

PRODUCT INFORMATION



IGF-2 (33-40) (trifluoroacetate salt)

Item No. 44047

Formal Name: (L-seryl-L-arginyl-L-valyl-L-seryl-L-arginyl-L-arginyl-L-seryl-L-arginine), trifluoroacetate salt

Synonyms: H-Ser-Arg-Val-Ser-Arg-Arg-Ser-Arg-OH, Insulin-like Growth Factor 2

Peptide Sequence: H-SRVSRRSR-OH

MF: C₃₈H₇₄N₂₀O₁₂ • XCF₃COOH

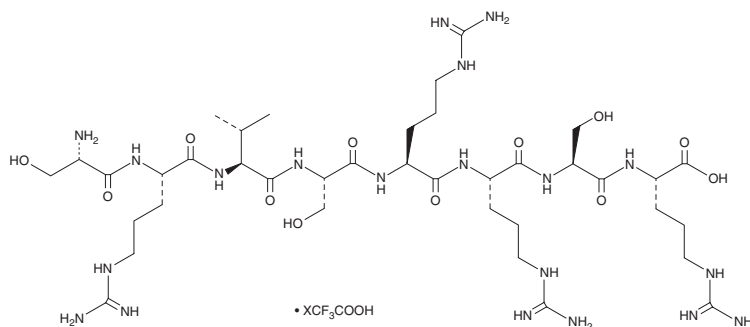
FW: 1,003.1

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

IGF-2 (33-40) (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

IGF-2 (33-40) is a peptide fragment of insulin-like growth factor 2 (IGF-2).¹ It has been used in the development of antibodies against IGF-2.²

References

1. Blyth, A.J., Kirk, N.S., and Forbes, B.E. Understanding IGF-II action through insights into receptor binding and activation. *Cells* **9(10)**, 2276 (2020).
2. Meyer-Schwickerath, R., Pfeiffer, A., Blum, W.F., et al. Vitreous levels of the insulin-like growth factors I and II, and the insulin-like growth factor binding proteins 2 and 3, increase in neovascular eye disease. Studies in nondiabetic and diabetic subjects. *J. Clin. Invest.* **92(6)**, 2620-2625 (1993).

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, 12/17/2025

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