

Pancreatic polypeptide (human) monoclonal antibody (6)

Pancreatic polypeptide (PP) is a 36 amino-acid residue peptide, with sequence and 3-D structural homology with NPY and PYY. PP is the least conserved of the three peptides and the first to be identified. It is almost exclusively expressed in the endocrine pancreas and released in response to cholinergic stimulation and food intake. PP has inhibitory effects on gall bladder secretion, intestinal motility and pancreatic secretion. PP may serve as a marker for certain gut endocrine tumors including carcinoid tumors.

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

Ordering Information

[Order Online »](#)

BPD-ABS-030-06-1	1mg
BPD-ABS-030-06-04	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	PP
Application	ELISA
Application Notes	Detects a band of ~4kDa by Western blot.
Clone	6
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.
Host	Mouse
Immunogen	Synthetic human Pancreatic Polypeptide.
Isotype	IgG1κ
Purity Detail	Protein A/G-affinity purified.
Species Reactivity	Human
UniProt ID	P01298
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-
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)