sheet**

Product Name: SHLP3 (Small humanin-like peptide 3)

Catalog Number: AS-65589 (0.1 mg) Lot Number: See label on vial

Sequence: NH<sub>2</sub>-Met-Leu-Gly-Tyr-Asn-Phe-Ser-Ser-Phe-Pro-Cys-Gly-Thr-lle-Ser-Ile-

Ala-Pro-Gly-Phe-Asn-Phe-Tyr-Arg-Leu-Tyr-Phe-Ile-Trp-Val-Asn-Gly-Leu-

Ala-Lys-Val-Val-Trp-COOH

H-MLGYNFSSFPCGTISIAPGFNFYRLYFIWVNGLAKVVW-OH

Molecular Weight: 4380.4

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: Recent advances in high-resolution sequencing have led to the discovery of

unique peptides derived from mitochondrial genome. <sup>1-2</sup> Currently 8 peptides are identified: humanin, mitochondrial open reading frame of the 12S tRNA-c (MOTS-c), and six small humanin-like peptides (SHLP1-6). <sup>1-2</sup> All of these

peptides are released into cytosol from mitochondria. SHLP3 shares

protective effects with Humanin, such as reduction in apoptosis, generation of reactive oxygen species and improving mitochondrial metabolism. In

addition, it may participate in age-related disease pathology<sup>1-2</sup>.

Related Products: Humanin, Cat# AS-60886

MOTS-c, Cat# AS-65587 SHLP-1, Cat# AS-65588 SHLP-4, Cat# AS-65590 SHLP-6, Cat# AS-65591

References:

1. Cobb L. J. et al., AGING 8 (4), 796-808 (2016).

2. Okada A. K. et al., Sci Reports 7 (7802), 1-10 (2017).

For Research Use Only