

Human NOV ELISA Kit (Nephroblastoma Overexpressed) ab193710

1 [References](#) 2 [图像](#)

概述

产品名称

检测方法

精确度

样品类型

检测类型

灵敏度

范围

回收率

实验步骤

种属反应性

产品概述

人NOV ELISA试剂盒(Nephroblastoma Overexpressed)

Colorimetric

批次内

样品	n	Mean	SD	CV%
Overall				< 10%

批次间

样品	n	Mean	SD	CV%
Overall				< 12%

Cell culture supernatant, Serum, Plasma

Sandwich (quantitative)

2 pg/ml

2 pg/ml - 30000 pg/ml

特定样本回收率

样品类型	平均%	范围
Serum	109.5	102% - 118%
Plasma	99.29	93% - 106%
Cell culture media	130.4	118% - 139%

One step assay

与反应: Human

Abcam's NOV (Nephroblastoma Overexpressed) Human ELISA Kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of human NOV in serum, plasma, and cell culture supernatant.

This assay employs an antibody specific for human NOV coated on a 96-well plate. Standards and samples are pipetted into the wells and the immobilized antibody captures NOV present in the samples. The wells are washed and biotinylated anti-human NOV 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 NOV 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.

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
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
700X HRP-Streptavidin Concentrate	1 x 200μl
Assay Diluent A	1 x 30ml
Biotinylated Human NOV detection antibody	2 vials
Human NOV Standards (lyophilized)	2 vials
Pre-coated Human NOV Microplate (12 strips x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

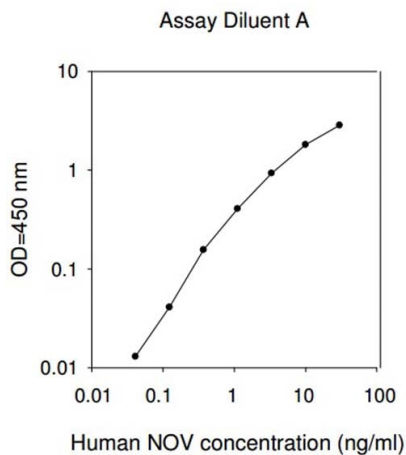
功能 Immediate-early protein likely to play a role in cell growth regulation.

组织特异性 Expressed in bone marrow, thymic cells and neuroblastoma. Increased expression in Wilms tumor of the stromal type.

序列相似性 Belongs to the CCN family.
Contains 1 CTCK (C-terminal cystine knot-like) domain.
Contains 1 IGFBP N-terminal domain.
Contains 1 TSP type-1 domain.
Contains 1 VWFC domain.

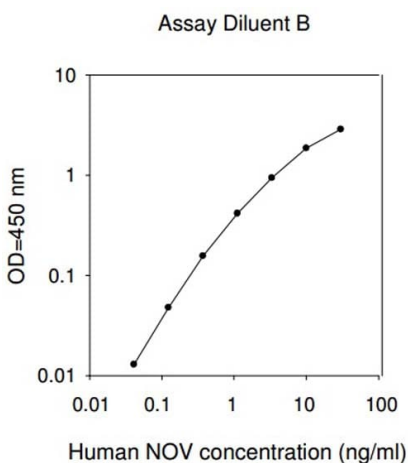
细胞定位 Secreted.

图片



Typical standard curve in Assay Diluent A

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



Typical standard curve in Assay Diluent B

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