

PRODUCT INFORMATION



GLP-1 (7-36) Acid (human, mouse, rat, porcine, bovine)

Item No. 41250

Formal Name: L-histidyl-L-alanyl-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 -glutamylglycyl-L-glutamyl-L-alanyl-L-alanyl-L-lysyl-L- α -glutamyl-L-phenylalanyl-L-iso-leucyl-L-alanyl-L-tryptophyl-L-leucyl-L-valyl-L-lysylglycyl-L-arginine, monoacetate

Synonyms: Beinaglutide, Glucagon-like Peptide 1 (7-36) Acid

Peptide Sequence: HAEGTFTSDVSSYLEGQAAKEFIAWLVKGR-OH

MF: $C_{149}H_{225}N_{39}O_{46} \cdot C_2H_4O_2$ • CH_3CO_2H

FW: 3,358.7

Purity: $\geq 98\%$

Supplied as: A solid

Storage: $-20^\circ C$

Stability: ≥ 4 years

H—His—Ala—Glu—Gly—Thr—Phe—Thr—Ser—Asp—Val—
Ser—Ser—Tyr—Leu—Glu—Gly—Gln—Ala—Ala—Lys—
Glu—Phe—Ile—Ala—Trp—Leu—Val—Lys—Gly—Arg—OH

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

Laboratory Procedures

Glucagon-like peptide 1 (GLP-1) (7-36) acid (human, mouse, rat, porcine, bovine) is supplied as a solid. A stock solution may be made by dissolving the GLP-1 (7-36) acid (human, mouse, rat, porcine, bovine) in the solvent of choice, which should be purged with an inert gas. GLP-1 (7-36) acid (human, mouse, rat, porcine, bovine) is sparingly soluble (1-10 mg/ml) in DMSO.

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 GLP-1 (7-36) acid (human, mouse, rat, porcine, bovine) can be prepared by directly dissolving the solid in aqueous buffers. GLP-1 (7-36) acid (human, mouse, rat, porcine, bovine) is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

GLP-1 (7-36) acid is a derivative of the hormone and GLP-1 receptor (GLP-1R) agonist GLP-1 (7-36) amide.^{1,2} It enhances insulin-induced phosphorylation of Akt in 3T3-L1 adipocytes when used at a concentration of 100 nM in the presence of a dipeptidyl peptidase 4 (DPP-4) inhibitor.² GLP-1 (7-36) acid (150 μ g/kg) reduces plasma levels of total cholesterol, triglycerides, free fatty acids, LDL, insulin, and leptin, as well as adipose tissue weight, in a mouse model of obesity induced by a high-fat diet. It decreases blood glucose levels, food intake, and gastric emptying in non-obese and *ob/ob* mice when administered at a dose of 3 mg/kg.¹

References

1. Fang, X., Du, Z., Duan, C., *et al.* Beinaglutide shows significantly beneficial effects in diabetes/obesity-induced nonalcoholic steatohepatitis in *ob/ob* mouse model. *Life Sci.* **270**:118966, (2021).
2. Zhang, F., Chen, Z., Wu, D., *et al.* Recombinant human GLP-1 beinaglutide regulates lipid metabolism of adipose tissues in diet-induced obese mice. *iScience* **24**(12), 103382 (2021).

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, 09/09/2025

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