

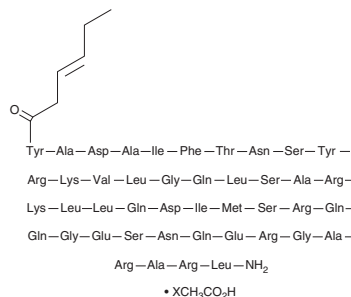
PRODUCT INFORMATION



Tesamorelin (acetate)

Item No. 42498

Formal Name: N-[(3E)-1-oxo-3-hexen-1-yl]-somatoliberin (human pancreatic islet), acetate
Synonyms: TH9507, N-hexanoyl-Tyr1 hGRF (1-44)
Peptide Sequence: N-hexanoyl-YADAIFTNSYRKVLGQLSARKLLQ DIMSRQQGESNQERGARARL-OH
MF: C₂₂₁H₃₆₆N₇₂O₆₇S • XC₂H₄O₂
FW: 5,135.9
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

Tesamorelin (acetate) is supplied as a solid. Aqueous solutions of tesamorelin (acetate) can be prepared by directly dissolving the solid in aqueous buffers. Tesamorelin (acetate) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Tesamorelin is a derivative of human growth hormone-releasing factor (1-44) (hGRF (1-44)) containing a hexanoyl adduct at the N-terminus, rendering it resistant to degradation by dipeptidyl peptidase 4 (DPP-4).¹ *In vivo*, tesamorelin (20 µg/kg twice daily) increases serum levels of insulin-like growth factor 1 (IGF-1) in growing pigs.² It also increases serum levels of growth hormone and IGF-1 in rats and dogs.¹ Formulations containing tesamorelin have been used in the treatment of excess abdominal fat in adults with HIV and lipodystrophy.

References

1. Ferdinandi, E.S., Brazeau, P., High, K., *et al.* Non-clinical pharmacology and safety evaluation of TH9507, a human growth hormone-releasing factor analogue. *Basic Clin. Pharmacol. Toxicol.* **100(1)**, 49-58 (2007).
2. Dubreuil, P. and Désévaux, C. Growth hormone-releasing factor: Structural modification or protection for more potent analogs. *Comb. Chem. High Throughput Screen.* **9(3)**, 171-174 (2006).

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, 02/10/2026

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