

# Gastrin I (rat), ELISA kit

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

The Gastrin I (rat), 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. This kit utilizes low sample volume and the ability to measure up to 37 samples in duplicate per 96-well plate.

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 »](#)

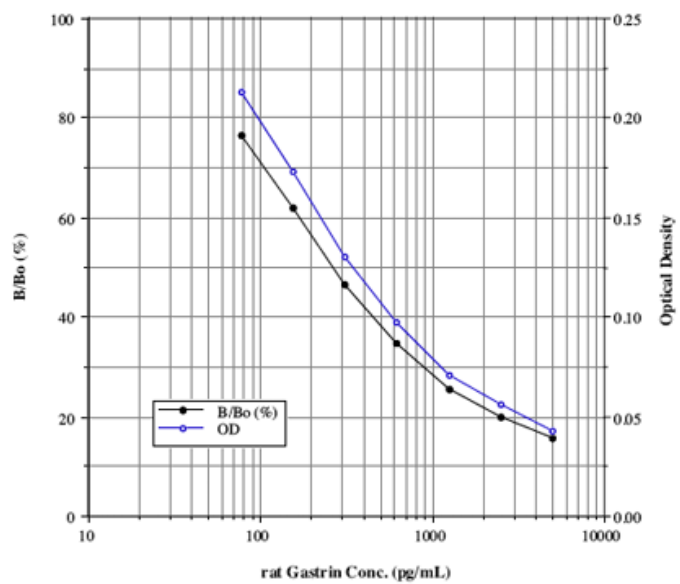
ADI-900-149

96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Efficient - only 50µl sample volume required
- Ultrasensitive - detect <80 pg/ml
- Higher throughput - assay up to 37 samples in duplicate per 96 well kit



## 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 rat 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 Rat Gastrin I (100%), Cholecystokinin 26-33 (CCK-8) (100%), human Gastrin I (30%), mouse Gastrin I (20.5%), Minigastrin (G13-I) (11%), Gastrin Tetrapeptide (CCK-4) (0.75%), Glucagon (0.5%) and <0.2%: Gastrin Inhibitory Polypeptide (GIP), Bombesin, Somatostatin-14, Gastrin Releasing Peptide (GRP)

Sensitivity 78.1 pg/ml (range 78.1 – 5,000 pg/ml)

Species Reactivity Rat

UniProt ID P04563

Wavelength 405 nm



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-  
fr@enzolifesciences.com](mailto:info-fr@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)