

PRODUCT INFORMATION

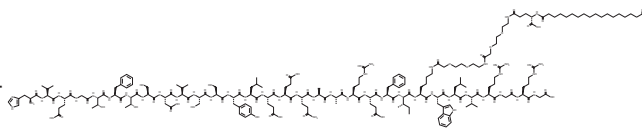


Ecnoglutide

Item No. 41272

CAS Registry No.: 2459531-73-6

Formal Name: L-histidyl-L-valyl-L- α -glutamylglycyl-L-threonyl-L-phenylalanyl-L-threonyl-L-seryl-L- α -aspartyl-L-valyl-L-seryl-L-seryl-L-tyrosyl-L-leucyl-L- α -glutamyl-L- α -glutamyl-L-glutaminy-L-alanyl-L-alanyl-L-arginyl-L- α -glutamyl-L-phenylalanyl-L-isoleucyl-N6-[N-(17-carboxy-1-oxoheptadecyl)-L- γ -glutamyl-2-[2-(2-aminoethoxy)ethoxy]acetyl-2-[2-(2-aminoethoxy)ethoxy]acetyl]-L-lysyl-L-tryptophyl-L-leucyl-L-valyl-L-arginylglycyl-L-arginyl-glycine



Synonym: XW003

Peptide Sequence: HVEGTFTSDVSSYLEEQAAREFIK-(AEEA-AEEA- $\{\gamma$ -Glu)-C18 diacid)-WLVRGRG
(AEEA = 2-(2-(2-aminoethoxy)ethoxy)acetic acid)

MF: C₁₉₄H₃₀₄N₄₈O₆₁

FW: 4,284.8

Purity: \geq 98%

Supplied as: A solid

Storage: -20°C

Stability: \geq 4 years

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

Laboratory Procedures

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

Description

Ecnoglutide is a cAMP-biased peptide glucagon-like peptide 1 receptor (GLP-1R) agonist.¹ It selectively induces cAMP accumulation over β -arrestin recruitment (EC_{50} s = 0.018 and 1,300 nM, respectively) without inducing GLP-1R receptor internalization (EC_{50} = >10,000 nM) in cell-based assays using human GLP-1R. *In vivo*, ecnoglutide (0.15 mg/kg) reduces fasting and post-prandial blood glucose levels in *db/db* mice. It also reduces body weight and food intake in *db/db* mice when administered at a dose of 0.05 mg/kg.

Reference

1. Guo, W., Xu, Z., Zou, H., *et al.* Discovery of ecnoglutide – A novel, long-acting, cAMP-biased glucagon-like peptide-1 (GLP-1) analog. *Mol. Metab.* **75:101762**, (2023).

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, 09/12/2024

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