

Rat ANP ELISA Kit (NPPA) ab108797

★★★★★ [1 Abreviews](#) [11 References](#) [画像数 2](#)

製品の概要

製品名	Rat ANP ELISA Kit (NPPA)										
検出方法	Colorimetric										
再現性	Intra-Assay (同時再現性)										
	<table border="1"> <thead> <tr> <th>サンプル</th> <th>N</th> <th>平均値</th> <th>SD</th> <th>CV%</th> </tr> </thead> <tbody> <tr> <td>Overall</td> <td></td> <td></td> <td></td> <td>6.3%</td> </tr> </tbody> </table>	サンプル	N	平均値	SD	CV%	Overall				6.3%
サンプル	N	平均値	SD	CV%							
Overall				6.3%							
	Inter-Assay (日差再現性)										
	<table border="1"> <thead> <tr> <th>サンプル</th> <th>N</th> <th>平均値</th> <th>SD</th> <th>CV%</th> </tr> </thead> <tbody> <tr> <td>Overall</td> <td></td> <td></td> <td></td> <td>10.7%</td> </tr> </tbody> </table>	サンプル	N	平均値	SD	CV%	Overall				10.7%
サンプル	N	平均値	SD	CV%							
Overall				10.7%							
サンプルの種類	Cell culture supernatant, Serum, Plasma, Tissue										
アッセイタイプ	Sandwich (quantitative)										
検出感度	= 6.1 pg/ml										
検出範囲	0.008 ng/ml - 2 ng/ml										
添加回収試験	102 %										
全工程の試験時間	5h 00m										
ステップ	Multiple steps standard assay										
種交差性	交差種: Rat 交差が予測される動物種: Pig 										

製品の概要
 Abcam's ANP (NPPA) Rat *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of ANP concentrations in cell culture supernatants, tissue-extract, serum and plasma.

An ANP specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an ANP 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 ANP captured in plate.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

試験プラットフォーム

Microplate

製品の特性

保存方法

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

内容	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
40X Biotinylated Rat ANP Antibody	1 x 150µl
ANP Microplate (12 x 8 well strips)	1 unit
ANP Standard	1 vial
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml

機能

Hormone playing a key role in cardiovascular homeostasis through regulation of natriuresis, diuresis, and vasodilation. Also plays a role in female pregnancy by promoting trophoblast invasion and spiral artery remodeling in uterus. Specifically binds and stimulates the cGMP production of the NPR1 receptor. Binds the clearance receptor NPR3.

関連疾患

Atrial standstill 2
Atrial fibrillation, familial, 6

配列類似性

Belongs to the natriuretic peptide family.

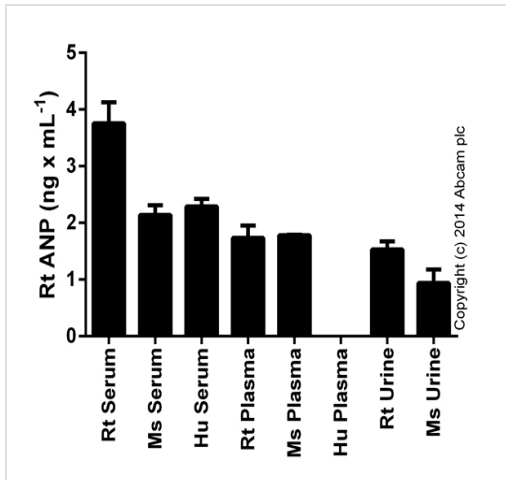
翻訳後修飾

Cleaved by CORIN upon secretion to produce the functional hormone.
Atrial natriuretic factor: Cleaved by MME. The cleavage initiates degradation of the factor and thereby regulate its activity.

細胞内局在

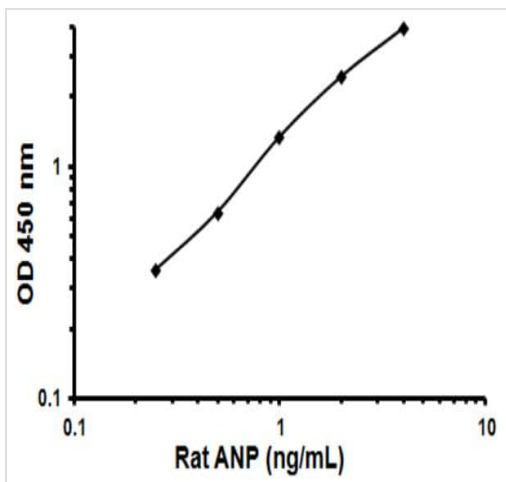
Secreted.

画像



Rat ANP measured in biological fluids, background signal subtracted (duplicates +/- SD).

Sandwich ELISA - ANP (NPPA) Rat ELISA kit (ab108797)



Representative Standard Curve using ab108797.

Typical Standard Curve

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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