

Chromogranin A polyclonal antibody

Chromogranin A (CgA) is a prohormone which is proteolytically processed in neurons prior to axonal transport. The differential expression of chromogranin A fragments in neuronal populations suggests distinct functional roles for each CgA-derived peptide. Immunohistochemical localization of CgA has been used as a marker for the identification and prognostic assessment of neuroendocrine tumours and CgA has also been reported to be a component of senile plaques in the brains of demented individuals.

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

Ordering Information

[Order Online »](#)

ENZ-ABS268-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CgA, CHGA, Parathyroid secretory protein 1

Application ELISA, IHC

Crossreactivity Expected reactivity due to sequence homology with bovine, chimpanzee, dog, horse, pig, mouse, rat.

Formulation Liquid. In PBS containing 0.09% sodium azide.

Host Sheep

Immunogen Synthetic peptide corresponding to aa 1-15 (L1PVNSPMNKGDTVM15) of native human chromogranin A, conjugated to KLH.

Species Reactivity Human

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)