Peptide YY (3-36) monoclonal antibody (4)

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-051-04-1	1mg
BPD-ABS-051-04-04	400µg

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, WB

Application NotesDetects a band of ~4kDa by Western blot.

Clone 4

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M

sodium chloride and 15mM sodium azide.

Host Mouse

Immunogen Synthetic peptide corresponding to aa 3-36 of human

Peptide YY coupled to carrier.

Isotype IgG2ak

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 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

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