

Chromogranin A (human) monoclonal antibody (PHE5) (Ready-to-Use)

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

Citations: 3

[View Online »](#)

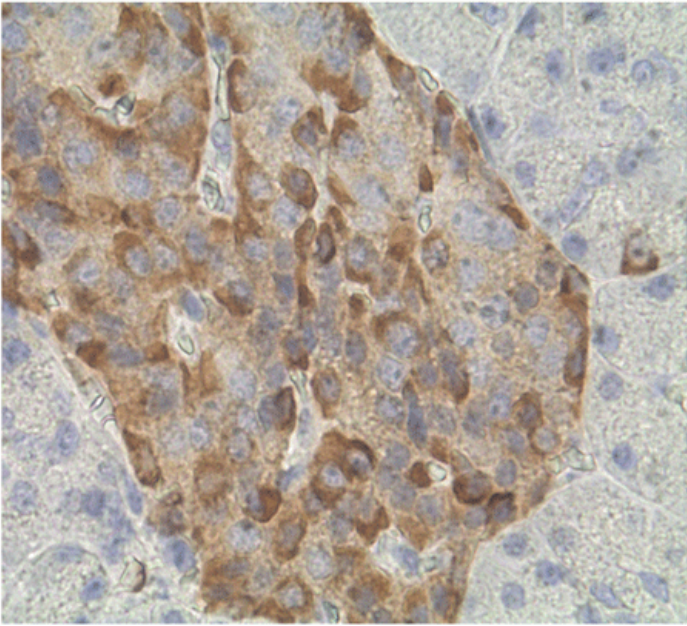
Ordering Information

[Order Online »](#)

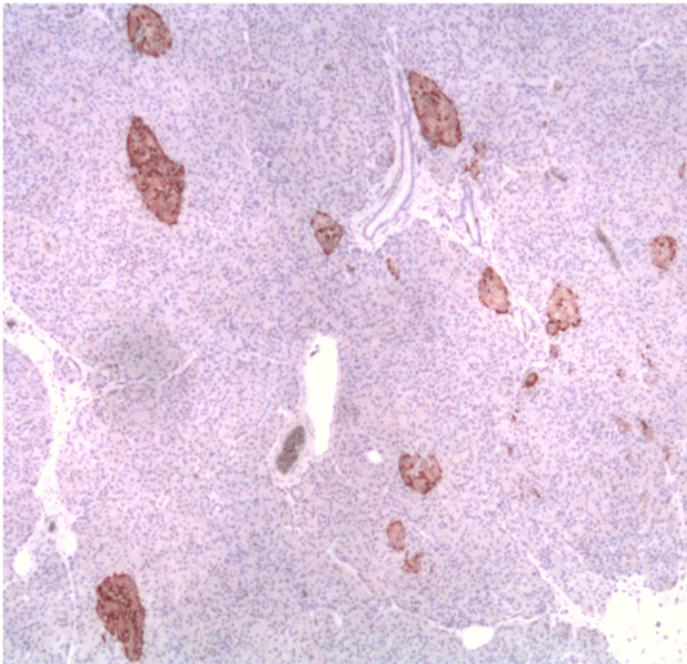
| | |
|-----------|-------------|
| ENZ-30932 | ~6 µg - 6ml |
|-----------|-------------|

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. |
| Long Term Storage | +4°C |
| Shipping | Blue Ice |

Regulatory Status RUO - Research Use Only

Product Details

| | |
|-------------------------------------|--|
| Alternative Name | Neuroendocrine, CgA |
| Application | IHC (PS) |
| Clone | PHE5 |
| Formulation | Liquid. <i>Ready-to-use</i> solution 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). 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](#)

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)