

Human Beta Nerve Growth Factor ELISA Kit (NGFB) ab99986

★★★★★ 6 Abreviews 12 References 画像数 1

製品の概要

製品名	Human Beta Nerve Growth Factor ELISA Kit (NGFB)
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 14 pg/ml
検出範囲	6.86 pg/ml - 5000 pg/ml
添加回収試験	> 100 %

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	106.7	90% - 122%
Serum	113.1	98% - 116%
Plasma	102.8	85% - 117%

ステップ	Multiple steps standard assay
種交差性	交差種: Human
製品の概要	Abcam's Beta Nerve Growth Factor (NGFB) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human Beta Nerve Growth Factor in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human Beta Nerve Growth Factor coated on a 96-well plate. Standards and samples are pipetted into the wells and Beta Nerve Growth Factor present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Beta Nerve Growth Factor antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of Beta Nerve Growth Factor bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Get results in 90 minutes with Human beta Nerve Growth Factor ELISA Kit ([ab193760](#)) from our

SimpleStep ELISA® range.

特記事項

NGFB concentration is very low in normal human serum samples and may be undetectable even with undiluted samples. Optimization may be required with urine samples.

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
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
800X HRP-Streptavidin Concentrate	1 x 200µl
Beta Nerve Growth Factor Microplate (12 x 8 wells)	1 unit
Beta Nerve Growth Factor Standard Human (recombinant) (lyophilized)	2 vials
Biotinylated anti-Human Beta Nerve Growth Factor	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

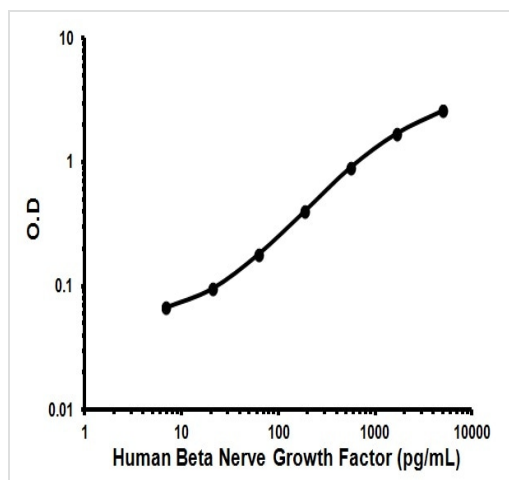
関連性

Nerve growth factor (NGF) is one of a family of neurotrophins that induce the survival and proliferation of neurons. In cell culture, NGF induces the formation of neurite projections and, in vivo, may stimulate the innervation of tissues. NGF plays a role in the repair, regeneration, and protection of neurons, and as such could serve as a therapeutic agent in neurodegenerative conditions such as Alzheimer's disease. NGF enhances survival, growth, neurotransmitter biosynthesis of sympathetic and sensory neurons; neurotrophic factor; cutaneous innervation; growth, differentiation and survival of B lymphocytes. It also has a possible role in allergy and tissue repair. NGF is found in the hypothalamus, pituitary, thyroid gland, testes, epididymis, vascular smooth muscle cells, fibroblasts, mast cells and eosinophils. NGF is upregulated by glutamate, vitamin D3, IL6, FGF basic, astrocyte specific IL1, TNF alpha, PDGF and TGF beta. It is downregulated by GABAergic neuronal activity, glucocorticoids and Schwann cell-specific TGF beta.

細胞内局在

Secreted.

画像



Representative Standard Curve using ab99986.

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