

Peptide YY (human) monoclonal antibody (1) (biotin conjugate)

Peptide tyrosine-tyrosine (PYY) amide is a 36 amino-acid. PYY is a member of the pancreatic polypeptide (PP) family of peptides (neuropeptide Y (NPY), PYY and PP). While NPY is localized in neurons (both in the central and peripheral nervous system) and PP in the pancreatic islets, PYY is found in both neurons and the gut. PYY is released in response to food intake from the same endocrine cells (L cells) in the intestinal mucosa as the glucagon-like peptides, and inhibits gall bladder secretion, gut motility and pancreatic secretion. These effects are similar to those of PP and overlap with the gut inhibitory activity of GLP-1.

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

Ordering Information

[Order Online »](#)

BPD-ABS-029-01B-015	150µg
---------------------	-------

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	PYY, peptide tyrosine-tyrosine amide
Application	ELISA
Application Notes	Detects a band of ~4kDa by Western blot.
Clone	1
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.14M sodium chloride and 15mM sodium azide.
Host	Mouse
Immunogen	Synthetic human Peptide YY.
Isotype	IgG1κ
Purity Detail	Protein A/G-affinity purified.
Species Reactivity	Human
UniProt ID	P10082
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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)