

# PRODUCT INFORMATION



## Glucose-dependent Insulinotropic Polypeptide (1-39) (porcine) (trifluoroacetate salt)

Item No. 36865

**Formal Name:** L-tyrosyl-L-alanyl-L- $\alpha$ -glutamylglycyl-L-threonyl-L-phenylalanyl-L-isoleucyl-L-seryl-L- $\alpha$ -aspartyl-L-tyrosyl-L-seryl-L-isoleucyl-L-alanyl-L-methionyl-L- $\alpha$ -aspartyl-L-lysyl-L-isoleucyl-L-arginyl-L-glutaminy-L-glutaminy-L- $\alpha$ -aspartyl-L-phenylalanyl-L-valyl-L-asparaginy-L-tryptophyl-L-leucyl-L-leucyl-L-alanyl-L-glutaminy-L-lysylglycyl-L-lysyl-L-lysyl-L-seryl-L- $\alpha$ -aspartyl-L-tryptophyl-L-lysyl-L-histidyl-L-asparagine, trifluoroacetate salt

**Synonyms:** Gastric Inhibitory Peptide (1-39), GIP (1-39)

**Peptide Sequence:** YAEGTFISDYSIAMDKIRQQDFVNWLLAQKGGKSDW  
KHN-OH

**MF:**  $C_{210}H_{316}N_{56}O_{61}S \cdot XCF_3COOH$

**FW:** 4,633.2

**Purity:**  $\geq 98\%$

**Supplied as:** A solid

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

**Stability:**  $\geq 4$  years

H-Tyr-Ala-Glu-Gly-Thr-Phe-Ile-Ser-Asp-Tyr-  
Ser-Ile-Ala-Met-Asp-Lys-Ile-Arg-Gln-Gln-  
Asp-Phe-Val-Asn-Trp-Leu-Leu-Ala-Gln-Lys-  
Gly-Lys-Lys-Ser-Asp-Trp-Lys-His-Asn-OH  
•  $XCF_3COOH$

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

### Laboratory Procedures

Glucose-dependent Insulinotropic Polypeptide (GIP) (1-39) (porcine) (trifluoroacetate salt) is supplied as a solid. A stock solution may be made by dissolving the GIP (1-39) (porcine) (trifluoroacetate salt) in water. We do not recommend storing the aqueous solution for more than one day.

### Description

GIP (1-39) is an endogenous 39-amino acid peptide hormone that induces insulin secretion.<sup>1</sup> It has been found in the upper intestines of pigs. GIP (1-39) (100 nM) enhances glucose-induced increases in insulin release in primary rat pancreatic  $\beta$ -cells to a greater extent than GIP (1-42).

### Reference

1. Xie, L., Lu, J., Östenson, C.-G., *et al.* GIP<sub>1-39</sub>, a novel insulinotropic peptide form and aspects on its mechanism of action. *Regul. Pept.* **121(1-3)**, 107-112 (2004).

#### 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