

Human NGF R ELISA Kit (CD271) ab155436

[1 References](#) [1 Image](#)

Overview

Product name Human NGF R ELISA Kit (CD271)

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Intra-assay				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
Inter-assay				< 12%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity < 80 pg/ml

Range 0.082 ng/ml - 20 ng/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	86.88	79% - 95%
Serum	82.75	75% - 93%
Plasma	84.17	76% - 93%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's NGF R (CD271) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay designed for the quantitative measurement of Human NGF R (CD271) in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Human NGF R coated on a 96-well plate. Standards

and samples are pipetted into the wells and NGF R present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human NGF R 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 NGF R bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Notes Optimization may be required with urine samples.
The standard is affinity purified.

Platform Microplate

Properties

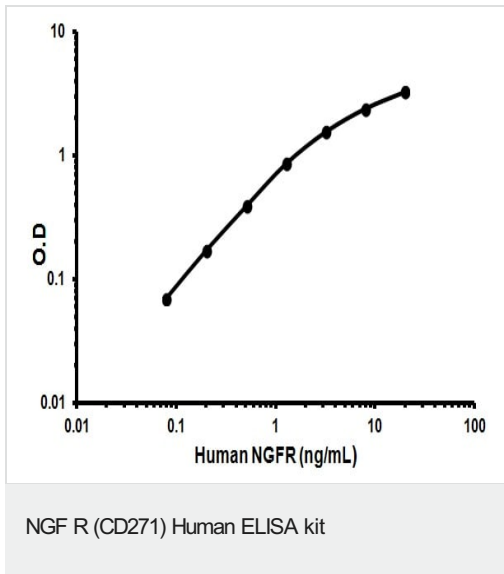
Storage instructions Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
500X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent C	1 x 30ml
Biotinylated anti-Human NGF R	2 vials
NGF R Microplate (12 x 8 wells)	1 unit
Recombinant Human NGF R Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Relevance Plays a role in the regulation of the translocation of GLUT4 to the cell surface in adipocytes and skeletal muscle cells in response to insulin, probably by regulating RAB31 activity, and thereby contributes to the regulation of insulin-dependent glucose uptake (By similarity). Low affinity receptor which can bind to NGF, BDNF, NT-3, and NT-4. Can mediate cell survival as well as cell death of neural cells. Necessary for the circadian oscillation of the clock genes ARNTL/BMAL1, PER1, PER2 and NR1D1 in the suprachiasmatic nucleus (SCN) of the brain and in liver and of the genes involved in glucose and lipid metabolism in the liver.

Cellular localization Membrane; Single-pass type I membrane protein.

Images



Representative standard curve using ab155436

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.com/abpromise> or contact our technical team.

Terms and conditions

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