

PRODUCT INFORMATION



Glucose-dependent Insulinotropic Polypeptide (3-42) (human) (trifluoroacetate salt)

Item No. 36908

Formal Name: L- α -glutamylglycyl-L-threonyl-L-phenylalanyl-L-isoleucyl-L-seryl-L- α -aspartyl-L-tyrosyl-L-seryl-L-isoleucyl-L-alanyl-L-methionyl-L- α -aspartyl-L-lysyl-L-isoleucyl-L-histidyl-L-glutamyl-L-glutamyl-L- α -aspartyl-L-phenylalanyl-L-valyl-L-asparaginyl-L-tryptophyl-L-leucyl-L-leucyl-L-alanyl-L-glutamyl-L-lysylglycyl-L-lysyl-L-lysyl-L-asparaginyl-L- α -aspartyl-L-tryptophyl-L-lysyl-L-histidyl-L-asparaginyl-L-isoleucyl-L-threonyl-L-glutamine, trifluoroacetate salt

H — Glu — Gly — Thr — Phe — Ile — Ser — Asp — Tyr — Ser — Ile —
Ala — Met — Asp — Lys — Ile — His — Gln — Gln — Asp — Phe —
Val — Asn — Trp — Leu — Leu — Ala — Gln — Lys — Gly — Lys —
Lys — Asn — Asp — Trp — Lys — His — Asn — Ile — Thr — Gln — OH
• XCF₃COOH

Synonyms: Gastric Inhibitory Peptide 1 (3-42), GIP-1 (3-42)

Peptide Sequence: EGTFFISDYSIAMDKIHQQDFVNWLLAQKGGKNDWKHNITQ-OH

MF: C₂₁₄H₃₂₄N₅₈O₆₃S • XCF₃COOH

FW: 4,749.3

Purity: ≥90%

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

Gastric inhibitory peptide 1 (GIP-1) (3-42) (human) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the GIP-1 (3-42) (human) (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

Description

GIP-1 (3-42) is a peptide fragment of the incretin hormone GIP (Item No. 25004) and a GIP receptor antagonist.¹ It is formed from GIP by serum dipeptidyl peptidase 4 (DDP-4). GIP-1 (3-42) (100 nM) reduces insulin secretion from BRIN-BD11 pancreatic cells. It increases plasma glucose levels and decreases plasma insulin levels in an *ob/ob* mouse model of diabetes when administered at a dose of 25 nmol/kg.

Reference

1. Gault, V.A., Parker, J.C., Harriott, P., *et al.* Evidence that the major degradation product of glucose-dependent insulinotropic polypeptide, GIP(3-42), is a GIP receptor antagonist *in vivo*. *J. Endocrinol.* **175**(2), 525-533 (2002).

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, 05/09/2024

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