

PRODUCT INFORMATION



Sauvagine (trifluoroacetate salt)

Item No. 34442

Formal Name: 5-oxo-L-prolylglycyl-L-prolyl-L-prolyl-L-isoleucyl-L-seryl-L-isoleucyl-L- α -aspartyl-L-leucyl-L-seryl-L-leucyl-L- α -glutamyl-L-leucyl-L-leucyl-L-arginyl-L-lysyl-L-methionyl-L-isoleucyl-L- α -glutamyl-L-isoleucyl-L- α -glutamyl-L-lysyl-L-glutamyl-L- α -glutamyl-L-lysyl-L- α -glutamyl-L-lysyl-L-glutamyl-L-glutamyl-L-alanyl-L-alanyl-L-asparaginy-L-asparaginy-L-arginyl-L-leucyl-L-leucyl-L-leucyl-L- α -aspartyl-L-threonyl-L-isoleucinamide, trifluoroacetate salt

MF: $C_{202}H_{346}N_{56}O_{63}S \cdot XCF_3COOH$

FW: 4,599.4

Purity: $\geq 95\%$

Supplied as: A solid

Storage: $-20^\circ C$

Stability: ≥ 4 years

H—Pyr—Gly—Pro—Pro—Ile—Ser—Ile—Asp—Leu—Ser—
Leu—Glu—Leu—Leu—Arg—Lys—Met—Ile—Glu—Ile—
Glu—Lys—Gln—Glu—Lys—Glu—Lys—Gln—Gln—Ala—
Ala—Asn—Asn—Arg—Leu—Leu—Leu—Asp—Thr—Ile—NH₂

• XCF₃COOH

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

Laboratory Procedures

Sauvagine (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the sauvagine (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Sauvagine is a neuropeptide hormone that has been found in frog (*P. sauvagei*) skin and member of the corticotropin-releasing factor (CRF) family, which includes mammalian CRF, urocortin (Item No. 24745), urocortin II (Item No. 24746), urocortin III (Item Nos. 24749 | 24750), and piscine urotensin 1.^{1,2} *In vivo*, sauvagine (0.5 $\mu g/kg$, s.c.) reduces systemic blood pressure in normotensive rats.² It inhibits anterior pituitary release of prolactin, thyroid-stimulating hormone (TSH), and growth hormone in rats. Sauvagine also increases plasma adrenocorticotrophic hormone (ACTH) levels in rats.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. Montecucchi, P.C., and Henschen, A. Amino acid composition and sequence analysis of sauvagine, a new active peptide from the skin of *Phyllomedusa sauvagei*. *Int. J. Pept. Protein Res.* **18(2)**, 113-120 (1981).
3. Rivier, C., Rivier, J., Lederis, K., et al. In vitro and in vivo ACTH-releasing activity of ovine CRF, sauvagine and urotensin I. *Regul. Pept.* **5(2)**, 139-143 (1983).

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