

Product Data Sheet

Product Name: MOTS-c

Catalog Number: AS-65587 (1 mg) Lot Number: See label on vial

Sequence: NH₂-Met-Arg-Trp-Gln-Glu-Met-Gly-Tyr-Ile-Phe-Tyr-Pro-Arg-Lys-Leu-Arg-

COOH

H-MRWQEMGYIFYPRKLR-OH

Molecular Weight: 2174.7

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 and associate with increased longevity and cell viability, reduced apoptosis, and other beneficial functions. ¹⁻³ MOTS-c was found to reduce insulin resistance,

decrease obesity, and promote homeostasis.3

Related Products: Humanin, Cat# AS-60886

SHLP-1, Cat# AS-65588 SHLP-3, Cat# AS-65589 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).

3. Lee C., et al., Cell Met 21, 443-454 (2015).