

PRODUCT INFORMATION



Proadrenomedullin (45-92) (human) (trifluoroacetate salt)

Item No. 27209

Formal Name: L- α -glutamyl-L-leucyl-L-arginyl-L-methionyl-L-seryl-L-seryl-L-seryl-L-tyrosyl-L-prolyl-L-threonylglycyl-L-leucyl-L-alanyl-L- α -aspartyl-L-valyl-L-lysyl-L-alanylglycyl-L-prolyl-L-alanyl-L-glutamyl-L-threonyl-L-leucyl-L-isoleucyl-L-arginyl-L-prolyl-L-glutamyl-L- α -aspartyl-L-methionyl-L-lysylglycyl-L-alanyl-L-seryl-L-arginyl-L-seryl-L-prolyl-L- α -glutamyl-L- α -aspartyl-L-seryl-L-seryl-L-prolyl-L- α -aspartyl-L-alanyl-L-alanyl-L-arginyl-L-isoleucyl-L-arginyl-L-valine, trifluoroacetate salt

H — Glu — Leu — Arg — Met — Ser — Ser — Ser — Tyr — Pro — Thr — Gly — Leu — Ala — Asp — Val — Lys — Ala — Gly — Pro — Ala — Gln — Thr — Leu — Ile — Arg — Pro — Gln — Asp — Met — Lys — Gly — Ala — Ser — Arg — Ser — Pro — Glu — Asp — Ser — Ser — Pro — Asp — Ala — Ala — Arg — Ile — Arg — Val — OH
• XCF₃COOH

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

MF: C₂₁₅H₃₅₉N₆₇O₇₃S₂ • XCF₃COOH

FW: 5,114.7

Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥4 years

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

Laboratory Procedures

Proadrenomedullin (45-92) (MR-proADM) (human) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the MR-proADM (human) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. MR-proADM (human) (trifluoroacetate salt) is soluble in the organic solvent formic acid a concentration of approximately 1 mg/ml.

Description

MR-proADM is a peptide that is comprised of amino acid residues 45-92 of preproadrenomedullin.¹ MR-proADM has no effect on precontracted isolated cat pulmonary arterial rings when used at concentrations up to 10 μ M, unlike adrenomedullin (1-52) (Item No. 24889) and adrenotensin, which induce relaxation and contraction, respectively.² Levels of MR-proADM are elevated in the plasma of patients with sepsis and those with cardiovascular disease and positively correlate with disease severity in patients with community-acquired pneumonia.^{1,3}

References

1. Morgenthaler, N.G., Struck, J., Alonso, C., *et al.* Measurement of midregional proadrenomedullin in plasma with an immunoluminometric assay. *Clin. Chem.* **51(10)**, 1823-1829 (2005).
2. Gumusel, B., Chang, J.-K., Hyman, A., *et al.* Adrenotensin: An ADM gene product with the opposite effects of ADM. *Life Sci.* **57(8)**, PL87-PL90 (1995).
3. Huang, D.T., Angus, D.C., Kellum, J.A., *et al.* Midregional proadrenomedullin as a prognostic tool in community-acquired pneumonia. *Chest* **136(3)**, 823-831 (2009).

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, 03/20/2023

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