

Product Data Sheet

Product Name: SHLP4 (Small humanin-like peptide 4)

Catalog Number: AS-65590 (1 mg) Lot Number: See label on vial

Sequence: NH₂-Met-Leu-Glu-Val-Met-Phe-Leu-Val-Asn-Arg-Gly-Lys-Ile-Cys-Arg-

Val-Pro-Phe-Thr-Phe-Phe-Asn-Leu-Ser-Leu-COOH

H-MLEVMFLVNRRGKICRVPFTFFNLSL-OH

Molecular Weight: 3131.85

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. 1-2 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). 1-2 All of these peptides are released into cytosol from mitochondria. SHLP4 may promote cell proliferation in certain cells and could be involved in reduced apoptosis 1-

Related Products: Humanin, Cat# AS-60886

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