

A new version of CISBIO Chromogranin A ELISA kit (#: CGA-ELISA-NG)

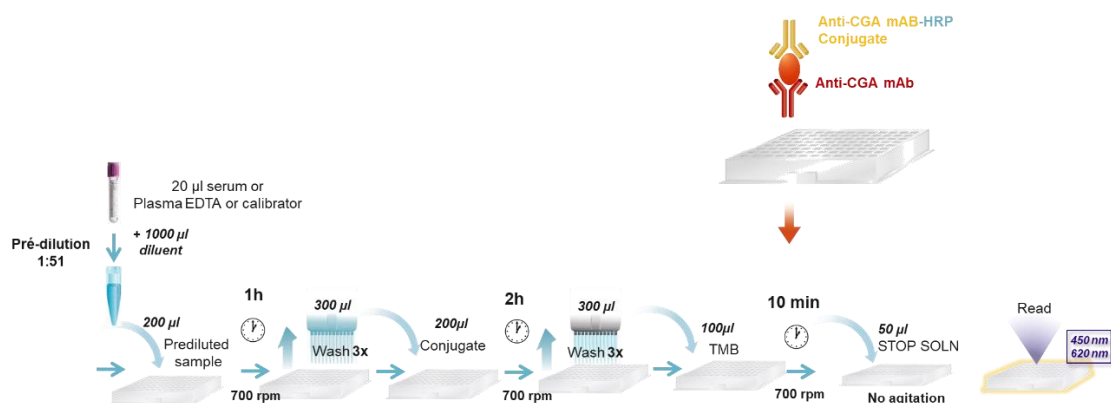
Introduction

Pioneered in the early 2000s by CISBIO, the chromogranin A assay (CGA) using RIA technology has provided a marker of choice for neuroendocrine tumors (NETs). In 2008, to follow the evolution of laboratory assays, a CGA ELISA kit was added to our diagnostic tools.

In 2022, a new version of this kit is available (CGA-ELISA-NG) and further improves the quality of the assays for this tumor marker. This version is now fully compatible with the new European regulatory rules of the IVDR 2017/746.

All other parameters have been optimized to best meet the requirements. Incubation times have been reduced, dilution and recovery tests have been improved, the robustness of the kit has been optimized. The following paragraphs support these improvements.

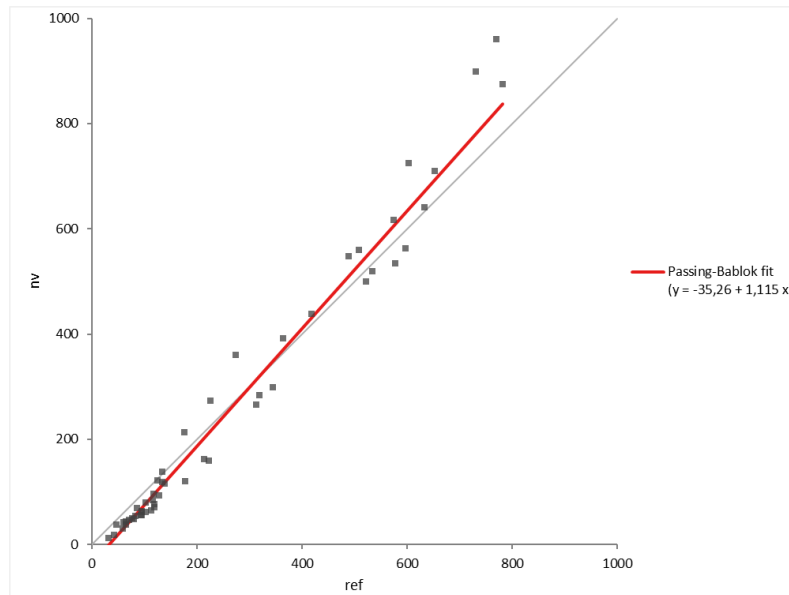
The new kit uses the same antibodies as the previous ones to maintain the same specificity. The antibodies are now produced in cell culture, thus reducing the use of animals.



Correlation between CGA-RIA-CT and CGA-ELISA-NG:

The experiment was performed on two batches of CGA-ELISA-NG and 2 batches of CGA-RIA-CT on 109 serum samples covering the whole calibration range.

