

PRODUCT INFORMATION



Adrenomedullin (1-12) (human) (trifluoroacetate salt)

Item No. 24890

Formal Name:	L-tyrosyl-L-arginyl-L-glutamyl-L-seryl-L-methionyl-L-asparaginyl-L-asparaginyl-L-phenylalanyl-L-glutamylglycyl-L-leucyl-L-arginine, trifluoroacetate salt	
Synonyms:	ADM, AM	H—Tyr—Arg—Gln—Ser—Met—Asn—Asn—Phe—Gln—Gly—
MF:	$C_{64}H_{100}N_{22}O_{19}S \cdot XCF_3COOH$	Leu—Arg—OH
FW:	1,513.7	
Purity:	≥95%	• XCF ₃ COOH
Supplied as:	A lyophilized powder	
Storage:	-20°C	
Stability:	≥4 years	

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

Laboratory Procedures

Adrenomedullin (1-12) (human) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the adrenomedullin (1-12) (human) (trifluoroacetate salt) in water. The solubility of adrenomedullin (1-12) (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 (1-12) is an N-terminal fragment of adrenomedullin (1-52) (Item No. 24889). Unlike adrenomedullin (1-52), adrenomedullin (1-12) has no effect on arterial pressure in cats and lacks antimicrobial activity.^{1,2}

References

1. Lippton, H., Chang, J.K., Hao, Q., *et al.* Adrenomedullin dilates the pulmonary vascular bed *in vivo*. *J. Appl. Physiol.* **76**(5), 2154-2156 (1994).
2. Allaker, R.P., Grosvenor, P.W., McAnerney, D.C., *et al.* Mechanisms of adrenomedullin antimicrobial action. *Peptides* **27**(4), 661-666 (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, 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