

PRODUCT INFORMATION



Kisspeptin-10 (mouse, rat)

Item No. 14036

CAS Registry No.: 478507-53-8

Formal Name: L-tyrosyl-L-asparaginyl-L-tryptophyl-L-asparaginyl-L-seryl-L-phenylalanyl-glycyl-L-leucyl-L-arginyl-L-tyrosinamide

Synonyms: Kp-10, YNWNFSGLRY-NH₂

MF: C₆₃H₈₃N₁₇O₁₅

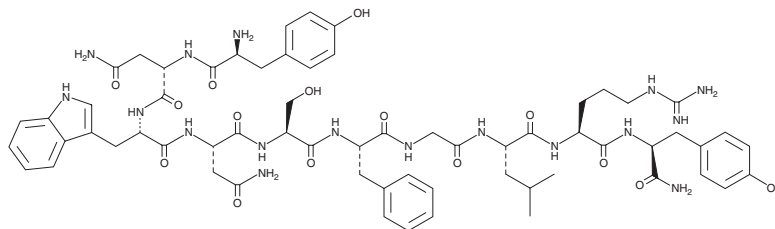
FW: 1,318.4

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

Kisspeptin-10 (mouse, rat) is supplied as a solid. A stock solution may be made by dissolving the kisspeptin-10 (mouse, rat) in water. We do not recommend storing the aqueous solution for more than one day.

Description

Kisspeptin-10 is a neuropeptide and ligand of the kisspeptin receptor, previously known as G protein-coupled receptor 54 (GPR54).¹ It corresponds to amino acids 110-119 of rodent kisspeptin-1 (KiSS1), a metastasis suppressor gene product. Kisspeptin-10 (10 μM) inhibits FGF2-induced migration of NIH3T3 cells expressing the mouse kisspeptin receptor. Intracerebroventricular administration of kisspeptin-10 increases serum follicle-stimulating hormone (FSH) and luteinizing hormone (LH) levels in adult male rats (ED₅₀s = 400 and 4 pmol/animal).²

References

1. Stafford, L.J., Xia, C., Ma, W., *et al.* Identification and characterization of mouse metastasis-suppressor KiSS1 and its G-protein-coupled receptor. *Cancer Res.* **62(19)**, 5399-5404 (2002).
2. Navarro, V.M., Castellano, J.M., Fernández-Fernández, R., *et al.* Effects of KiSS-1 peptide, the natural ligand of GPR54, on follicle-stimulating hormone secretion in the rat. *Endocrinology* **146(4)**, 1689-1697 (2005).

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, 04/18/2023

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