

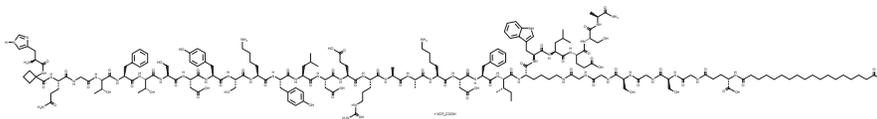
PRODUCT INFORMATION



Survodutide (trifluoroacetate salt)

Item No. 44293

Synonym: BI 456906
MF: $C_{192}H_{289}N_{47}O_{61} \cdot XCF_3COOH$
FW: 4,231.7
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Survodutide (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the survodutide (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Survodutide (trifluoroacetate salt) is soluble (≥ 10 mg/ml) in DMSO.

Description

Survodutide is a peptide agonist of the glucagon receptor (GCGR) and glucagon-like peptide 1 receptor (GLP-1R).¹ It induces cAMP production in CHO-K1 cells expressing human GCGR or GLP-1R (EC_{50} s = 0.52 and 0.33 nM, respectively). *In vivo*, survodutide (30 and 100 nmol/kg) reduces food intake in wild-type but not *Glp1r*^{-/-} mice. It reduces body weight, plasma triglyceride and cholesterol levels, and hepatic alanine transaminase (ALT) and aspartate aminotransferase (AST) levels in a mouse model of diet-induced obesity. It also increases energy expenditure independent of locomotor activity and body weight changes in the same model.

Reference

1. Zimmermann, T., Thomas, L., Baader-Pagler, T., *et al.* BI 456906: Discovery and preclinical pharmacology of a novel GCGR/GLP-1R dual agonist with robust anti-obesity efficacy. *Mol. Metab.* **66:101633**, (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, 10/16/2025

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