

Human Neurogranin ELISA Kit (NRGN) ab277083

重组 SimpleStep ELISA

5 图像

概述

产品名称 人Neurogranin ELISA试剂盒(NRGN)

检测方法 Colorimetric

批次内				
样品	n	Mean	SD	CV%
Plasma	8			7.2%

批次间				
样品	n	Mean	SD	CV%
Plasma	3			14.9%

样品类型 Cell culture supernatant, Serum, EDTA Plasma, Cit plasma

检测类型 Sandwich (quantitative)

灵敏度 1.4 pg/ml

范围 9.375 pg/ml - 600 pg/ml

特定样本回收率		
样品类型	平均%	范围
Cell culture supernatant	76	72% - 78%
Serum	86	80% - 93%
EDTA Plasma	96	88% - 103%
Cit plasma	104	98% - 109%

检测时间 1h 30m

实验步骤 One step assay

种属反应性 与反应: Human

产品概述

Human Neurogranin ELISA kit (ab277083) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Human Neurogranin protein in human serum, plasma - citrate, plasma EDTA, and cell culture supernatant samples. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human Neurogranin with 1.4 pg/mL sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

-Single-wash protocol reduces assay time to 90 minutes or less

-High sensitivity, specificity and reproducibility from superior antibodies

-Fully validated in biological samples

-96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate (**ab203359**) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

说明

Neurogranin (NRGN), encoded by the *NRGN* gene, is a postsynaptic protein kinase that is highly expressed in the brain and regulates the availability of calmodulin; neurogranin binds to calmodulin in the absence of calcium. Phosphorylation by protein kinase C has been observed to decrease the calmodulin binding activity. Studies have previously shown that *NRGN* expression is regulated during development by thyroid hormones and retinoids for proper development of the central nervous system. ab277083 has been developed to specifically recognize full length neurogranin.

平台

Pre-coated microplate (12 x 8 well strips)

性能

存放说明

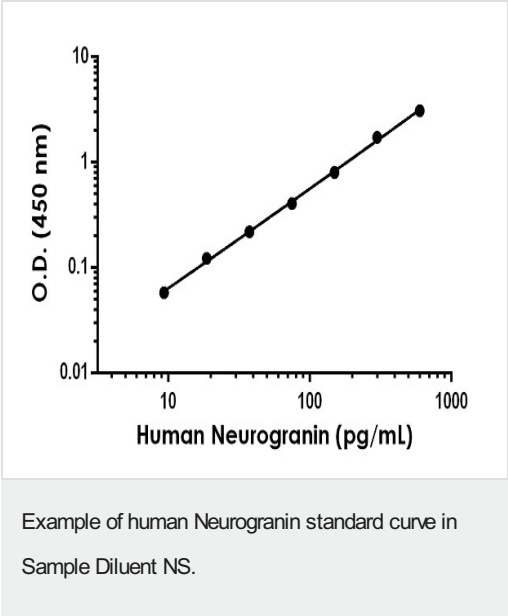
Store at +4°C. Please refer to protocols.

组件	1 x 96 tests
10X Human Neurogranin (NRGN) Capture Antibody	1 x 600µl
10X Human Neurogranin (NRGN) Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 4BI	1 x 6ml
Human Neurogranin (NRGN) Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit

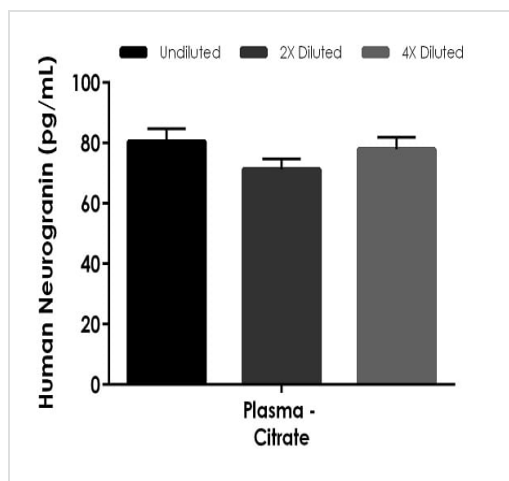
组件	1 x 96 tests
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

功能	Acts as a "third messenger" substrate of protein kinase C-mediated molecular cascades during synaptic development and remodeling. Binds to calmodulin in the absence of calcium.
组织特异性	In the cerebral cortex, found in the cell bodies of neurons in layers II-VI, and in apical and basal dendrites of pyramidal neurons. Is not found in the dendrites in patients with Alzheimer disease.
序列相似性	Belongs to the neurogranin family. Contains 1 collagen-like domain. Contains 1 IQ domain.

图片

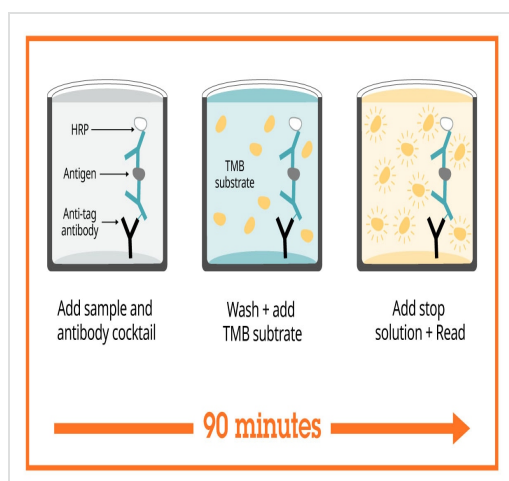


The Neurogranin standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



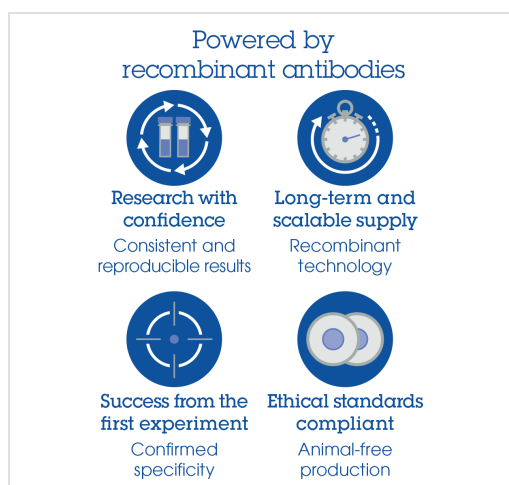
Interpolated concentrations of native Neurogranin in human plasma samples.

The concentrations of Neurogranin were measured in duplicates, interpolated from the Neurogranin standard curves and corrected for sample dilution. Undiluted samples are as follows: plasma (citrate) 100% and plasma (heparin) 12.5%. The interpolated dilution factor corrected values are plotted (mean \pm SD, $n=2$). The mean Neurogranin concentration was determined to be 77 pg/mL in plasma (citrate).



SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.

Sandwich ELISA - Human Neurogranin ELISA Kit
(NRGN) (ab277083)



To learn more about the advantages of recombinant antibodies see [here](#).

Sandwich ELISA - Human Neurogranin ELISA Kit
(NRGN)

**Get more done with
SimpleStep ELISA**



Easy to use
Single-wash 90-minute
protocol



Flexible
Matched antibody pairs
available



Precision antibodies
High sensitivity, specificity
and reproducibility



Scalable
Now in 10-pack and
384-well formats

Sandwich ELISA - Human Neurogranin ELISA Kit
(NRGN) (ab277083)

To learn more about the advantages of SimpleStep ELISA® kits
see [here](#).

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors