



Product Data Sheet

Product Name:	Amylin (1-37), rat	
Catalog Number:	AS-60253-05 (0.5 mg) AS-60253-1 (1 mg)	Lot Number: See label on vial
Sequence:	H-Lys-Cys-Asn-Thr-Ala-Thr-Cys-Ala-Thr-Gln-Arg-Leu-Ala-Asn-Phe-Leu-Val-Arg-Ser-Ser-Asn-Asn-Leu-Gly-Pro-Val-Leu-Pro-Pro-Thr-Asn-Val-Gly-Ser-Asn-Thr-Tyr-NH ₂ (Disulfide bridge: 2-7) (3-letter code) KCNTATCATQRLANFLVRSSNNLGPVLPPTNVGSNTY-NH ₂ (Disulfide bridge: 2-7) (1-letter code)	
Molecular Weight:	3920.5	
Peptide Purity:	>95%	
Appearance:	Lyophilized white powder	

Peptide Reconstitution: Amylin (1-37) is freely soluble in water.

Storage: Amylin (1-37) 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.

Reference: Cooper, GJ. *Endo. Rev.* **15**, 163 (1994).

Related Products:

Name	Cat #	Size
Amylin (8-37), rat ATQRLANFLVRSSNNLGPVLPPTNVGSNTY-NH ₂	AS-21749	1 mg
Amylin (1-37), human KCNTATCATQRLANFLVHSSNCFGAILSSTNVGSNTY (Disulfide bridge: 2-7)	AS-60804	1 mg
Amylin (1-37), human, amide KCNTATCATQRLANFLVHSSNCFGAILSSTNVGSNTY-NH ₂ (Disulfide bridge: 2-7)	AS-60254-1	1 mg

For Research Use Only