

Product Data Sheet

Product Name: Peptide YY (3-36), human

Catalog Number: AS-24405 (0.5 mg) Lot Number: See label on vial

AS-24406 (1 mg)

Sequence: H-IIe-Lys-Pro-Glu-Ala-Pro-Gly-Glu-Asp-Ala-Ser-Pro-Glu-Glu-Leu-Asn-

Arg-Tyr-Ala-Ser-Leu-Arg-His-Tyr-Leu-Asn-Leu-Val-Thr-Arg-Gln-Arg-

Tyr-NH2 (3-letter code)

IKPEAPGEDASPEELNRYYASLRHYLNLVTRQRY-NH2 (1-letter code)

Molecular Weight: 4049.5

Peptide Purity: >95%

Appearance: Lyophilized white powder

Peptide Reconstitution: Peptide YY is freely soluble in water.

Storage: Peptide YY 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: This peptide, a PYY (3-36), a Y2R agonist, is released from the gastrointestinal tract postprandially in proportion to the calorie content of a meal and can inhibit food intake. Ref: Grandt, D. et al. *Regul. Peptides* **40**, 161 (1992); Batterham, R. et al. *Nature.* **418**, 650 (2002).

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.

Synthetic human PYY[3–36] and acyl-ghrelin were synthesized in-house or purchased from Anaspec, Inc. (San Jose, CA), respectively. The purity and identity of the peptides were confirmed by mass spectroscopy. Peptide stock solutions in deionized water were aliquoted into polypropylene microcentrifuge tubes, lyophilized, and frozen (–20 C) until freshly redissolved in physiological saline on each experimental day-<u>Adams, Sean H. et al. *Endocrinology*</u>. **145**, 4967 (2004).

Published Citations:

Adams, Sean H. et al. *Endocrinology*. **145**, 4967 (2004).

For Research Use Only