

PRODUCT INFORMATION

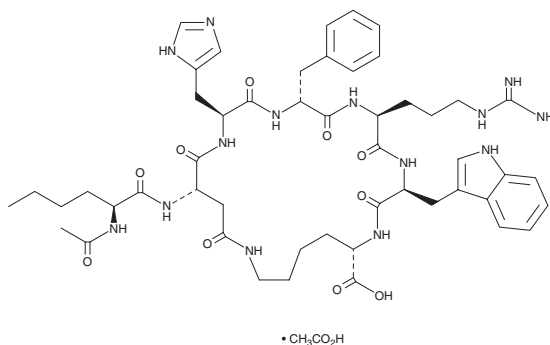


PT-141 (acetate)

Item No. 35175

CAS Registry No.: 1607799-13-2
Formal Name: N-acetyl-L-norleucyl-L- α -aspartyl-L-histidyl-D-phenylalanyl-L-arginyl-L-tryptophyl-L-lysine (2 \rightarrow 7)-lactam, monoacetate

Synonym: Bremelanotide Acetate
MF: C₅₀H₆₈N₁₄O₁₀ • C₂H₄O₂
FW: 1,085.2
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

PT-141 (acetate) is supplied as a solid. A stock solution may be made by dissolving the PT-141 (acetate) in the solvent of choice, which should be purged with an inert gas. PT-141 (acetate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of PT-141 (acetate) in these solvents is approximately 10 mg/ml.

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 PT-141 (acetate) can be prepared by directly dissolving the solid in aqueous buffers. The solubility of PT-141 (acetate) in PBS (pH 7.2) is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

PT-141 is a cyclic peptide agonist of melanocortin receptor 4 (MC4R) and an analog of α -melanocyte stimulating hormone (α -MSH; Item No. 29923).¹ It binds to MC3R and MC4R and increases intracellular cAMP levels in HEK293 cells expressing MC4R. It increases dopamine release in the medial preoptic area (mPOA) of the hypothalamus in female rats.² PT-141 (100 and 200 $\mu\text{g}/\text{kg}$) increases the number of solicitations, an appetitive sexual behavior, by ovariectomized estrogen- and progesterone-primed female rats. Intranasal administration of PT-141 (50 $\mu\text{g}/\text{kg}$) induces penile erections in male rats.¹ Formulations containing PT-141 have been used in the treatment of hypoactive sexual desire disorder (HSDD) in premenopausal women.

References

1. Molinoff, P.B., Shadiack, A.M., Earle, D., *et al.* PT-141: A melanocortin agonist for the treatment of sexual dysfunction. *Ann. N.Y. Acad. Sci.* **994**(1), 96-102 (2003).
2. Pfau, J., Giuliano, F., and Gelez, H. Bremelanotide: An overview of preclinical CNS effects on female sexual function. *J. Sex Med.* **4**(Suppl 4), 269-279 (2007).

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, 10/06/2021

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