

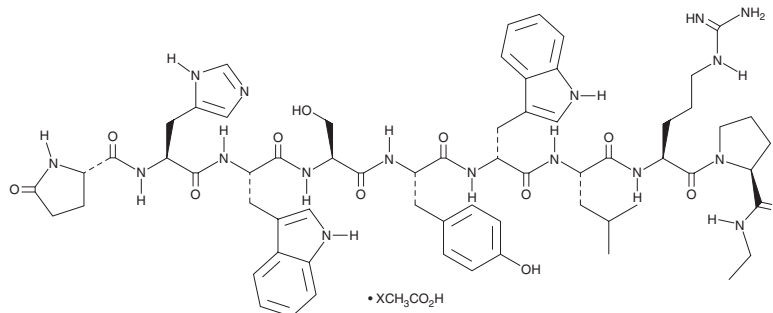
PRODUCT INFORMATION



Deslorelin (acetate)

Item No. 34041

CAS Registry No.: 1642579-30-3
Formal Name: 6-D-tryptophan-9-(N-ethyl-L-prolinamide)-1-9-luteinizing hormone-releasing factor (swine), acetate
Synonyms: D-Trp⁶-Pro⁹-des-Gly¹⁰-GnRH ethylamide, pGlu-His-Trp-Ser-Tyr-D-Trp-Leu-Arg-Pro-NH₂
MF: C₆₄H₈₃N₁₇O₁₂ • XC₂H₄O₂
FW: 1,282.5
Purity: ≥98%
UV/Vis.: λ_{max}: 219 nm
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

Deslorelin (acetate) is supplied as a solid. A stock solution may be made by dissolving the deslorelin (acetate) in the solvent of choice, which should be purged with an inert gas. Deslorelin (acetate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of deslorelin (acetate) in ethanol is approximately 10 mg/ml and approximately 5 mg/ml in DMSO and DMF.

Description

Deslorelin is a peptide agonist of gonadotropin-releasing hormone (GnRH).¹ It binds to the GnRH receptor (GnRHR) with a K_i value of 0.1 nM. Deslorelin decreases testicular volume, semen volume, sperm motility, and total sperm count, as well as increases the percentage of abnormal sperm, in dogs when administered via a subcutaneous implant at doses of 3, 6, and 12 mg/kg.² It decreases disease severity and plasma levels of estradiol, 17-hydroxyprogesterone, and androstenedione in ferrets with adrenocortical disease (ACD) when administered at doses ranging from 3 to 4.41 mg/kg per implant.³ Deslorelin (4.7 mg/implant per animal) suppresses ovulation and estrus in cats.⁴ Formulations containing deslorelin have been used in the management of ACD in ferrets and ovulation regulation and artificial insemination in horses.

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. Junaidi, A., Williamson, P.E., Martin, G.B., *et al.* Dose-response studies for pituitary and testicular function in male dogs treated with the GnRH superagonist, deslorelin. *Reprod. Domest. Anim.* **44**(5), 725-734 (2009).
3. Wagner, R.A., Piché, C.A., Jöchle, W., *et al.* Clinical and endocrine responses to treatment with deslorelin acetate implants in ferrets with adrenocortical disease. *Am. J. Vet. Res.* **66**(5), 910-914 (2005).
4. Ackermann, C.L., Volpato, R., Destro, F.C., *et al.* Ovarian activity reversibility after the use of deslorelin acetate as a short-term contraceptive in domestic queens. *Theriogenology* **78**(4), 817-822 (2012).

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, 11/16/2022

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