

# Gastric inhibitory peptide (human) monoclonal antibody (32)

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-054-26-1	1mg
BPD-ABS-054-26-04	400µg

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Short Term Storage** +4°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** GIP

**Application** ELISA

**Clone** 32

**Formulation** Liquid. In 0.01 M phosphate buffer, pH 7.4, containing 0.5 M NaCl and 15 mM sodium azide.

**Host** Mouse

**Immunogen** Synthetic human gastric inhibitory polypeptide (1-42)

**Purity Detail** Protein A purified.

**Species Reactivity** Human

**Worry-free Guarantee** This antibody is covered by our [Worry-Free Guarantee](#).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)