

PRODUCT INFORMATION

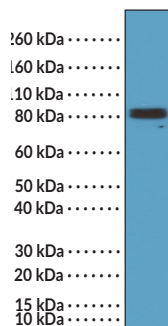
Chromogranin A (N-Term) Rabbit Monoclonal Antibody (Clone RM385)

Item No. 32312

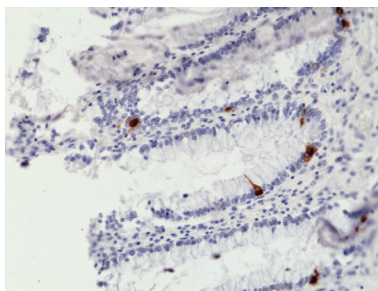
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	CGA, Parathyroid Secretory Protein 1, Pituitary Secretory Protein 1
Immunogen:	Peptide from the N-terminal region of human chromogranin A
Cross Reactivity:	(+) Chromogranin A
Species Reactivity:	(+) Human
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM385
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:1,000-1:10,000 for IHC and 1:1,000-1:2,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

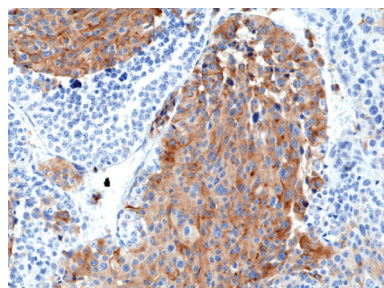
Images



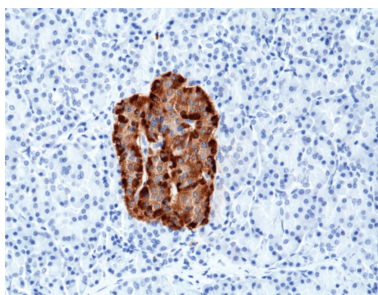
WB of human brain tissue lysate using Chromogranin A (N-Term) Rabbit Monoclonal Antibody (Clone RM385) at a dilution of 1:1,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human colon tissue using Chromogranin A (N-Term) Rabbit Monoclonal Antibody (Clone RM385) at a dilution of 1:10,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded medullary thyroid carcinoma tissue using Chromogranin A (N-Term) Rabbit Monoclonal Antibody (Clone RM385) at a dilution of 1:2,500.



Immunohistochemical staining of formalin-fixed and paraffin-embedded normal human pancreas tissue using Chromogranin A (N-Term) Rabbit Monoclonal Antibody (Clone RM385) at a dilution of 1:2,500.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 02/21/2024

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Chromogranin A is a glycoprotein and member of the chromogranin/secretogranin family of neuroendocrine secretory proteins that is encoded by *CHGA* in humans.¹⁻³ It is expressed in endocrine, neuroendocrine, and immune cells and localized to secretory granules. Chromogranin A is composed of multiple peptide domains and is proteolytically cleaved by prohormone convertases, plasmin, kallikrein, and cathepsin L to generate various bioactive peptides, including paeceastatin, vasostatin-1, catestatin, and serpinin, which have roles in insulin sensitivity, vasodilation, inflammation, and cardiac contractility, respectively. Plasma levels of chromogranin A are increased in patients with hypertension, inflammatory bowel disease, rheumatoid arthritis, and severe sepsis, as well as carcinoid, pancreatic islet, or thyroid tumors.³ Cayman's Chromogranin A (N-Term) Rabbit Monoclonal Antibody (Clone RM385) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Di Giacinto, P., Rota, F., Rizza, L., *et al.* Chromogranin A: From laboratory to clinical aspects of patients with neuroendocrine tumors. *Int. J. Endocrinol.* 8126087 (2018).
2. Eissa, N., Hussein, H., Hendy, G.N., *et al.* Chromogranin-A and its derived peptides and their pharmacological effects during intestinal inflammation. *Biochem. Pharmacol.* **152**, 315-326 (2018).
3. Mahata, S.K. and Corti, A. Chromogranin A and its fragments in cardiovascular, immunometabolic, and cancer regulation. *Ann. N. Y. Acad. Sci.* **1455(1)**, 34-58 (2019).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM