

## **Product Data Sheet**

Product Name:	SHLP1 (Small humanin-like peptide 1)	
Catalog Number:	AS-65588 (1 mg)	Lot Number: See label on vial
Sequence:	NH <sub>2</sub> -Met-Cys-His-Trp-Ala-Gly-Gly-Ala-Ser-Asn-Thr-Gly-Asp-Ala-Arg-Gly- Asp-Val-Phe-Gly-Lys-Gln-Ala-Gly-COOH H-MCHWAGGASNTGDARGDVFGKQAG-OH	
Molecular Weight:	2393.6	
Peptide Purity:	Peak Area by HPLC ≥95%	
Storage:	This peptide is shipped at ambient temperally ophilized peptide at –20°C or lower. Record aliquoted and stored at –20°C or lower.	
Description:	Recent advances in high-resolution sequer unique peptides derived from mitochondria are identified: humanin, mitochondrial oper c (MOTS-c), and six small humanin-like pe peptides are released into cytosol from mit- increased longevity and cell viability, reduc beneficial functions. <sup>1-2</sup>	I genome. <sup>1-2</sup> Currently 8 peptides n reading frame of the 12S tRNA- ptides (SHLP1-6). <sup>1-2</sup> All of these ochondria and associate with
Related Products:	Humanin, Cat# AS-60886 MOTS-c, Cat# AS-65587 SHLP-3, Cat# AS-65589 SHLP-4, Cat# AS-65590 SHLP-6, Cat# AS-65591	
References:	<ol> <li>Cobb L. J. et al., AGING 8 (4), 796-808</li> <li>Okada A. K. et al., Sci Reports 7 (7802)</li> </ol>	

For Research Use Only