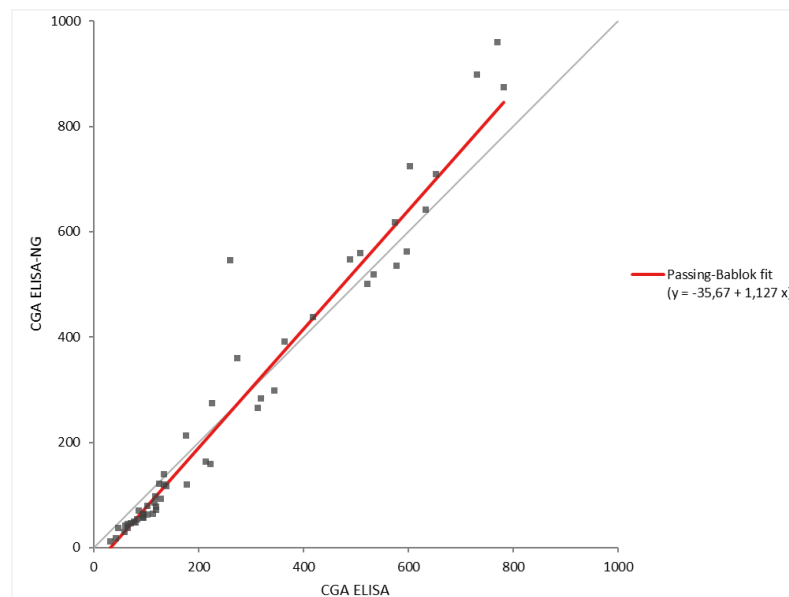
The results below show a good correlation between the kits with a slope close to 1.0 which will facilitate the switch from previous ELISA version to the CGA-ELISA-NG version.



Correlation - r | 0,988

Correlation between CGA-ELISA and CGA-ELISA-NG:

Similar correlation was performed using 53 samples covering the whole range of concentrations. The little gap on the slope is linked to the adjustment of the new kit to obtain optimal dilution performances along the calibration curve. Nevertheless, the correlation is good.



Correlation - r | 0,977

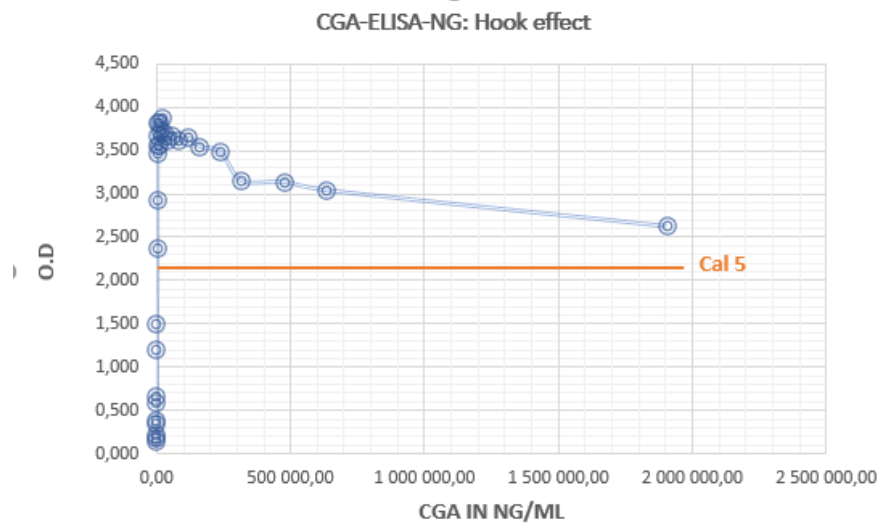
Essential analytical characteristics of the CGA-ELISA-NG kit:

Hook effect:

In this experiment, CGA antigen is diluted in normal human serum to cover a range of CGA concentration from 2 million ng/mL to 0 ng/mL.

We can see on the graph below that up to 2 million ng/mL, no hook effect appears making this kit perfectly compatible with very high serum concentrations of CGA avoiding problems of false

negatives. In the opinion of clinicians, the highest values that can be encountered are below one million ng/mL (L. Chardon HCL Lyon 2021 internal meeting).



Recovery and Dilution:

To control the similarity of the assay response between exogenous CGA and serum CGA, recovery tests were performed. The results show a perfect similarity of the assays whatever the origin of the antigen hence meeting Cisbio internal standards in force (90%-110%).

	Recoveries		
	[CGA] ng /mL	% Recovery Min	% Recovery Max
Sample 1	200	98,1%	105,2%
Sample 2	401	100,8%	106,2%
Sample 3	770	97,5%	100,1%
Sample 4	170	96,4%	105,2%

As NETs can be associated with very high concentrations of CGA, out-of-range samples must be diluted. To ensure that dilutions do not compromise the accuracy of the results provided, cascade dilution tests were performed (1/2; 1/4, 1/8, 1/16).

