

Chromogranin A (human) monoclonal antibody (PHE5)

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

Citations: 3

[View Online »](#)

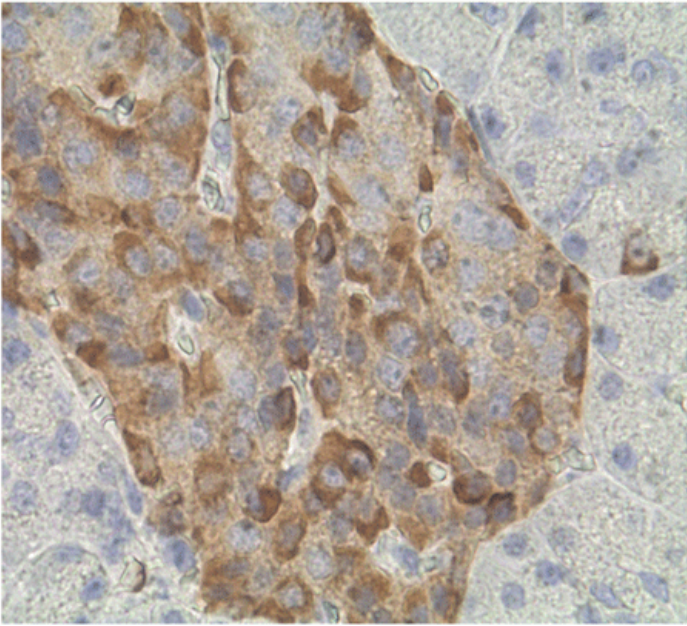
Ordering Information

[Order Online »](#)

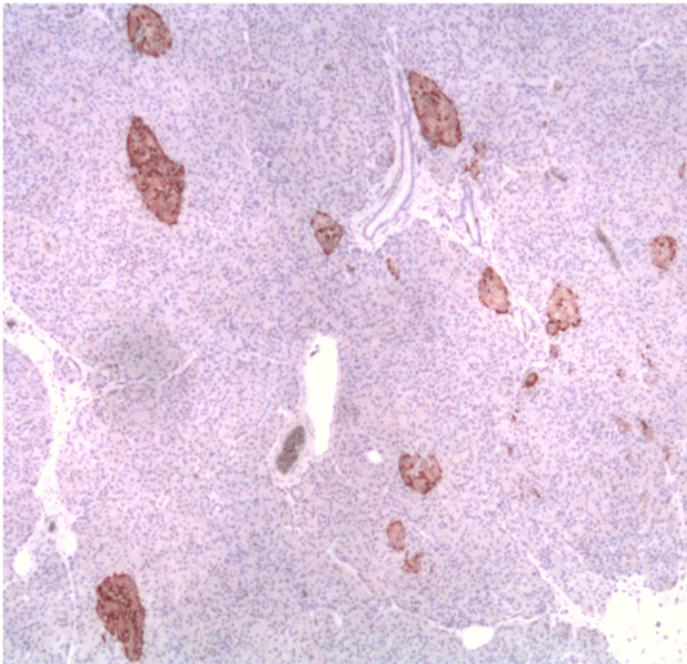
ENZ-C34932	~10 µg - 0.5ml
------------	----------------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of human endocrine pancreas tissue stained with chromogranin A, mAb (PHE5). Original magnifications: 5x and 63x, respectively.



Immunohistochemistry analysis of human endocrine pancreas tissue stained with chromogranin A, mAb (PHE5). Original magnifications: 5x and 63x, respectively.

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Neuroendocrine, CgA
Application	IHC (PS)
Clone	PHE5
Formulation	Liquid. In PBS containing 0.1% sodium azide and 1% BSA.
Host	Mouse
Immunogen	Human pheochromocytoma.
Isotype	IgG1
Purity Detail	Affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (paraffin sections; this antibody can be used for staining of paraffin-embedded tissue sections fixed in neutral buffered formalin and is compatible with most commonly used histological fixatives. Choose a good quality enzyme-linked anti-mouse IgG1 antibody for detection). A 20- to 50-fold dilution of the antibody is suggested. Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually.
Source	Purified from hybridoma tissue culture supernatant.
Species Reactivity	Human

Specificity

PHE5 is specific to chromogranin A, a highly acidic protein associated with the “dense core” granules of many cells of the neural and neuroendocrine system. It reacts with neuroendocrine cells of the gut, pancreatic islet and adrenal medulla, but does not react with any cells in the brain or with melanoma cells. The antibody stains carcinoid tumors, pheochromocytomas and paragangliomas. It is nonreactive with carcinomas, lymphomas and sarcomas.

UniProt ID

P10645

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)