

PRODUCT INFORMATION



IGRP (206-214) (mouse) (trifluoroacetate salt)

Item No. 44049

Formal Name: L-valyl-L-tyrosyl-L-leucyl-L-lysyl-L-threonyl-L-asparaginyl-L-valyl-L-phenylalanyl-L-leucine, trifluoroacetate salt

Synonyms: Islet-specific Glucose-6-phosphatase Catalytic Subunit-related Protein (206-214), Val-Tyr-Leu-Lys-Thr-Asn-Val-Phe-Leu-OH

Peptide Sequence: VYLKTNVFL-OH

MF: C₅₄H₈₅N₁₁O₁₃ • XCF₃COOH

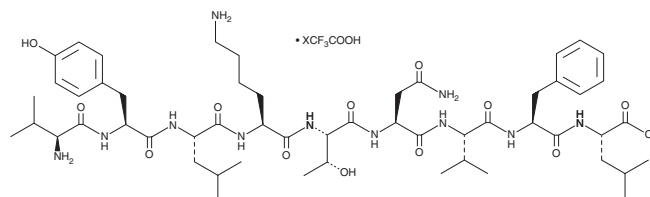
FW: 1,096.3

Purity: ≥95%

Supplied as: A crystalline 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

IGRP (206-214) (mouse) (trifluoroacetate salt) is supplied as a crystalline solid. A stock solution may be made by dissolving the IGRP (206-214) (mouse) (trifluoroacetate salt) in the solvent of choice, which should be purged with an inert gas. IGRP (206-214) (mouse) (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) in DMSO.

IGRP (206-214) (mouse) (trifluoroacetate salt) is slightly soluble (0.1-1 mg/ml) 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

Islet-specific glucose-6-phosphatase catalytic subunit-related protein (IGRP) (206-214) is an autoantigenic peptide for β -cells.¹ It promotes mouse T cell-induced lysis of RMA-S cells expressing the MHC class I receptor in a concentration-dependent manner. IGRP (206-214) is bound in a tetramer assay by CD8⁺ T cells isolated from pancreatic islets but not those isolated from peripheral blood in non-obese diabetic mice.

Reference

1. Lieberman, S.M., Evans, A.M., Han, B., *et al.* Identification of the β cell antigen targeted by a prevalent population of pathogenic CD8⁺ T cells in autoimmune diabetes. *Proc. Natl. Acad. Sci. USA* **100**(14), 8384-8388 (2003).

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, 09/30/2025

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