

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](#).

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com