

# PRODUCT INFORMATION



## Kisspeptin-10 (zebrafish) (trifluoroacetate salt)

Item No. 35652

**Formal Name:** L-tyrosyl-L-asparaginyl-L-leucyl-L-asparaginyl-L-seryl-L-phenylalanyl-glycyl-L-leucyl-L-arginyl-L-tyrosinamide, trifluoroacetate salt

**Synonyms:** Kp-10, Tyr-Asn-Leu-Asn-Ser-Phe-Gly-Leu-Arg-Tyr-NH<sub>2</sub>

**Peptide Sequence:** YNLNSFGLRY-NH<sub>2</sub>

**MF:** C<sub>58</sub>H<sub>84</sub>N<sub>16</sub>O<sub>15</sub> • XCF<sub>3</sub>COOH

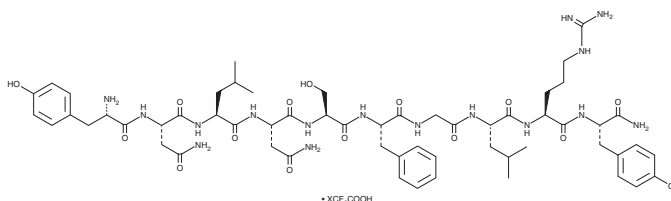
**FW:** 1,245.4

**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 (zebrafish) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the kisspeptin-10 (zebrafish) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Kisspeptin-10 (zebrafish) (trifluoroacetate salt) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of kisspeptin-10 (zebrafish) (trifluoroacetate salt) in these solvents is approximately 20 and 10 mg/ml, respectively.

### Description

Kisspeptin-10 is a neuropeptide expressed in zebrafish brains, ovaries, and testes and is involved in gonadal development and steroid expression.<sup>1</sup> It increases levels of the mRNA encoding luteinizing hormone (Lh) receptor, as well as progesterin receptor a (Pra) and Prb, in isolated zebrafish ovarian follicles when used at a concentration of 10 ng/ml. Kisspeptin-10 increases luteinizing hormone serum levels in female, but not male, goldfish when administered at doses of 0.1 and 1 µg/g.<sup>2</sup>

### References

1. Hatef, A., Rajeswari, J.J., and Unniappan, S. Kisspeptin stimulates oocyte maturation, and food deprivation modulates the abundance of kisspeptin system in zebrafish gonads. *Aquacult. Fish.* **7(5)**, 484-493 (2022).
2. Li, S., Zhang, Y., Liu, Y., et al. Structural and functional multiplicity of the kisspeptin/GPR54 system in goldfish (*Carassius auratus*). *J. Endocrinol.* **201(3)**, 407-418 (2009).

#### 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, 10/31/2025

#### 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