

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-023-02-04	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

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-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)