The results show very good dilution performances with a recovery close to 1.0, a slope interpolation close to 1.0 and a correlation coefficient also close to 1.0

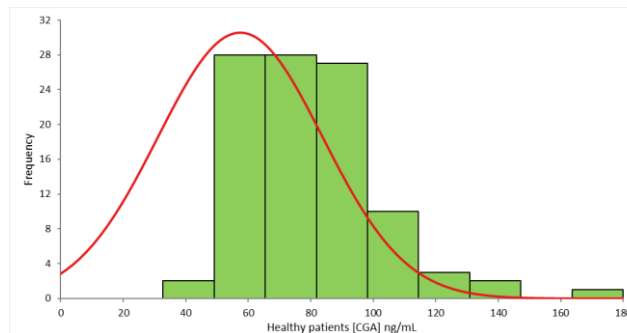
	Dilutions			
	[CGA] ng /mL	Average % Recovery	Interp. Slope	Correl Coef
Sample 1	615	93,8%	1,003	0,999
Sample 2	405	98,7%	1,006	1,000
Sample 3	845	96,0%	1,000	1,000
Sample 4	902	88,5%	0,999	0,999

Biotin interferences:

Due to the resurgence of high serum biotin levels as a result of treatments or dietary supplements, the kit has been evaluated and no significant interference up to biotin concentrations of 600 ng/mL was shown

Normal Values:

Based on a serum library of 101 healthy samples, the distribution of normal values is shown below. The 95th percentile value that can be considered as a cut-off value is estimated at 101 ng/mL



Clinical validation of the kit:

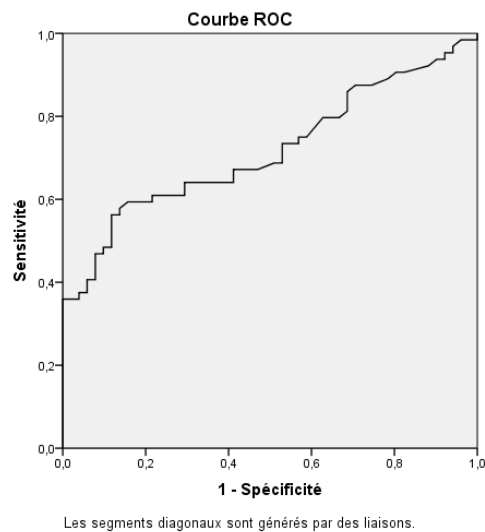
A retrospective study was carried out in partnership with the medical biology laboratory of the Hospices Civils de Lyon under the aegis of Pr Thomas Walter, specialist in neuroendocrine tumors of the digestive system (Report available on request).

The table below provides the precise type of populations studied (positive and negative groups) in terms of pathology or treatments administered.

Variables	NEN patients, n=229		« Control », n=100		p
	n available	n available	n available	n available	
Female, n (%)	229	107 (47%)	100	52 (52%)	0.38
Age in years, median (range)	229	65 (19-93)	100	61 (18-85)	<0.001
Chronic gastritis, n (%)	229	3 (1%)	100	3 (3%)	0.29
Chronic congestive heart failure, n (%)	229	0 (0%)	100	1 (1%)	0.13
Renal deficiency, n (%)	222	39 (18%)	93	17 (18%)	0.88
Inflammatory syndrome, n (%)	147	57 (39%)	63	42 (67%)	<0.001
Concurrent PPI, n (%)	229	44 (19%)	100	40 (40%)	<0.001
MEN1 syndrome, n (%)	227	14 (6%)	100	0 (0%)	<0.001
Concurrent somatostatin analogs, n (%)	229	123 (54%)	100	0 (0%)	<0.001
Biological parameters, median (range)					
CRP in mg/mL	147	3 (1-162)	63	11 (1-130)	0.003
GFR (CKD-EPI) in mL/mn	221	88 (7 to>90)	92	>90 (8 to>90)	0.56
Median CgA by RIA in ng/mL (range)	229	176 (20-431,846)	23	102 (35-2,365)	0.58
Median CgA by ELISA (range)	229	155 (0-227,586)	100	82 (18-1,995)	0.09
CgA ELISA	229		100		<0.001
< 1 ULN (≤ 94 ng/mL)		87 (38%)		56 (56%)	
1-2 ULN (95-188 ng/mL)		33 (14%)		20 (20%)	
> 2 ULN (>188 ng/mL)		109 (48%)		24 (24%)	

NEN, neuroendocrine neoplasm; GFR, glomerular filtration rate; CRP, C-reactive protein; CgA, chromogranin A; PPI, proton pump inhibitors.

