

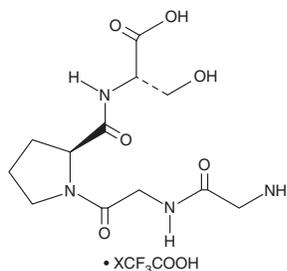
PRODUCT INFORMATION



Exendin-4 (29-32) (trifluoroacetate salt)

Item No. 45083

Formal Name: glycyglycyl-L-prolyl-L-serine, trifluoroacetate salt
Synonym: Gly-Gly-Pro-Ser-OH
Peptide Sequence: GGPS-OH
MF: C₁₂H₂₀N₄O₆ • XCF₃COOH
FW: 316.3
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

Exendin-4 (29-32) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the exendin-4 (29-32) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. Exendin-4 (29-32) (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of exendin-4 (29-32) (trifluoroacetate salt) can be prepared by directly dissolving the solid in aqueous buffers. Exendin-4 (29-32) (trifluoroacetate salt) is sparingly soluble (1-10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Exendin-4 (29-32) is a peptide fragment corresponding to residues 29-32 of exendin-4 (Item No. 11096), a peptide agonist of the glucagon-like peptide 1 receptor (GLP-1R).

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, 02/25/2026

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