

PRODUCT INFORMATION



Bofanglutidum (trifluoroacetate salt)

Item No. 42549

Synonym: HEGTFTSDVSYLEGQAAK-
(AEEA-AEEA-γGlu-C22 diacid)-
EFIAWLVRGRG-OH (AEEA =
2-(2-(2-aminoethoxy)ethoxy)acetic
acid)

MF: C₁₈₉H₂₉₅N₄₅O₅₉ • XCF₃COOH

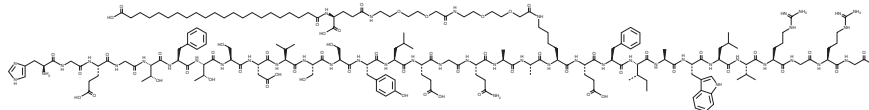
FW: 4,141.7

Purity: ≥95%

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

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

Description

Bofanglutidum is a glucagon-like peptide 1 receptor (GLP-1R) agonist.¹

Reference

1. Gan, Z., Chen, W., Zhang, Y., et al. Long-acting GLP-1 compound. *Gan & Lee Pharmaceuticals Co., Ltd.* PCT/CN2020/141057 (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, 01/27/2025

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