

PRODUCT INFORMATION



CJC-1295 (acetate)

Item No. 32704

Formal Name: L-tyrosyl-D-alanyl-L- α -aspartyl-L-alanyl-L-isoleucyl-L-phenylalanyl-L-threonyl-L-glutamyl-L-seryl-L-tyrosyl-L-arginyl-L-lysyl-L-valyl-L-leucyl-L-alanyl-L-glutamyl-L-leucyl-L-seryl-L-alanyl-L-arginyl-L-lysyl-L-leucyl-L-leucyl-L-glutamyl-L- α -aspartyl-L-isoleucyl-L-leucyl-L-seryl-L-argininamide, acetate

Synonyms: CJC-1295-no DAC, GHRH (1-29)-NH₂, Modified GRF (1-29)

MF: C₁₅₂H₂₅₂N₄₄O₄₂ • XC₂H₄O₂

FW: 3,367.9

Purity: ≥95%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥4 years

H—Tyr—D-Ala—Asp—Ala—Ile—Phe—Thr—Gln—Ser—Tyr—

Arg—Lys—Val—Leu—Ala—Gln—Leu—Ser—Ala—Arg—

Lys—Leu—Leu—Gln—Asp—Ile—Leu—Ser—Arg—NH₂

• XCH₃CO₂H

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

Laboratory Procedures

CJC-1295 (acetate) is supplied as a crystalline solid. A stock solution may be made by dissolving the CJC-1295 (acetate) in the solvent of choice, which should be purged with an inert gas. CJC-1295 (acetate) is soluble in the organic solvent DMSO. The solubility of CJC-1295 (acetate) in DMSO is approximately 1 mg/ml. CJC-1295 (acetate) is slightly soluble in ethanol and dimethyl formamide.

Description

CJC-1295 is a synthetic peptide derivative of growth hormone-releasing hormone (GHRH).¹ Unlike CJC-1295-DAC, CJC-1295 does not contain a drug affinity complex (DAC) moiety and has a shorter half-life than the DAC-conjugated version in the blood.² CJC-1295 increases the secretion of growth hormone in primary rat anterior pituitary cells when used from picomolar to micromolar concentrations and into rat blood when administered subcutaneously at a dose of 1 μ mol/kg.²

References

1. Bridon, D.P., Boudjellab, N., Leger, R., *et al.* Long lasting growth hormone releasing factor derivatives. *Conjuchem Inc.* **WO02/062844A2** (2002).
2. Teichman, S.L., Neale, A., Lawrence, B., *et al.* Prolonged stimulation of growth hormone (GH) and insulin-like growth factor I secretion by CJC-1295, a long-acting analog of GH-releasing hormone, in healthy adults. *J. Clin. Endocrinol. Metab.* **91**(3), 799-805 (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, 08/06/2024

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