

PRODUCT INFORMATION



Urocortin II (human) (trifluoroacetate salt)

Item No. 24746

CAS Registry No.: 398001-88-2

Formal Name: L-isoleucyl-L-valyl-L-leucyl-L-seryl-L-leucyl-L- α -aspartyl-L-valyl-L-prolyl-L-isoleucylglycyl-L-leucyl-L-leucyl-L-glutamyl-L-isoleucyl-L-leucyl-L-leucyl-L- α -glutamyl-L-glutamyl-L-alanyl-L-arginyl-L-alanyl-L-arginyl-L-alanyl-L-alanyl-L-arginyl-L- α -glutamyl-L-glutamyl-L-alanyl-L-threonyl-L-threonyl-L-asparagyl-L-alanyl-L-arginyl-L-isoleucyl-L-leucyl-L-alanyl-L-arginyl-L-valylglycyl-L-histidyl-L-cysteine, trifluoroacetate salt

H-Ile-Val-Leu-Ser-Leu-Asp-Val-Pro-Ile-Gly-Leu-Leu-Gln-Ile-Leu-Leu-Glu-Gln-Ala-Arg-Ala-Arg-Ala-Ala-Arg-Glu-Gln-Ala-Thr-Thr-Asn-Ala-Arg-Ile-Leu-Ala-Arg-Val-Gly-His-Cys-NH₂

MF: C₁₉₄H₃₃₉N₆₃O₅₄S • XCF₃COOH

FW: 4,450.3

• XCF₃COOH

Purity: \geq 95%

Supplied as: A lyophilized powder

Storage: -20°C

Stability: \geq 4 years

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

Laboratory Procedures

Urocortin II (human) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the urocortin II (human) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Urocortin II (human) (trifluoroacetate salt) is soluble in the organic solvent formic acid at a concentration of approximately 1 mg/ml.

Description

Urocortin II is a neuropeptide hormone and member of the corticotropin-releasing factor (CRF) family which includes mammalian CRF, urocortin (Item No. 24745), urocortin III (Item Nos. 24749 | 24750), frog sauvagine, and piscine urotensin I.¹ Human urocortin II shares 34, 43, and 37-40% sequence homology with rat and human CRF, human urocortin (Item No. 24744), and human urocortin III (Item No. 24748). Urocortin II increases cell shortening and accelerates relaxation of rabbit ventricular myocytes in a time- and concentration-dependent manner. *In vivo*, urocortin reduces arterial blood pressure in normotensive and spontaneously hypertensive rats *via* peripheral CRF₂ receptor agonism. It induces dose-dependent tachycardia and hypotension in rats when administered at doses of 3 and 30 pmol/kg.² Urocortin (10 and 20 μ g/kg) also reduces the visceral pain response to colorectal distension in conscious rats and delays gastric emptying in mice.^{1,3}

References

1. Adão, R., Santos-Ribeiro, D., Rademaker, M.T., *et al.* Urocortin 2 in cardiovascular health and disease. *Drug Discov. Today* **20(7)**, 906-914 (2015).
2. Gardiner, S.M., March, J.E., Kemp, P.A., *et al.* Regional hemodynamic actions of selective corticotropin-releasing factor type 2 receptor ligands in conscious rats. *J. Pharmacol. Exp. Ther.* **312(1)**, 53-60 (2005).
3. Million, M., Wang, L., Wang, Y., *et al.* CRF₂ receptor activation prevents colorectal distension induced visceral pain and spinal ERK1/2 phosphorylation in rats. *Gut* **55(2)**, 172-181 (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, 11/29/2022

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