

Human NUCB2 ELISA Kit ab213812

1 References 1 Image

Overview

Product name Human NUCB2 ELISA Kit

Detection method Colorimetric

Precision Intra-assay

Sample	n	Mean	SD	CV%
1	16	58pg/ml	2.55	= 4.4%
2	16	278pg/ml	17.51	= 7.2%
3	16	993pg/ml	57.2	= 5.6%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	62pg/ml	3.41	= 5.5%
2	24	286pg/ml	20.59	= 7.2%
3	24	1045pg/ml	74.19	= 7.1%

Sample type Cell culture supernatant, Serum

Assay type Sandwich (quantitative)

Sensitivity < 10 pg/ml

Range 31.2 pg/ml - 2000 pg/ml

Assay time 3h 30m

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Human

Product overview The Human NUCB2 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213812) is designed for the quantitative detection of Human NUCB2 in cell culture supernatants and serum.

The ELISA kit is based on standard sandwich enzyme-linked immunosorbent assay technology. A monoclonal antibody from mouse specific for NUCB2 has been pre-coated onto 96-well plates.

Standards (Expression system for standard: E.coli; Immunogen sequence: V25-L106) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for NUCB2 is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was added and unbound conjugates were washed away with PBS or TBS buffer. HRP substrate TMB was used to visualize HRP enzymatic reaction. TMB was catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the Human NUCB2 amount of sample captured in plate.

**Notes** NUCB2 is a naturally occurring protein molecule produced by the brains of mammals. It is responsible for regulating appetite and production of body fat. NUCB2 is a metabolic polypeptide encoded in the N-terminal region of the protein precursor, Nucleobindin2 (NUCB2). Recombinant human NUCB2 is a 9.7 kDa protein containing 82 amino acid residues. Originally identified as a hypothalamic neuropeptide, NUCB2 is also expressed in other areas of the brain, and in pancreatic islets of Langerhans, gastric endocrine cells and adipocytes. NUCB2 plays a role in hunger and energy regulation in an independent manner.

**Platform** Pre-coated microplate (12 x 8 well strips)

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	Identifier	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Adhesive Plate Seal		4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-Human NUCB2 coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti-Human NUCB2 Antibody		1 x 100µl
Lyophilized Recombinant Human NUCB2 Standard		2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml
Wash Buffer (25X)		1 x 20ml

**Function** Calcium-binding protein. May have a role in calcium homeostasis.

**Tissue specificity** Predominantly expressed in spleen, testis and normal stomach.

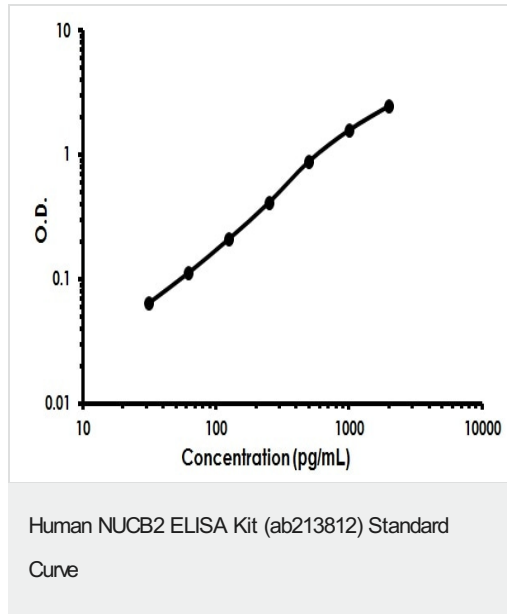
**Sequence similarities** Belongs to the nucleobindin family.

Contains 2 EF-hand domains.

## Cellular localization

Membrane. Cytoplasm. Secreted.

## Images



**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.com/abpromise> or contact our technical team.

## Terms and conditions

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