

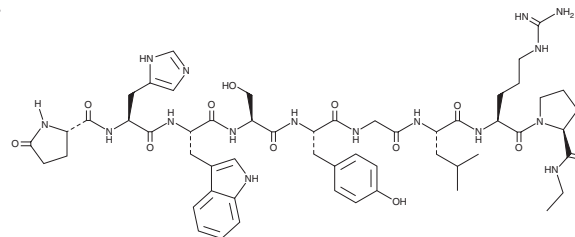
PRODUCT INFORMATION



Fertirelin

Item No. 42967

CAS Registry No.: 38234-21-8
Formal Name: 9-(N-ethyl-L-prolinamide)-1-9-luteinizing hormone-releasing factor (swine)
Synonyms: pGlu-His-Trp-Ser-Tyr-Gly-Leu-Arg-Pro-ethylamide, pGlu-HWSYGLRP-ethylamide, TAP-031, U-9689
MF: $C_{55}H_{76}N_{16}O_{12}$
FW: 1,153.3
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.

Laboratory Procedures

Fertirelin is supplied as a solid. A stock solution may be made by dissolving the fertirelin in the solvent of choice, which should be purged with an inert gas. Fertirelin 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 fertirelin can be prepared by directly dissolving the solid in aqueous buffers. Fertirelin is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Fertirelin is a peptide agonist of the gonadotropin-releasing hormone receptor (GnRHR) and a derivative of GnRH.¹ It binds to GnRHR in CHO cells expressing the human receptor ($K_i = 0.7$ nM).¹ *In vivo*, fertirelin (200 $\mu\text{g}/\text{animal}$) increases serum levels of luteinizing hormone and induces luteinization of follicular cysts in dairy cows.²

References

1. Nederpelt, I., Georgi, V., Schiele, F., *et al.* Characterization of 12 GnRH peptide agonists - a kinetic perspective. *Br. J. Pharmacol.* **173**(1), 128-141 (2016).
2. Osawa, T., Nakao, T., Kimura, M., *et al.* Fertirelin and buserelin compared by LH release, milk progesterone and subsequent reproductive performance in dairy cows treated for follicular cysts. *Theriogenology* **44**(6), 835-847 (1995).

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, 04/10/2025

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