

PRODUCT INFORMATION



Exendin-4 (29-39) (trifluoroacetate salt)

Item No. 45084

Formal Name: glycyglycyl-L-prolyl-L-seryl-L-serylglycyl-L-alanyl-L-prolyl-L-prolyl-L-prolyl-L-serine, trifluoroacetate salt

Peptide Sequence: GGPSSGAPPPS-OH

MF: $C_{38}H_{59}N_{11}O_{15} \cdot XCF_3COOH$

FW: 910.0

Purity: $\geq 98\%$

Supplied as: A solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years

H—Gly—Gly—Pro—Ser—Ser—Gly—Ala—Pro—Pro—Pro—Ser—OH
• XCF_3COOH

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

Laboratory Procedures

Exendin-4 (29-39) is supplied as a solid. Aqueous solutions of exendin-4 (29-39) can be prepared by directly dissolving the solid in aqueous buffers. Exendin-4 (29-39) is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Exendin-4 (29-39) is a peptide fragment corresponding to residues 29-39 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, 02/27/2026

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