Chromogranin A (human) monoclonal antibody (PHE5)

This antibody is covered by our Worry-Free Guarantee.

Citations: 3

View Online »

Ordering Information

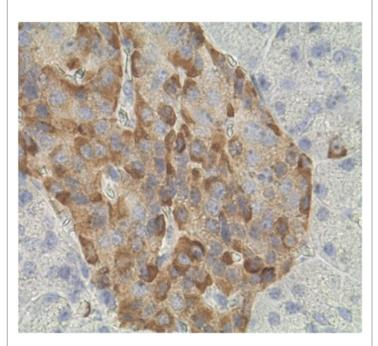
Order Online »

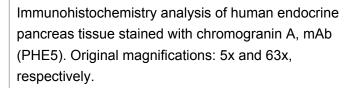
ENZ-C34932

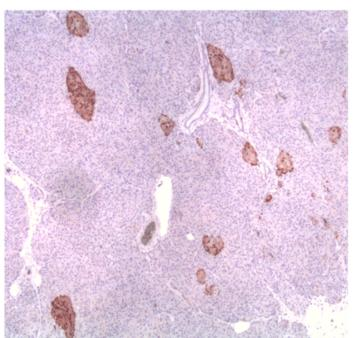
~10 µg - 0.5ml

Manuals, SDS & CofA

View Online »







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

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.

lgG1

Purity Detail Affinity purified.

Recommendation Dilutions/Conditions Immunohistochemistry (paraffin sections; this antibody can

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

be used for staining of paraffin-embedded tissue sections

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



info-