

PRODUCT INFORMATION



KISS1-305 (trifluoroacetate salt)

Item No. 41456

Formal Name: D-tyrosyl-3-(4-pyridinyl)-D-alanyl-L-asparaginyl-L-seryl-L-phenylalanyl-2-azaglycyl-L-leucyl-N-[imino(methylamino)methyl]-L-ornithyl-L-phenylalaninamide, trifluoroacetate salt

Synonym: D-Tyr-D-Pya(4)-Asn-Ser-Phe-azaGly-Leu-Arg(Me)-Phe-NH₂

Peptide Sequence: yxNSFZLBF-NH₂ (x = D-3-(4-Pyridinyl)-alanine; Z = 2-Azaglycine; B = N-methylarginine)

MF: C₅₆H₇₆N₁₆O₁₂ • XCF₃COOH

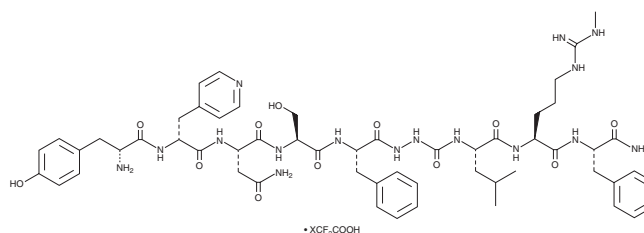
FW: 1,165.3

Purity: ≥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.

Laboratory Procedures

KISS1-305 (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the KISS1-305 (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. KISS1-305 (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in acetonitrile and 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 KISS1-305 (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. KISS1-305 (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

KISS1-305 is a peptide agonist of the kisspeptin receptor.¹ It induces calcium mobilization in a reporter assay using CHO cells expressing the kisspeptin receptor (EC₅₀ = 2.6 nM for the human receptor). KISS1-305 reduces plasma levels of testosterone below the limit of detection in male rats. It increases the number of ovulated oocytes in horse chorionic gonadotropin-stimulated immature female rats when administered at a dose of 100 nmol/kg.²

References

1. Asami, T., Nishizawa, N., Matsui, H., *et al.* Physicochemically and pharmacokinetically stable nonapeptide KISS1 receptor agonists with highly potent testosterone-suppressive activity. *J. Med. Chem.* **57**(14), 6105-6115 (2014).
2. Asami, T., Nishizawa, N., Matsui, H., *et al.* Design, synthesis, and biological evaluation of novel investigational nonapeptide KISS1R agonists with testosterone-suppressive activity. *J. Med. Chem.* **56**(21), 8298-8307 (2013).

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, 09/05/2024

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