

PRODUCT INFORMATION



Amylin (8-37) (human) (trifluoroacetate salt)

Item No. 24894

Formal Name: L-alanyl-L-threonyl-L-glutamyl-L-arginyl-L-leucyl-L-alanyl-L-asparagyl-L-phenylalanyl-L-leucyl-L-valyl-L-histidyl-L-seryl-L-seryl-L-asparagyl-L-asparagyl-L-phenylalanylglycyl-L-alanyl-L-isoleucyl-L-leucyl-L-seryl-L-seryl-L-threonyl-L-asparagyl-L-valylglycyl-L-seryl-L-asparagyl-L-threonyl-L-tyrosine, trifluoroacetate salt

Synonyms: IAPP (8-37) (human), Islet Amyloid Polypeptide (8-37) (human)

MF: $C_{138}H_{215}N_{41}O_{46} \cdot XCF_3COOH$

FW: 3,184.4

Purity: $\geq 95\%$

Supplied as: A lyophilized powder

Storage: $-20^{\circ}C$

Stability: ≥ 4 years

H — 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 — NH_2

• XCF_3COOH

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Amylin (8-37) (human) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the amylin (8-37) (human) (trifluoroacetate salt) in water. The solubility of amylin (8-37) (human) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Amylin (8-37) is a peptide fragment of amylin (Item Nos. 24274 | 24275).¹ It inhibits osteoblast proliferation, but not bone resorption, induced by full-length human amylin and does not act as an agonist at amylin receptors. Amylin (8-37) induces formation of polymorphic fibrils containing coiled and laterally-associated sheet structures.² It is cytotoxic to RINm5F islet β -cells *in vitro* when used at a concentration of 25 μM , an effect that is inversely correlated with mature fibril content.³

References

1. Cornish, J., Callon, K.E., Lin, C.Q., *et al.* Dissociation of the effects of amylin on osteoblast proliferation and bone resorption. *Am. J. Physiol.* **274**(5 Pt 1), E827-833 (1998).
2. Goldsbury, C., Goldie, K., Pellaud, J., *et al.* Amyloid fibril formation from full-length and fragments of amylin. *J. Struct. Biol.* **130**(2-3), 352-362 (2000).
3. Konarkowska, B., Aitken, J.F., Kistler, J., *et al.* The aggregation potential of human amylin determines its cytotoxicity towards islet β -cells. *FEBS J.* **273**(15), 3614-3624 (2006).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/18/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM