

PRODUCT INFORMATION



CJC-1295-DAC (trifluoroacetate)

Item No. 31497

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-arginyl-N-[3-(2,5-dihydro-2,5-dioxo-1H-pyrrol-1-yl)-1-oxopropyl]-L-lysineamide, trifluoroacetate

Synonyms: DAC-GRF

Peptide Sequence: YaDAIFTQSYRKVLAQLSARKLLQDILSRL-NH₂

MF: C₁₆₅H₂₆₉N₄₇O₄₆ • XCF₃COOH

FW: 3,647.3

Purity: $\geq 98\%$

Supplied as: A solid

Storage: -20°C

Stability: ≥ 4 years

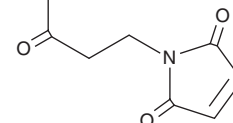
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

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-Lys-NH₂

• XCF₃COOH



Laboratory Procedures

CJC-1295-DAC (trifluoroacetate) is supplied as a solid. A stock solution may be made by dissolving the CJC-1295-DAC (trifluoroacetate) in the solvent of choice, which should be purged with an inert gas. CJC-1295-DAC (trifluoroacetate) is sparingly soluble (1-10 mg/ml) in DMSO.

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 CJC-1295-DAC (trifluoroacetate) can be prepared by directly dissolving the solid in aqueous buffers. CJC-1295-DAC (trifluoroacetate) 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

CJC-1295-DAC is a synthetic peptide derivative of growth hormone-releasing hormone (GHRH) that contains a maleimidopropionic acid (MPA) moiety, which functions as a drug affinity complex (DAC).¹ The DAC binds serum albumin, increasing the stability of the bioactive peptide in rat plasma. CJC-1295-DAC increases plasma GH levels in rats and increases pituitary GH levels, serum insulin-like growth factor (IGF-1) levels, body weight, and body length in GHRH knockout mice.^{1,2}

References

1. Jetté, L., Léger, R., Thibaut, K., *et al.* Human growth hormone-releasing factor (hGRF)₁₋₂₉-albumin bioconjugates activate the GRF receptor on the anterior pituitary in rats: Identification of CJC-1295 as a long-lasting GRF analog. *Endocrinology* **146**(7), (2005).
2. Alba, M., Fintini, D., Sagazio, A., *et al.* Once-daily administration of CJC-1295, a long-acting growth hormone-releasing hormone (GHRH) analog, normalizes growth in the GHRH knockout mouse. *Am. J. Physiol. Endocrinol. Metab.* **291**(6), E1290-1294 (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, 07/31/2026

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