

Gastrin I (human), ELISA kit

The Gastrin I (human), ELISA kit is a colorimetric competitive enzyme immunoassay kit

The Gastrin I (human), EIA kit is a colorimetric competitive enzyme immunoassay kit with results in 5 hours. Absorbance is read at 405 nm. Ready-to-use liquid color-coded reagents are provided to reduce error in the lab.

Gastrins are a family of sequence-related carboxyamidated peptides produced by endocrine G cells of the antrum mucosa, and are potent stimulators of gastric acid secretion. Antral distension, partially digested proteins, amino acids, and vagal stimulation resulting from smelling, tasting, chewing or swallowing food all contribute to gastrin release from G cell storage. In addition, caffeine, alcohol, hypoglycemia, antacids and elevated calcium levels also stimulate gastrin release. Increased serum gastrin levels are associated with duodenal ulcers, Helicobacter pylori infections, colorectal carcinomas, and other tumors and cancerous lesions.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

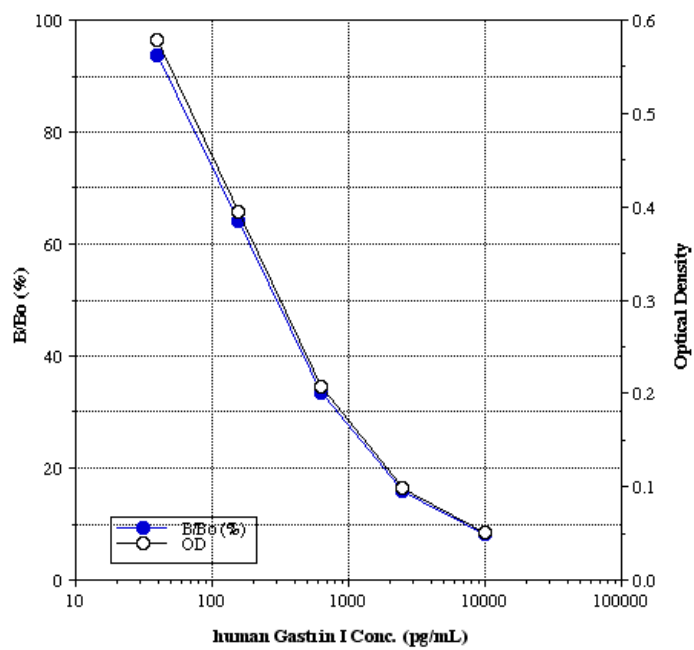
ADI-900-026

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Safe - non-radioactive reagents
- Ready-to-use - color coded reagents and pre-coated plates reduce errors
- highly sensitive
- Measure up to 39 samples in duplicate in 5 hours



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	Colorimetric detection, ELISA
Application Notes	For the quantitative determination of human Gastrin I in culture supernatants, plasma, and serum.
Assay Time	5 hours
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Contents	GxR IgG Microtiter plate, Conjugate, Antibody, Assay buffer, Wash buffer concentrate, Standard, pNpp Substrate, Stop solution
Crossreactivity	Gastrin I (1-17) (100%), Minigastrin (G13-I) (74.6%), rat Gastrin I (70.7%), Gastrin II (G-17-II, sulfated) (9.30%), Cholecystokinin 26-33 (CCK-8) (2.67%), Gastrin Tetrapeptide (CCK-4) (1.60%), Big Gastrin (G34-I) (0.80%) and <0.001%: Gastrin Releasing Peptide (GRP), Gastrin Inhibitory Polypeptide (GIP), Glucagon, Bombesin, Pancreatic Polypeptide, Vasoactive Intestinal Peptide (VIP), Somatostatin-14
Sensitivity	7.27 pg/ml (range 39.1 – 10,000 pg/ml)
Species Reactivity	Human
UniProt ID	P01350
Wavelength	405 nm



Last modified: May 29, 2024

ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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