

PRODUCT INFORMATION

GnRH (free acid; trifluoroacetate salt)

Item No. 24762

Formal Name:	10-glycine-luteinizing hormone-releasing factor (swine), trifluoroacetate salt	
Synonyms:	Gonadorelin, Gonadotropin-releasing Hormone, LHRH	H—Glp—His—Trp—Ser—Tyr—Gly—Leu—Arg—Pro—Gly—OH
MF:	C ₅₅ H ₇₄ N ₁₆ O ₁₄ • XCF ₃ COOH	
FW:	1,183.3	
Purity:	≥95%	• XCF ₃ COOH
Supplied as:	A lyophilized powder	
Storage:	-20°C	
Stability:	≥4 years	

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

Laboratory Procedures

GnRH (free acid; trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the GnRH (free acid; trifluoroacetate salt) in water. The solubility of GnRH (free acid; trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

GnRH is a peptide hormone produced in the hypothalamus that is secreted in a pulsatile fashion and stimulates the release of luteinizing hormone and follicle-stimulating hormone.¹ The free acid form of GnRH is biologically inactive.²

References

1. Conn, P.M. and Crowley, W.F., Jr. Gonadotropin-releasing hormone and its analogs. *Annu. Rev. Med.* **45**, 391-405 (1994).
2. Meidan, R. and Koch, Y. Binding of luteinizing hormone-releasing hormone analogues to dispersed rat pituitary cells. *Life Sci.* **28(17)**, 1961-1967 (1981).

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, 11/10/2022

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