

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 »](#)

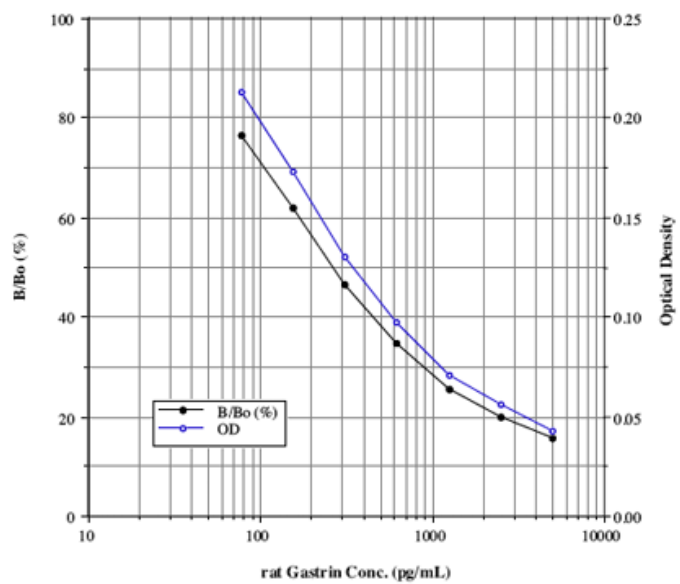
ADI-900-149

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

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