

Rat BNP 45 ELISA Kit **ab108816**

★★★★★ [1 Abreviews](#) [23 References](#) [2 Images](#)

Overview

Product name	Rat BNP 45 ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			4.9%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			7.1%
Sample type	Serum, Plasma			
Assay type	Sandwich (quantitative)			
Sensitivity	= 7.9 ng/ml			
Range	15.625 pg/ml - 1000 pg/ml			
Assay duration	Multiple steps standard assay			
Species reactivity	<b>Reacts with:</b> Rat			
Product overview	Abcam's BNP 45 rat <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of BNP 45 in plasma, serum, tissue extract, and cell culture supernatants.			

A BNP 45 specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a BNP 45 specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of BNP 45 captured in plate.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid

**repeated freeze-thaw cycles.**

**Notes**

There is significant cross-reactivity between BNP32 and BNP45.

**Platform**

Microplate

**Properties**

**Storage instructions**

Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
70X Biotinylated Rat BNP 45 Antibody	1 x 90µl
BNP 45 Microplate (12 x 8 well strips)	1 unit
BNP 45 Standard (Lyophilized)	1 vial
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml

**Function**

Cardiac hormone which may function as a paracrine antifibrotic factor in the heart. Also plays a key role in cardiovascular homeostasis through natriuresis, diuresis, vasorelaxation, and inhibition of renin and aldosterone secretion. Specifically binds and stimulates the cGMP production of the NPR1 receptor. Binds the clearance receptor NPR3.

**Tissue specificity**

Brain and also in atria, but at much lower levels than ANP.

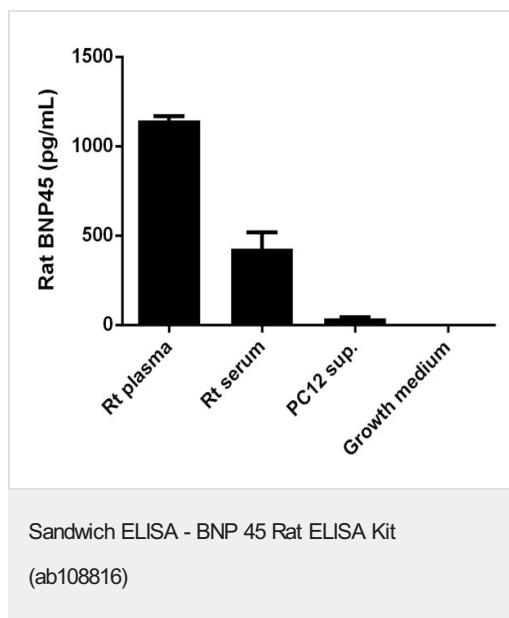
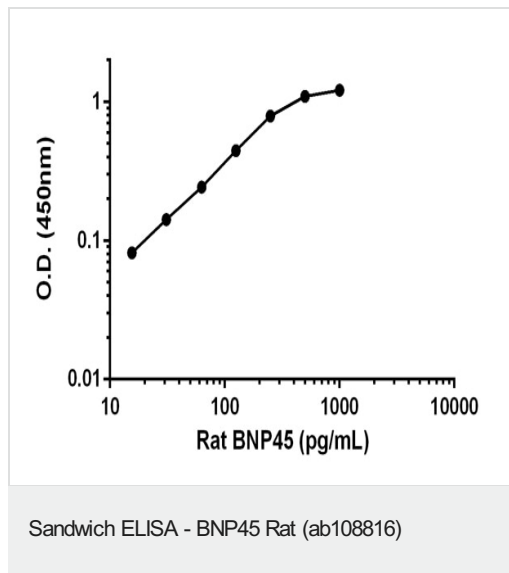
**Sequence similarities**

Belongs to the natriuretic peptide family.

**Cellular localization**

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