

PRODUCT INFORMATION

IGF-1 (30-41) (trifluoroacetate salt)

Item No. 44013

Formal Name:	((S)-2-((S)-3-((S)-2-((S)-2-((S)-2-((S)-2-((S)-2-(2-aminoacetamido)-3-(4-hydroxyphenyl)propanamido)acetamido)-3-hydroxypropanamido)-3-hydroxypropanamido)-5-hydroxypropanamido)-5-guanidinopentanamido)-5-guanidinopentanamido)-2-oxobutanoyl)pyrrolidin-1-yl)-L-glutaminy-L-threonine, trifluoroacetate salt	H—Gly—Tyr—Gly—Ser—Ser—Ser—Arg—Arg—Ala—Pro—
Synonyms:	IGF-1 (78-89), IGF-1c, Insulin-like Growth Factor-1 (30-41), Insulin-like Growth Factor-1 (78-89)	Gln—Thr—OH • XCF ₃ COOH
Peptide Sequence:	GYGSSRRAPQT-OH	
MF:	C ₅₁ H ₈₃ N ₁₉ O ₁₉ • XCF ₃ COOH	
FW:	1,266.3	
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

IGF-1 (30-41) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the IGF-1 (30-41) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. IGF-1 (30-41) (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in DMSO.

IGF-1 (30-41) (trifluoroacetate salt) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Insulin-like growth factor 1 (IGF-1) (30-41) is a peptide fragment of the polypeptide hormone IGF-1 (Item No. 33981).¹ It has been used as a starting material in the synthesis of β -sheet mimics of IGF-1.

Reference

1. Shang, Y., Zhi, D., Feng, G., *et al.* Supramolecular nanofibers with superior bioactivity to insulin-like growth factor-I. *Nano Lett.* **19**(3), 1560-1569 (2019).

WARNING

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, 11/03/2025

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