

Product datasheet

Human BNP ELISA Kit ab193694

1 图像

概述

产品名称 人BNP ELISA试剂盒

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
Overall				< 10%

批次间

样品	n	Mean	SD	CV%
Overall				< 12%

样品类型 Cell culture supernatant, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 14 pg/ml

范围 14 pg/ml - 1000 pg/ml

回收率 特定样本回收率

样品类型	平均%	范围
Serum	78.12	68% - 83%
Plasma	80.96	67% - 107%
Cell culture media	94.31	83% - 105%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's BNP Human ELISA Kit (ab193694) is an in vitro enzyme-linked immunosorbent assay for the quantitative measurement of Human BNP in serum, plasma and cell culture supernatants. This assay employs an antibody specific for Human BNP coated on a 96-well plate. Standards and samples are pipetted into the wells and the immobilized antibody captures BNP present in

the samples. The wells are washed and biotinylated anti-Human BNP antibody is added. After washing away any unbound biotinylated antibody, an HRP-conjugated streptavidin is pipetted to the wells. After incubation, the wells are again washed, followed by the addition of a TMB substrate solution to the wells. Color will develop in proportion to the amount of BNP bound in each well. Addition of the Stop Solution will change the color from blue to yellow, and the intensity of the color is measured at 450 nm.

经测试应用

适用于: Sandwich ELISA

平台

Pre-coated microplate (12 x 8 well strips)

性能

存放说明

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

组件	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer Concentrate	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated Human BNP Detection Antibody (lyophilized)	2 vials
Human BNP Standards (lyophilized)	2 vials
Pre-coated Human BNP Microplate (12 strips x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

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.

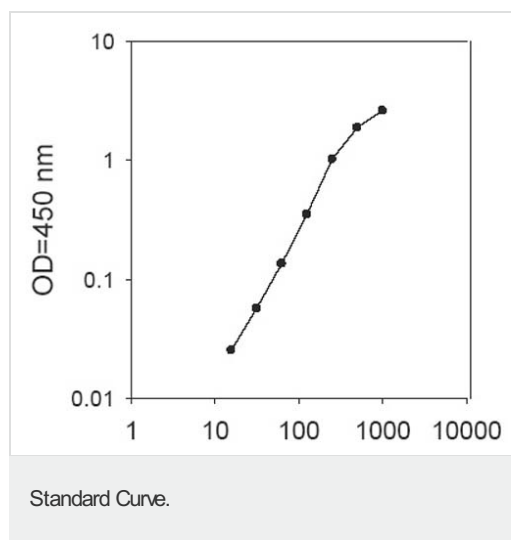
应用

Our [Abpromise guarantee](#) covers the use of **ab193694** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

应用	Ab评论	说明
Sandwich ELISA		Use at an assay dependent concentration.

图片



Standard curve using ab193694 BNP Human ELISA Kit.

Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

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