

## **Product Data Sheet**

Product Name: SHLP6 (Small humanin-like peptide 6)

Catalog Number: AS-65591 (1 mg) Lot Number: See label on vial

Sequence: NH<sub>2</sub>-Met-Leu-Asp-Gln-Asp-Ile-Pro-Met-Val-Gln-Pro-Leu-Leu-Lys-Val-Arg-

Leu-Phe-Asn-Asp-COOH

H-MLDQDIPMVQPLLKVRLFND-OH

Molecular Weight: 2385.9

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. SHLP6 has been shown to increase apoptosis with potential cancer targeting roles and also

may be involved in aging. 1-2

Related Products: Humanin, Cat# AS-60886

MOTS-c, Cat# AS-65587 SHLP-1, Cat# AS-65588 SHLP-3, Cat# AS-65589 SHLP-4, Cat# AS-65590

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