

PRODUCT INFORMATION



GLP-2 (1-34) (human) (trifluoroacetate salt)

Item No. 24756

Formal Name: L-histidyl-L-alanyl-L- α -aspartylglycyl-L-seryl-L-phenylalanyl-L-seryl-L- α -aspartyl-L- α -glutamyl-L-methionyl-L-asparaginyl-L-threonyl-L-isoleucyl-L-leucyl-L- α -aspartyl-L-asparaginyl-L-leucyl-L-alanyl-L-alanyl-L-arginyl-L- α -aspartyl-L-phenylalanyl-L-isoleucyl-L-asparaginyl-L-tryptophyl-L-leucyl-L-isoleucyl-L-glutamyl-L-threonyl-L-lysyl-L-isoleucyl-L-threonyl-L- α -aspartyl-L-arginine, trifluoroacetate salt

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

Synonym: Glucagon-like Peptide 2 (1-34)
MF: C₁₇₁H₂₆₆N₄₈O₅₆S • XCF₃COOH
FW: 3,922.3
Purity: \geq 95%
Supplied as: A lyophilized powder
Storage: -20°C
Stability: \geq 4 years

• XCF₃COOH

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

Laboratory Procedures

GLP-2 (1-34) (human) (trifluoroacetate salt) is supplied as a lyophilized powder. A stock solution may be made by dissolving the GLP-2 (1-34) (human) (trifluoroacetate salt) in water. The solubility of GLP-2 (1-34) (human) (trifluoroacetate salt) in water is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

GLP-2 (1-34) (human) is a synthetic form of the human peptide GLP-2 (Item No. 24414) that contains the full human sequence with the addition of an arginine residue at the C-terminal end.¹ It increases survival of myenteric neurons cultured alone (EC₅₀ = 0.35 nM) and when co-cultured with mast cells to induce cell death.² It does not stimulate insulin release from isolated rat pancreatic islets in the presence of glucose when used at concentrations ranging from 2.5 to 250 nM.¹ GLP-2 (1-34) (human) also does not affect formation of adenomas in a transgenic model of intestinal tumorigenesis using Apc^{Min/+} mice.³

References

- Schmidt, W.E., Siegel, E.G., and Creutzfeldt, W. Glucagon-like peptide-1 but not glucagon-like peptide-2 stimulates insulin release from isolated rat pancreatic islets. *Diabetologia* **28(9)**, 704-707 (1985).
- Voss, U., Sand, E., Hellström, P.M., *et al.* Glucagon-like peptides 1 and 2 and vasoactive intestinal peptide are neuroprotective on cultured and mast cell co-cultured rat myenteric neurons. *BMC Gastroenterol.* **12**, 30 (2012).
- Koehler, J.A., Harper, W., Barnard, M., *et al.* Glucagon-like peptide-2 does not modify the growth or survival of murine or human intestinal tumor cells. *Cancer Res.* **68(19)**, 7897-7904 (2008).

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, 11/17/2022

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