

Human CNTF ELISA Kit ab99994

1 References 画像数 2

製品の概要

製品名	Human CNTF ELISA Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 0.2 ng/ml
検出範囲	0.823 ng/ml - 200 ng/ml
添加回収試験	> 92 %

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	95.45	84% - 104%
Plasma	93.28	83% - 104%

ステップ	Multiple steps standard assay
種交差性	交差種: Human 非交差種: Rat

**製品の概要**

Abcam's CNTF Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human CNTF in plasma and cell culture supernatants. (Human CNTF concentration is pretty low in normal plasma, it may not be detected in this assay).

This assay employs an antibody specific for Human CNTF coated on a 96-well plate. Standards and samples are pipetted into the wells and CNTF present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human CNTF 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 CNTF bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

We have not been able to detect endogenous Human CNTF in normal serum with ab99994, only in serum spiked with Human CNTF.

**特記事項** Optimization may be required with urine samples.

**試験プラットフォーム** Microplate

## 製品の特性

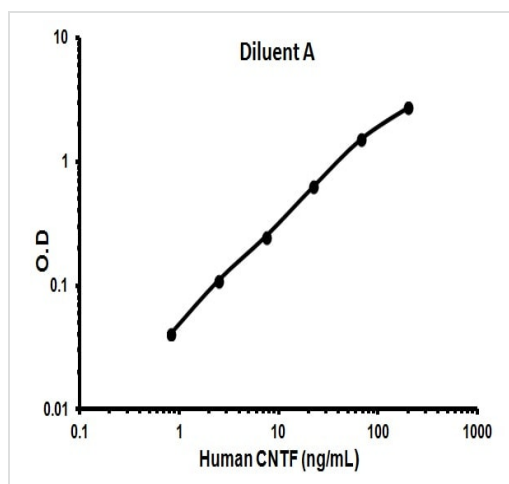
**保存方法** Store at -20°C. Please refer to protocols.

内容	1 x 96 tests
20X Wash Buffer	1 x 25ml
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human CNTF	2 vials
CNTF Microplate (12 strips x 8 wells)	1 unit
Recombinant Human CNTF Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**関連性** Ciliary neurotrophic factor (CNTF) is expressed in glial cells within the central and peripheral nervous systems. CNTF stimulates gene expression, cell survival or differentiation in a variety of neuronal cell types such as sensory, sympathetic, ciliary and motor neurons. CNTF itself lacks a classical signal peptide sequence of a secreted protein, but is thought to convey its cytoprotective effects after release from adult glial cells by some mechanism induced by injury. In addition to its neuronal actions, CNTF also acts on non neuronal cells such as glia, hepatocytes, skeletal muscle, embryonic stem cells and bone marrow stromal cells.

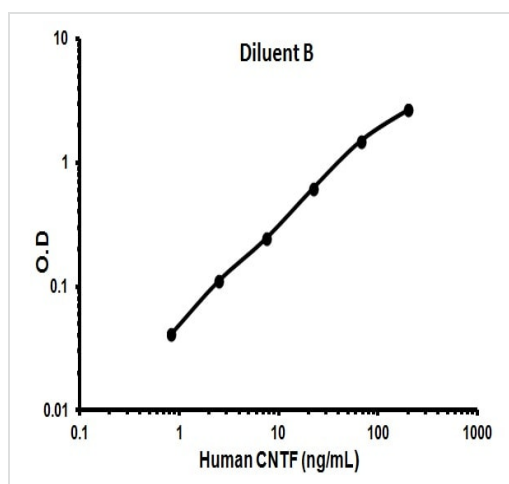
**細胞内局在** Cytoplasmic

## 画像



Representative Standard Curve using ab99994

Typical Standard Curve



Representative Standard Curve using ab99994

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