

Vasoactive intestinal peptide monoclonal antibody (2)

Vasoactive intestinal peptide (VIP) is a 28-amino-acid peptide belonging to the secretin-glucagon family of peptide hormones. Its sequence is identical in man, ox, pig and rat. VIP is widely distributed in the central and peripheral nervous systems, acting as a neurotransmitter or neuromodulator. Peripherally it is involved in vaso- and bronchodilatation, smooth muscle relaxation and stimulation of secretion. Ectopic endocrine tumors secreting VIP (VIPomas) lead to raised plasma levels of VIP and are associated with watery diarrhea, hypokalemia and achlorhydria syndrome. VIP shows high sequence homology with peptide histidine-methionine (PHM) in man or peptide histidine-serine (PHI) in pigs or rats, present in the same precursor, and with the neuropeptide pituitary adenylate cyclase activating polypeptide (PACAP).

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

Ordering Information

[Order Online »](#)

BPD-ABS-023-02-1	1mg
BPD-ABS-023-02-04	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 VIP

Application ELISA, WB

Application Notes Detects a band of ~3kDa by Western blot.

Clone 2

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.

Host Mouse

Immunogen Vasoactive intestinal peptide.

Isotype IgG1κ

Purity Detail Protein A/G-affinity purified.

Species Reactivity Bovine, Human, Porcine, Rat

UniProt ID P01282

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



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