



## Product Data Sheet

<b>Product Name:</b>	Peptide YY, human	
<b>Catalog Number:</b>	AS-24401 (0.5 mg) AS-24402 (1 mg)	<b>Lot Number:</b> See label on vial
<b>Sequence:</b>	H-Tyr-Pro-Ile-Lys-Pro-Glu-Ala-Pro-Gly-Glu-Asp-Ala-Ser-Pro-Glu-Glu-Leu-Asn-Arg-Tyr-Tyr-Ala-Ser-Leu-Arg-His-Tyr-Leu-Asn-Leu-Val-Thr-Arg-Gln-Arg-Tyr-NH <sub>2</sub> (3-letter code) YPIKPEAPGEDASPEELNRYYYASLRHYLNLVTRQRY-NH <sub>2</sub> (1-letter code)	
<b>Molecular Weight:</b>	4309.8	
<b>Peptide Purity:</b>	>95%	
<b>Appearance:</b>	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:** PYY is a 36-amino acid regulatory gut hormone present in endocrine cells in the intestine. This peptide acts both as an endocrine and a paracrine agent. Ref: Tatemoto, K. et al. *BBRC* **157**, 713 (1988); Halldén, G. and G. Aponte, *J. Biol. Chem.* **272**, 12591 (1997); Murphy, KG. and SR. Bloom, *Exp. Physiol.* **89**, 507 (2004); Batterham, R. et al. *Nature* **418**, 650 (2002); Gribble, F. et al. *Diabetes* **52**,1147 (2003); Batterham, R. et al. *N. Engl. J. Med.* **349**, 941 (2003).

### Related Products:

<b>Name</b>	<b>Cat #</b>	<b>Size</b>
Peptide YY (3-36), human	AS-24405	0.5 mg
IKPEAPGEDASPEELNRYYYASLRHYLNLVTRQRY-NH <sub>2</sub>	AS-24406	1mg

*For Research Use Only*