



## Product Data Sheet

---

**Product Name:** Amylin (1-37), human

**Catalog Number:** AS-60804 (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-His-Ser-Ser-Asn-Asn-Phe-Gly-Ala-Ile-Leu-Ser-Ser-Thr-Asn-Val-Gly-Ser-Asn-Thr-Tyr-OH (Disulfide bridge: 2-7) (3-letter code)  
KCNTATCATQRLANFLVHSSNFGAILSSTNVGSNTY  
(Disulfide bridge: 2-7) (1-letter code)

**Molecular Weight:** 3904.5

**% Peak Area by HPLC:** ≥ 95

**Appearance:** Lyophilized white powder

**Peptide Reconstitution:** Amylin (1-37) peptide is freely soluble in H<sub>2</sub>O.

**Storage:** Amylin (1-37) 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:** This peptide is a major constituent of protein deposits identified in the Islets of Langerhans of patients with noninsulin-dependent diabetes mellitus. Ref: Jhamandas, J. et al. *J. Neurophysiol.* **89**, 2923 (2003).

*For Research Use Only*