



Product Data Sheet

Product Name: Ghrelin, human
Catalog Number: AS-24157 (1 mg) Lot Number: See label on vial
AS-24158 (5 mg)
Sequence: H-Gly-Ser-[Ser(n-octanoyl)]-Phe-Leu-Ser-Pro-Glu-His-Gln- Arg-Val-Gln-
Gln-Arg-Lys-Glu-Ser-Lys-Lys-Pro-Pro-Ala- Lys-Leu-Gln-Pro-Arg-OH
(3-letter code)
GS-S(n-octanoyl)-FLSPEHQRVQQRKESKKPPAKLQPR (1-letter code)
Molecular Weight: 3371.9
% Peak Area by HPLC: ≥ 95

Appearance: Lyophilized white powder

Peptide Reconstitution: Using H₂O, reconstitute by adding 100 µl to 1 mg Ghrelin peptide.

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: Ghrelin is 28 amino acids, in which the Serine 3 residue is n-octanoylated. This is an endogenous ligand specific for growth hormone secretagogue receptor (GHS-R). Ref: Kojima, M. et al. *Nature* **402**, 656 (1999).

Related Products:

Name	Cat #	Size
Biotin-Ghrelin, human (Biotin-GS-S(n-octanoyl)-FLSPEHQRVQQRKESKKPPAKLQPR)	AS-29391	1 mg
[Des-octanoyl]-Ghrelin, human (GSSFLSPEHQRVQQRKESKKPPAKLQPR)	AS-61177	1 mg
Ghrelin, rat (GS-S(n-octanoyl)-FLSPEHQKAQQRKESKKPPAKLQPR)	AS-24159 AS-24160	1 mg 5 mg
[Des-octanoyl]-Ghrelin, rat (GSSFLSPEHQKAQQRKESKKPPAKLQPR)	AS-60981	1 mg

For Research Use Only