

PRODUCT INFORMATION



Adrenomedullin (1-52) (porcine) (trifluoroacetate salt)

Item No. 27406

Formal Name: adrenomedullin (porcine), trifluoroacetate salt
Synonyms: ADM, AM
Peptide Sequence: YRQSMNNFQGLRSFGCRFGTCTVQKLAHQIYQFTDK
DKDGVAPRSKISPQGY-NH₂ (Cys16 and 21 bridge)
MF: C₂₆₂H₄₀₃N₇₉O₇₆S₃ • XCF₃COOH
FW: 5,971.8
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

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

Description

Adrenomedullin (1-52) is a peptide that corresponds to amino acids 1-52 of the porcine adrenomedullin sequence and contains a disulfide bond between cysteine residues at positions 16 and 21. It contains a glycine at position 40 instead of an asparagine residue as in the human adrenomedullin (1-52) (Item No. 24889) sequence.¹ It has been found in the porcine adrenal medulla cortex, pancreatic islet acinus and duct, and the anterior and posterior pituitary.²

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. Washimine, H., Asada, Y., Kitamura, K., *et al.* Immunohistochemical identification of adrenomedullin in human, rat, and porcine tissue. *Histochem. Cell. Biol.* **103(4)**, 251-254 (1995).

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