

PRODUCT INFORMATION



Adrenomedullin (22-52) (human) (trifluoroacetate salt)

Item No. 24892

Formal Name:	L-threonyl-L-valyl-L-glutaminy-L-lysyl-L-leucyl-L-alanyl-L-histidyl-L-glutaminy-L-iso-leucyl-L-tyrosyl-L-glutaminy-L-phenylalanyl-L-threonyl-L- α -aspartyl-L-lysyl-L- α -aspartyl-L-lysyl-L- α -aspartyl-L-asparaginy-L-valyl-L-alanyl-L-prolyl-L-arginyl-L-seryl-L-lysyl-L-iso-leucyl-L-seryl-L-prolyl-L-glutaminyglycyl-L-tyrosinamide, trifluoroacetate salt	H—Thr—Val—Gln—Lys—Leu—Ala—His—Gln—Ile—Tyr—Gln—Phe—Thr—Asp—Lys—Asp—Lys—Asp—Asn—Val—Ala—Pro—Arg—Ser—Lys—Ile—Ser—Pro—Gln—Gly—Tyr—NH ₂
Synonyms:	ADM, AM	
MF:	C ₁₅₉ H ₂₅₂ N ₄₆ O ₄₈ • XCF ₃ COOH	• XCF ₃ COOH
FW:	3,576.0	
Purity:	≥90%	
Supplied as:	A lyophilized powder	
Storage:	-20°C	
Stability:	≥2 years	

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

Laboratory Procedures

Adrenomedullin (22-52) (human) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the adrenomedullin (22-52) (human) (trifluoroacetate salt) in water. The solubility of adrenomedullin (22-52) (human) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Adrenomedullin (22-52) is a C-terminal fragment of adrenomedullin (1-52) (Item No. 24889).¹ *In vitro*, adrenomedullin (22-52) reduces basal corticosterone production in a mixture of rat adrenocortical and adrenomedullary cells.² It also reverses increases in ACTH-stimulated corticosterone production induced by adrenomedullin (1-52). Adrenomedullin (22-52) (0.5 and 5 μ g/kg/min) has no effect on basal regional cerebral blood flow but reverses increases in regional cerebral blood flow induced by rat adrenomedullin in rats.³ Unlike adrenomedullin (1-52), adrenomedullin (22-52) has no effect on mesenteric arterial perfusion pressure in cats.¹

References

1. Tian, Q., Zhao, D., Tan, D.Y., et al. Vasodilator effect of human adrenomedullin (13-52) on hypertensive rats. *Can. J. Physiol. Pharmacol.* **73(7)**, 1065-1069 (1995).
2. Ziolkowska, A., Budzynska, K., Trejter, M., et al. Effects of adrenomedullin and its fragment 22-52 on basal and ACTH-stimulated secretion of cultured rat adrenocortical cells. *Int. J. Mol. Med.* **11(5)**, 613-615 (2003).
3. Dogan, A., Suzuki, Y., Koketsu, N., et al. Intravenous infusion of adrenomedullin and increase in regional cerebral blood flow and prevention of ischemic brain injury after middle cerebral artery occlusion in rats. *J. Cereb. Blood Flow Metab.* **17(1)**, 19-25 (1997).

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