

Product Data Sheet

Product Name: Ghrelin, rat

Catalog Number: AS-24159 (1 mg) Lot Number: See label on vial

AS-24160 (5 mg)

Sequence: H-Gly-Ser-[Ser(n-octanoyl)]-Phe-Leu-Ser-Pro-Glu-His-Gln-Lys-Ala-Gln-

Gln-Arg-Lys-Glu-Ser-Lys-Lys-Pro-Pro-Ala-Lys-Leu-Gln-Pro-Arg-OH

(3-letter code)

GS-S(n-octanoyl)-FLSPEHQKAQQRKESKKPPAKLQPR (1-letter code)

Molecular Weight: 3314.8

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Ghrelin peptide is freely soluble in H₂O.

Storage: Ghrelin 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: Endogenous growth-hormone releasing peptide with novel regulatory mechanism.

Additional Information: Listed below are relevant information that may provide a guideline on how to use this product. End users will have to adapt to their own specific applications.

After arrival (2 wk), rats were randomly assigned to undergo twice daily (8:00 PM and 8:00 AM) subcutaneous injections of rat ghrelin (n = 8, 200 ng/injection; AnaSpec, San Jose, CA) or vehicle (water, n = 8). The ghrelin dose was chosen to be lower but in the same order of magnitude of the one previously reported to markedly increase body weight without affecting food intake in mice, and lack of food intake effects were verified in preliminary short-term experiments-Barazzoni, R. et al. *Am J Physiol Endocrinol Metab* **288**, 228 (2004).

Rat ghrelin, purchased from AnaSpec, Inc. (San Jose, CA, USA) was dissolved in distilled water (1mg/ml), and stored at -20°C until the time of preparation for administration. Immediately before administration by intraperitoneal (i.p.) injection, ghrelin was diluted again with 0.9% saline to the final concentration of 0.1 mg/ml-Liu, Y. et al. *Chinese J. Physio.* 49, 244 (2006).

Cells were treated with lipopolysaccharide (LPS) (0–1 μ g/mL) in the presence or absence of n-octonylated ghrelin (AnaSpec) dissolved in PBS at concentrations ranging from 0 to 1 μ mol/L-Waseem, T. et al. *Surgery* **143**, 334 (2008).

Published Citations:

Barazzoni, R. et al. *Am J Physiol Endocrinol Metab* **288**, 228 (2004). Liu, Y. et al. *Chinese J. Physio.* **49**, 244 (2006). Waseem, T. et al. *Surgery* **143**, 334 (2008). Yang, J. et al. *Cell* **132**, 387 (2008). Granado, M. et al. *Mole. & Cell.* **309**, 67 (2009). Xu, J. et al. *Neurosci. Let.* **453**, 58 (2009).

Related Products:

Name	Cat # Si	ze
[Des-octanoyl]-Ghrelin, rat	AS-60981	1 mg
GSSFLSPEHQKAQQRKESKKPPAKLQPR		
Ghrelin, human	AS-24157	1 mg
GS-S(n-octanoyl)-FLSPEHQRVQQRKESKKPPAKLQPR	AS-24158	5 mg

For Research Use Only