

PRODUCT INFORMATION

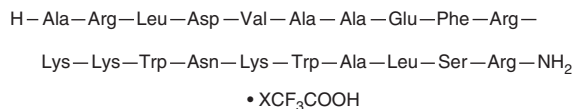


Proadrenomedullin (N-20) (bovine, porcine) (trifluoroacetate salt)

Item No. 27407

Formal Name: L-alanyl-L-arginyl-L-leucyl-L- α -aspartyl-L-valyl-L-alanyl-L-alanyl-L- α -glutamyl-L-phenylalanyl-L-arginyl-L-lysyl-L-lysyl-L-tryptophyl-L-asparaginyl-L-lysyl-L-tryptophyl-L-alanyl-L-leucyl-L-seryl-L-argininamide, trifluoroacetate salt

Synonyms: Mid-Regional proADM, Mid-Regional Proadrenomedullin, MR-proADM, MR-proAM, proADM, ProAM



MF: C₁₁₂H₁₇₈N₃₆O₂₆ • XCF₃COOH

FW: 2,444.8

Purity: \geq 95%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years

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

Laboratory Procedures

Proadrenomedullin (N-20) (bovine, porcine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the proadrenomedullin (N-20) (bovine, porcine) (trifluoroacetate salt) in water. The solubility of proadrenomedullin (N-20) (bovine, porcine) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Proadrenomedullin (N-20) is an endogenous peptide consisting of the 20 N-terminal amino acids of proadrenomedullin, which also corresponds to amino acids 22-41 of preproadrenomedullin.¹⁻³ Proadrenomedullin (N-20) has been found in porcine adrenal medulla.³ The porcine and bovine proadrenomedullin (N-20) amino acid sequence is similar to human proadrenomedullin (N-20) (Item No. 27102) but contains a substitution at position 7.

References

1. Kitamura, K., Kangawa, K., Kojima, M., *et al.* Complete amino acid sequence of porcine adrenomedullin and cloning of cDNA encoding its precursor. *FEBS Lett.* **338(3)**, 306-310 (1994).
2. Barker, S., Wood, E., Clark, A.J., *et al.* Cloning of bovine preproadrenomedullin and inhibition of its basal expression in vascular endothelial cells by staurosporine. *Life Sci.* **62(16)**, 1407-1415 (1998).
3. Kitamura, K., Kangawa, K., Ishiyama, Y., *et al.* Identification and hypotensive activity of proadrenomedullin N-terminal 20 peptide (PAMP). *FEBS Lett.* **351(1)**, 35-37 (1994).

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, 03/21/2023

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