

PRODUCT INFORMATION



Angiotensinogen (1-13) (human) (trifluoroacetate salt)

Item No. 24988

Formal Name: L- α -aspartyl-L-arginyl-L-valyl-L-tyrosyl-L-
isoleucyl-L-histidyl-L-prolyl-L-phenylalanyl-
L-histidyl-L-leucyl-L-valyl-L-isoleucyl-L-
histidine, trifluoroacetate salt
MF: C₇₉H₁₁₆N₂₂O₁₇ • XCF₃COOH
FW: 1,645.9
Purity: \geq 95%
Supplied as: A lyophilized powder
Storage: -20°C
Stability: \geq 4 years

H—Asp—Arg—Val—Tyr—Ile—His—Pro—Phe—His—Leu—
Val—Ile—His—OH
• XCF₃COOH

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

Laboratory Procedures

Angiotensinogen (1-13) (human) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the angiotensinogen (1-13) (human) (trifluoroacetate salt) in water. The solubility of angiotensinogen (1-13) (human) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Angiotensinogen (1-13) is a synthetic peptide precursor of angiotensin I (Item No. 24737) that corresponds to amino acids 1-13 of endogenous human angiotensinogen.¹ It is cleaved by renin at the Leu-Val peptide bond to release angiotensin I, which is then converted to angiotensin II (Item No. 17150) by angiotensin converting enzyme (ACE) and exerts vasoconstrictive properties. Angiotensinogen (1-13) has been used in the production of anti-angiotensinogen monoclonal antibodies.²

References

1. Tewksbury, D.A., Dart, R.A., and Travis, J. The amino terminal amino acid sequence of human angiotensinogen. *Biochem. Biophys. Res. Commun.* **99(4)**, 1311-1315 (1981).
2. Sawa, H., Kawaguchi, H., Mochizuki, N., *et al.* Distribution of angiotensinogen in diseased human hearts. *Mol. Cell Biochem.* **132(1)**, 15-23 (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, 12/05/2022

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