

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](#).

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

Last modified: May 29, 2024



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)