

PRODUCT INFORMATION

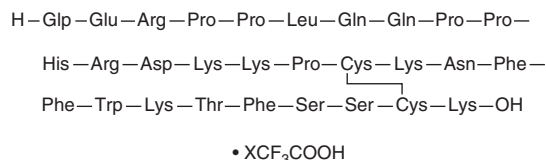


Cortistatin-29 (rat) (trifluoroacetate salt)

Item No. 24740

CAS Registry No.: 1815618-17-7

Formal Name: 5-oxo-L-prolyl-L- α -glutamyl-L-arginyl-L-prolyl-L-prolyl-L-leucyl-L-glutamyl-L-glutamyl-L-prolyl-L-prolyl-L-histidyl-L-arginyl-L- α -aspartyl-L-lysyl-L-lysyl-L-prolyl-L-cysteinyl-L-lysyl-L-asparaginyl-L-phenylalanyl-L-phenylalanyl-L-tryptophyl-L-lysyl-L-threonyl-L-phenylalanyl-L-seryl-L-seryl-L-cysteinyl-L-lysine, cyclic (17 $\rightarrow$ 28)-disulfide, trifluoroacetate salt



MF: C₁₆₁H₂₄₀N₄₆O₄₁S₂ • XCF₃COOH

FW: 3,540.0

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

Cortistatin-29 (rat) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the cortistatin-29 (rat) (trifluoroacetate salt) in water. The solubility of cortistatin-29 (rat) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Cortistatin-29 is a neuropeptide that is structurally similar to somatostatin-28.¹ It is produced by cleavage of precortistatin to procortistatin, which is cleaved at dibasic amino acids to form cortistatin-29 and cortistatin-14 as well as other partial cleavage products. Cortistatin mRNA is expressed in the human brain and in interneurons of the rat hippocampus and cerebral cortex.^{1,2} Cortistatin-29 binds to somatostatin (SST) receptors with IC₅₀ values of 2.8, 7.1, 0.2, 3, and 13.7 nM for SST₁₋₅, respectively.² Cortistatin-29 is found at similar levels as cortistatin-14 in mouse AtT20 cells but is secreted at a lower level. Cortistatin-29 corresponds to residues 85-112 of the rat peptide sequence.

References

1. de Lecea, L., Criado, J.R., Prospero-Garcia, O., *et al.* A cortical neuropeptide with neuronal depressant and sleep-modulating properties. *Nature* **381**(6579), 242-245 (1996).
2. Spier, A.D. and de Lecea, L. Cortistatin: A member of the somatostatin neuropeptide family with distinct physiological functions. *Brain Res. Brain Res. Rev.* **33**(2-3), 228-241 (2000).

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, 11/17/2022

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