

PRODUCT INFORMATION



Eloralintide (acetate)

Item No. 45770

Formal Name: L-γ-glutamyl-S-(mercaptomethyl)-L-cysteinyl-L-asparaginyl-L-threonyl-L-alanyl-L-threonyl-L-cysteinyl-L-alanyl-L-threonylglycyl-L-ornithyl-L-leucyl-L-alanyl-L-α-glutamyl-α-methyl-L-phenylalanyl-L-leucyl-L-valyl-L-arginyl-L-seryl-L-seryl-L-asparaginyl-N-methyl-L-asparaginyl-L-phenylalanylglycyl-L-prolyl-N-[N-(19 carboxy-1-oxononadecyl)-L-γ-glutamyl-L-γ-glutamyl]-L-lysyl-L-leucyl-L-prolyl-L-prolyl-L-threonyl-L-α-glutamyl-L-valylglycyl-L-seryl-L-asparaginyl-L-threonyl-L-tyrosinamide, cyclic (2→7)-thioether, acetate

Peptide Sequence: γE-[CNTATC]ATG-Orn-LAE-α-Me-L-Phe-LVRSSN-N-Me-Asn-FGPK(γE-C20 diacid)-LPPTEVGSNTY-NH₂

MF: C₂₀₁H₃₁₉N₄₉O₆₅S₂ • XC₂H₄O₂

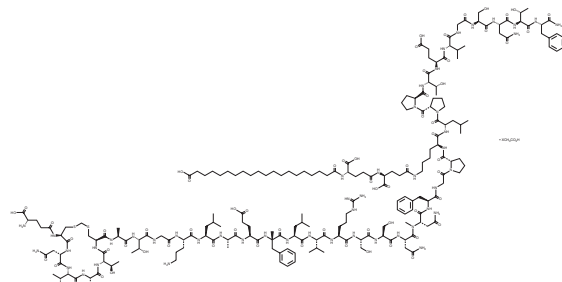
FW: 4,526.2

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

Eloralintide (acetate) is supplied as a solid. A stock solution may be made by dissolving the eloralintide (acetate) in the solvent of choice, which should be purged with an inert gas. Eloralintide (acetate) is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Eloralintide is an agonist of the amylin 1 (AMY₁) receptor.¹ It selectively activates the AMY₁ receptor over the calcitonin and AMY₃ receptors in cAMP activation assays using UM-UC-3 cells expressing the human receptors (EC₅₀s = 23.9, 291, and 253.8 pM, respectively). Eloralintide reduces food intake and induces a loss of up to 12.3% of body weight, primarily through reductions in fat mass, in diet-induced obese (DIO) rats at 3-100 nmol/kg per animal.

Reference

1. Briere, D.A., Qu, H., Lansu, K., *et al.* Eloralintide (LY3841136), a novel amylin receptor agonist for the treatment of obesity: From discovery to clinical proof of concept. *Mol. Metab.* **102**, 102271 (2025).

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, 07/02/2026

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