

Product Data Sheet

Product Name: Nesfatin-1 (24-53), human

Catalog Number: AS-65571 (0.5 mg) Lot Number: See label on vial

Sequence: Pro-Asp-Thr-Gly-Leu-Tyr-Tyr-Asp-Glu-Tyr-Leu-Lys-Gln-Val-Ile-Asp-Val-Leu-

Glu-Thr-Asp-Lys-His-Phe-Arg-Glu-Lys-Leu-Gln-Lys-NH2

PDTGLYYDEYLKQVIDVLETDKHFREKLQK-NH2

Molecular Weight: 3685.3

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: Nesfatin-1 was recently identified as an anorexigenic peptide and is

derived from Nucleobindin2 protein. Nesfatin-1 reduces body weight gain, food and water intake in rodents. In humans high levels of Nesfatin-1 in plasma are associated with overweight individuals. In addition, Nesfatin-1

has been linked to anxiety and stress.

References: 1. Feijóo-Bandín S., et.al., *Endocrinology* **154 (12)**, 4757-4767 (2013)

2. Vas S., et.al., PLOS 8 (4), 1-10 (2013)

3. Ramanjaneya M., et.al., *Endocrinology* **151 (7)**, 3169-3180 (2010)

For Research Use Only