

PRODUCT INFORMATION



GLP-1 (7-36) amide (trifluoroacetate salt)

Item No. 15069

| | | |
|---------------------|--------------------------------------|---|
| Synonym: | Glucagon-like Peptide 1 (7-36) amide | H—His—Ala—Glu—Gly—Thr—Phe—Thr—Ser—Asp—Val— |
| FW: | 3,297.6 | Ser—Ser—Tyr—Leu—Glu—Gly—Gln—Ala—Ala—Lys— |
| Supplied as: | A crystalline solid | Glu—Phe—Ile—Ala—Trp—Leu—Val—Lys—Gly—Arg—NH ₂ |
| Storage: | -20°C | |
| Stability: | ≥4 years | • XCF ₃ COOH |

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

Laboratory Procedures

GLP-1 (7-36) amide (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the GLP-1 (7-36) amide (trifluoroacetate salt) in water. The solubility of GLP-1 (7-36) amide (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

GLP-1 (7-36) amide is a peptide hormone released from intestinal L-cells upon nutrient consumption. It binds the GLP-1 receptor ($pIC_{50} = 8.5$) in the pancreas and displays various antidiabetic effects by potentiating glucose-induced secretion of insulin from pancreatic β -cells, increasing insulin expression, inhibiting β -cell apoptosis, promoting β -cell neogenesis, reducing glucagon secretion, delaying gastric emptying, promoting satiety, and increasing peripheral glucose disposal.¹

Reference

1. Donnelly, D. The structure and function of the glucagon-like peptide-1 receptor and its ligands. *Br. J. Pharmacol.* **166**(1), 27-41 (2012).

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, 10/24/2022

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