

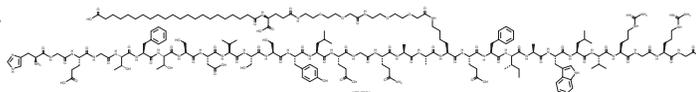
PRODUCT INFORMATION



Bofanglutide (trifluoroacetate salt)

Item No. 44596

Formal Name: 1 [8-glycine,26-[N⁶-[(22S)-22,44-dicarboxy-1,10,19,24-tetraoxo-3,6,12,15-tetraoxa-9,18,23-triazatetracont-1-yl]-L-lysine],34-arginine],7-37-glucagon-like peptide, trifluoroacetate salt



Synonym: GZR18

MF: C₁₈₉H₂₉₅N₄₅O₅₉ • XCF₃COOH

FW: 4,141.7

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

Bofanglutide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the bofanglutide (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Bofanglutide (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

Bofanglutide is a derivative of glucagon-like peptide 1 (GLP-1).¹ It binds to the GLP-1 receptor (GLP-1R) in HEK293 cells expressing the human receptor (IC₅₀ = 3.67 nM) and enhances glucose-stimulated insulin secretion in isolated mouse pancreatic β-cells when used at concentrations of 30 and 90 nM. Bofanglutide decreases non-fasting blood glucose levels in *db/db* mice (ED₅₀ = 3.024 nmol/kg). It reduces food and water intake and lowers serum triglyceride levels in the same mice.

Reference

1. Zhang, M., Zhang, Y., Peng, X., et al. GZR18, a novel long-acting GLP-1 analog, demonstrated positive *in vitro* and *in vivo* pharmacokinetic and pharmacodynamic characteristics in animal models. *Eur. J. Pharmacol.* **928**, 175107 (2022).

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, 02/16/2026

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