



Product Data Sheet

| | | |
|--------------------------|--|-------------------------------|
| Product Name: | Nesfatin-1 (24-53), mouse | |
| Catalog Number: | AS-65572 (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-Glu-Val-Leu-Glu-Thr-Asp-Pro-His-Phe-Arg-Glu-Lys-Leu-Gln-Lys-NH2 PDTGLYYDEYLKQVIEVLETDPHFREKLQK-NH2 | |
| Molecular Weight: | 3672.3 | |
| 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 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 was linked to the anxiety and stress. | |
| Related Products: | Nesfatin-1, human, Cat# AS-65571 | |
| References: | <ol style="list-style-type: none">1. Feijóo-Bandín S., et.al., <i>Endocrinology</i> 154 (12), 4757-4767 (2013)2. Vas S., et.al., <i>PLOS</i> 8 (4), 1-10 (2013)3. Ramanjaneya M., et.al., <i>Endocrinology</i> 151 (7), 3169-3180 (2010) | |

For Research Use Only