Gastric inhibitory peptide (3-42) (human) monoclonal antibody (25)

Human gastric inhibitory polypeptide (GIP) is a 42 amino acid peptide belonging to the glucagon-secretin family of peptide hormones. It is secreted by endocrine cells in the duodenal mucosa and stimulates glucose-dependent insulin secretion as well as GLP-1 release from more distal endocrine (L) cells in the intestinal mucosa. GIP shows amino-acid sequence similarities to glucagon, GLP-1 and GLP-2 (from approximately 50% identity for glucagon to 30% identity for GLP-2).

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BPD-ABS-057-25-1	1mg
BPD-ABS-057-25-04	400μg

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name GIP

Application ELISA

Clone 25

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M

sodium chloride and 15mM sodium azide.

GenBank ID 2695

Host Mouse

Immunogen Synthetic human gastric inhibitory polypeptide

Purity Detail Protein A purified.

Species Reactivity Human

Specificity The epitope lies within the N-terminal region of GIP (3-42)

P09681 **UniProt ID**

This antibody is covered by our Worry-Free Guarantee **Worry-free Guarantee**

Last modified: May 29, 2024