At the end of this study, a ROC curve was performed excluding patients with at least 1 false positive risk factor (heart failure, renal failure, PPI, chronic gastritis and NEM 1) and those on somatostatin analog treatments (risk of false negative). A total of 64 NET patients and 51 Controls were taken into consideration. Other classical parameters were also calculated (Specificity, Sensitivity, AUC, PPV, NPV



- Power of the study: > 80%
- Significance of the study: $p < 0,05$
- Specificity=0,78
- Sensitivity=0,61
- AUC=0,72 (0,62-0,81)
- PPV=0,59
- NPV=0,80

These data confirm those of the international recommendations demonstrating the interest of the assay in the follow-up of GEP-NET (prognostic biomarker and early marker of response to a given treatment)

Bibliography:

Tsai et al,

The prognostic and predictive role of CGA in gastroenteropancreatic neuroendocrine tumors- A single center experience
Forntiers in oncology, 11:article 741096, 2021

Dam et al,

Prospective study of CGA as a predictor of progression in patients with pancreatic, small intestine and unknown primary neuroendocrine tumors
Neuroendocrinology 110: 217-224, 2020

Zhang C. et al,

Serum chromogranin A for the diagnosis of gastroenteropancreatic neuroendocrine neoplasms and its association with tumour expression.
Oncology Letters 17: 1497-1504, 2019

Jun E et al.

Diagnostic value of chromogranin A in pancreatic neuroendocrine tumors depends on tumor size: A prospective observational study from a single institute.
Surgery. 2017 Jul;162(1):120-30

Rogowski W et al.

Baseline chromogranin A and its dynamics are prognostic markers in gastroenteropancreatic neuroendocrine tumors.

Future Oncol. 2017;13(12):1069-79

Cheng Y et al.

Serum chromogranin A levels for the diagnosis and follow-up of well-differentiated non-functioning neuroendocrine tumors.

Tumour biology. 2016; 37(3):2863-9

d'Herbomez M et al.

Biomarkers of neuroendocrine tumors.

Ann Biol Clin. 2016; 74(6):669-79.

Erickson JA et al.

A chromogranin A ELISA absent of an apparent high-dose hook effect observed in other chromogranin A ELISAs.

Clin Chim Acta. 2016; 452:120-3

Gut P et al.

Chromogranin A - unspecific neuroendocrine marker. Clinical utility and potential diagnostic pitfalls.

Arch Medical Sci: AMS. 2016; 12(1):1-9

Kim M et al.

The Role of Plasma Chromogranin A as Assessment of Treatment Response in Non-functioning Gastroenteropancreatic Neuroendocrine Tumors.

Cancer Res treat. 2016; 48(1):153-61

Lyubimova NV et al.

Chromogranin As a Biochemical Marker of Neuroendocrine Tumors.

Bull Exp Biol Med. 2016; 160(5):702-4

Shanahan MA et al.

Chromogranin A predicts survival for resected pancreatic neuroendocrine tumors.

J Surg Res. 2016; 201(1):38-43

Glinicki P et al.

Comparison of chromogranin A (CgA) levels in serum and plasma (EDTA2K) and the respective reference ranges in healthy males.

Endokrynol Pol. 2015; 66(1):53-6.

Hallet J et al.

Exploring the rising incidence of neuroendocrine tumors: a population-based analysis of epidemiology, metastatic presentation, and outcomes.

Cancer. 2015; 121(4):589-97

Han X et al.

The value of serum chromogranin A as a predictor of tumor burden, therapeutic

response, and nomogram-based survival in well-moderate nonfunctional pancreatic neuroendocrine tumors with liver metastases.

Eur J Gastroenterol Hepatol. 2015; 27(5):527-35.

Kim M et al.

The Role of Plasma Chromogranin A as Assessment of Treatment Response in Non-functioning Gastroenteropancreatic Neuroendocrine Tumors.

Cancer Res Treat. 2016; 48(1):153-61

Attwood SE et al.

Long-term safety of proton pump inhibitor therapy assessed under controlled, randomised clinical trial conditions: data from the SOPRAN and LOTUS studies.

Aliment Pharmacol Ther. 2015; 41(11):1162-74.

Rehfeld JF.

Chromogranin A in gastrinomas: Promises and pitfalls.

Clin Chim Acta. 2015; 15;446:15-20.

P Glinicki et al.

Chromogranin A (CgA): structure, biological function, pre-analytical, analytical, and clinical aspects of its measurement in blood

Postepy Nauk Medycznych. 2014; XXVII(12):847-51.

Piotr Glinicki et al.

Comparison of chromogranin A levels in serum and plasma (EDTA2K) and the respective reference ranges in healthy males

Endocrine Abstracts. 2014; (35):532

Hijioka M et al.

Serum chromogranin A is a useful marker for Japanese patients with pancreatic neuroendocrine tumors.
Cancer Sci. 2014; 105(11):1464-71.

Modlin IM et al.

Neuroendocrine tumor biomarkers: current status and perspectives.
Neuroendocrinology. 2014; 100(4):265-77.

Onal IK et al.

Chromogranin A as a marker of disease activity in inflammatory bowel disease.
Scand J Gastroenterol. 2014; 49(12):1501-2.

Pedersen L et al.

Preanalytical factors of importance for measurement of Chromogranin A.
Clin Chim Acta. 2014; 436:41-4.