

Rat BNP 32 ELISA Kit ab108815

10 References 1 Image

Overview

Product name	Rat BNP 32 ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			4.9%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			10.7%
Sample type	Cell culture supernatant, Serum, Plasma, Tissue			
Assay type	Sandwich (quantitative)			
Sensitivity	= 18 pg/ml			
Range	0.031 ng/ml - 2 ng/ml			
Recovery	= 103 %			
Assay time	5h 00m			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Rat			
Product overview	Abcam's BNP 32 Rat <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of BNP 32 in plasma, serum, tissue extract, and cell culture supernatants.			

A BNP 32 specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a BNP 32 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 32 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
50X Biotinylated Rat BNP 32 Antibody	1 x 120µl
BNP 32 Microplate (12 x 8 well strips)	1 unit
BNP 32 Standard	1 vial
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml

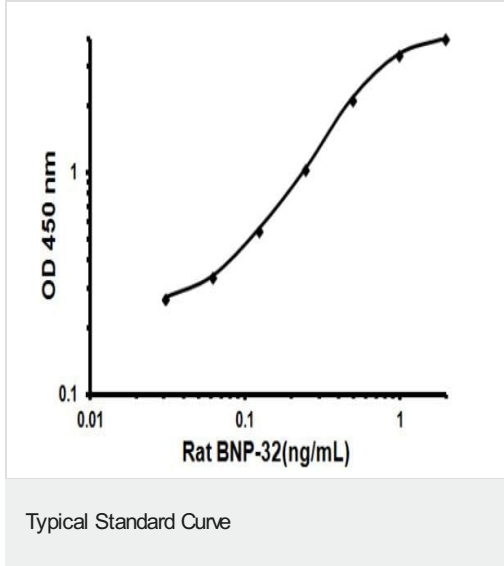
Function Cardiac hormone which may function as a paracrine antifibrotic factor in the heart. Also plays a key role in cardiovascular homeostasis through natriuresis, diuresis, vasorelaxation, and inhibition of renin and aldosterone secretion. Specifically binds and stimulates the cGMP production of the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



Representative Standard Curve using ab108815.

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