

# PRODUCT INFORMATION



## CTX-1211 (trifluoroacetate salt)

Item No. 42081

**Formal Name:** N2-acetyl-L-arginyl-L-cysteinyl-D-alanyl-L-arginyl-D-phenylalanyl-L-arginyl-L-tryptophyl-L-cysteinamide, cyclic (2→8)-disulfide, trifluoroacetate salt

**Synonym:** Acetyl-Arg-cyclo(Cys-D-Ala-Arg-D-Phe-Arg-Trp-Cys)-NH<sub>2</sub>

**MF:** C<sub>49</sub>H<sub>73</sub>N<sub>19</sub>O<sub>9</sub>S<sub>2</sub> • XCF<sub>3</sub>COOH

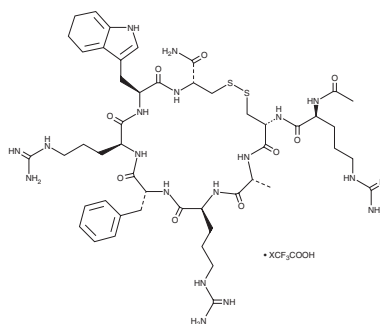
**FW:** 1,136.4

**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

CTX-1211 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the CTX-1211 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. CTX-1211 (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in ethanol.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of CTX-1211 (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. CTX-1211 (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

### Description

CTX-1211 is a peptide agonist of melanocortin receptors.<sup>1</sup> It increases intracellular cAMP levels in cells expressing melanocortin receptor 1 (MC1R), MC3R, or MC4R (EC<sub>50</sub>s = 6.2, 3.9, and 2.7, nM, respectively). CTX-1211 (1 mg/kg) sensitizes male and female mice to reductions in food intake and body weight induced by liraglutide (Item No. 24727) but with a lower effect on food intake in female mice.<sup>2</sup>

### References

1. Sharma, S. Peptide compositions. *Rhythm Metabolic Inc.* **WO 2014/144260 A1** (2014).
2. Dahir, N.S., Gui, Y., Wu, Y., *et al.* Subthreshold activation of the melanocortin system causes generalized sensitization to anorectic agents in mice. *J. Clin. Invest.* **134(14)**, e178250 (2024).

#### 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, 10/30/2024

#### 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