

Rat BNP 45 ELISA Kit ab108816

★★★★★ 1 Abreviews 23 References 2 图像

概述

产品名称 大鼠BNP 45 ELISA试剂盒

检测方法 Colorimetric

批次内				
样品	n	Mean	SD	CV%
Overall				4.9%

批次间				
样品	n	Mean	SD	CV%
Overall				7.1%

样品类型 Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 = 7.9 ng/ml

范围 15.625 pg/ml - 1000 pg/ml

实验步骤 Multiple steps standard assay

种属反应性 与反应: Rat

产品概述

Abcam's BNP 45 rat *in vitro* 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.

说明	There is significant cross-reactivity between BNP32 and BNP45.
平台	Microplate

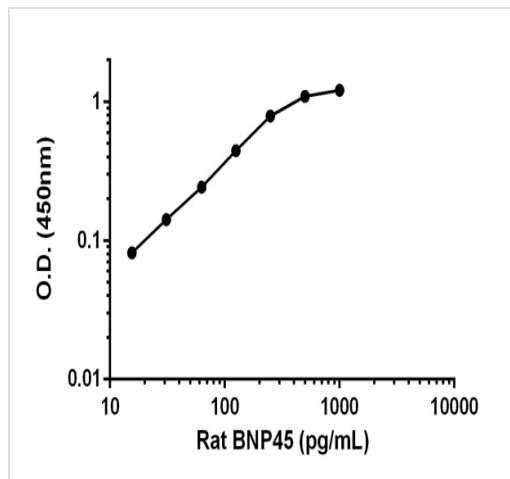
性能

存放说明 Store at -20°C. Please refer to protocols.

组件	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

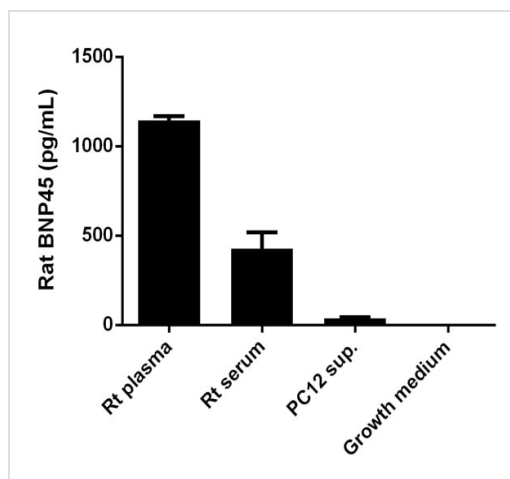
功能	Cardiac hormone which may function as a paracrine antifibrotic factor in the heart. Also plays a key role in cardiovascular homeostasis through natriuresis, diuresis, vasorelaxation, and inhibition of renin and aldosterone secretion. Specifically binds and stimulates the cGMP production of the NPR1 receptor. Binds the clearance receptor NPR3.
组织特异性	Brain and also in atria, but at much lower levels than ANP.
序列相似性	Belongs to the natriuretic peptide family.
细胞定位	Secreted.

图片



Standard curve : mean of duplicates (+/-SD) with background readings subtracted

Sandwich ELISA - BNP45 Rat (ab108816)



Rat BNP45 measured in biological fluids showing quantity (pg) per mL of tested sample

Sandwich ELISA - BNP 45 Rat ELISA Kit (ab108816)

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