

Peptide Histidine-Methionine (human) monoclonal antibody (3)

Human peptide histidine-methionine (PHM) is a 27-amino-acid peptide belonging to the glucagon-secretin family of peptide hormones. PHM is produced from the same prohormone as VIP, with which it shows considerable sequence homology. The VIP/PHM prohormone is produced by neurons in the central and peripheral nervous systems, often together with acetylcholine. In some situations, the peptides reinforce and prolong the effects of neurotransmitter release. Like VIP, PHM is a vasodilator and stimulator of intestinal fluid secretion. High plasma levels of PHM are diagnostic of the corresponding peptide-producing tumor (VIPoma), associated with the watery diarrhea-hypokalemia-achlorhydria syndrome.

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

Ordering Information

[Order Online »](#)

BPD-ABS-020-03-1	1mg
BPD-ABS-020-03-04	400µg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	PHM
Application	ELISA, IHC
Clone	3
Crossreactivity	Cross-reacts with PACAP (approximately 12%; incomplete, non-parallel inhibition). ABS 020-03 does not cross-react with GLP-1, GLP-2, VIP or glucagon.
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.
GenBank ID	7432
Host	Mouse
Immunogen	Synthetic human peptide histidine-methionine.
Isotype	IgG2a
Purity Detail	Protein A/G-purified.
Source	From culture supernatant.
Species Reactivity	Human
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)