

Product datasheet

Mouse NOV ELISA Kit (CCN3) α b205570

Recombinant SimpleStep ELISA[®]

[1 References](#) [7 Images](#)

Overview

Product name Mouse NOV ELISA Kit (CCN3)

Detection method Colorimetric

Precision

Intra-assay

| Sample | n | Mean | SD | CV% |
|---------|---|------|----|------|
| Overall | 5 | | | 8.4% |

Inter-assay

| Sample | n | Mean | SD | CV% |
|---------|---|------|----|------|
| Overall | 3 | | | 5.9% |

Sample type

Cell culture supernatant, Plasma, Cell culture extracts, Tissue Extracts

Assay type

Sandwich (quantitative)

Sensitivity

51 pg/ml

Range

130 pg/ml - 8000 pg/ml

Recovery

Sample specific recovery

| Sample type | Average % | Range |
|--------------------------|-----------|-------------|
| Cell culture supernatant | 111 | 107% - 116% |
| Hep Plasma | 85 | 85% - 85% |
| EDTA Plasma | 88 | 86% - 88% |
| Cit plasma | 104 | 103% - 106% |
| Mouse brain extract | 113 | 108% - 117% |

Assay time

1h 30m

Assay duration

One step assay

Species reactivity**Reacts with:** Mouse**Product overview**

Mouse NOV ELISA Kit (CCN3) (ab205570) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of NOV (CCN3) protein in cell culture supernatant, tissue extracts, plasma, and cell culture extracts. It uses our proprietary SimpleStep ELISA® technology. Quantitate Mouse NOV (CCN3) with 51 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate (**ab203359**) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

Notes

NOV is a member of the CCN family of extracellular matrix associated signaling proteins. As a secreted protein, NOV regulates cell survival, adhesion, proliferation, differentiation and migration. In humans, NOV expression is associated with metastasis and poor prognosis.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

Platform

Pre-coated microplate (12 x 8 well strips)

Properties**Storage instructions**

Store at +4°C. Please refer to protocols.

| Components | 1 x 96 tests |
|--|--------------|
| 10X Mouse NOV Capture Antibody | 1 x 600µl |
| 10X Mouse NOV Detector Antibody | 1 x 600µl |
| 10X Wash Buffer PT (ab206977) | 1 x 20ml |
| 50X Cell Extraction Enhancer Solution (ab193971) | 1 x 1ml |
| 5X Cell Extraction Buffer PTR (ab193970) | 1 x 10ml |
| Antibody Diluent 4BR | 1 x 6ml |
| Mouse NOV Lyophilized Recombinant Protein | 2 vials |

| Components | 1 x 96 tests |
|---|--------------|
| Plate Seals | 1 unit |
| Sample Diluent 75BS | 1 x 20ml |
| Sample Diluent NS (ab193972) | 1 x 50ml |
| SimpleStep Pre-Coated 96-Well Microplate (ab206978) | 1 unit |
| Stop Solution | 1 x 12ml |
| TMB Development Solution | 1 x 12ml |

Function

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

Tissue specificity

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

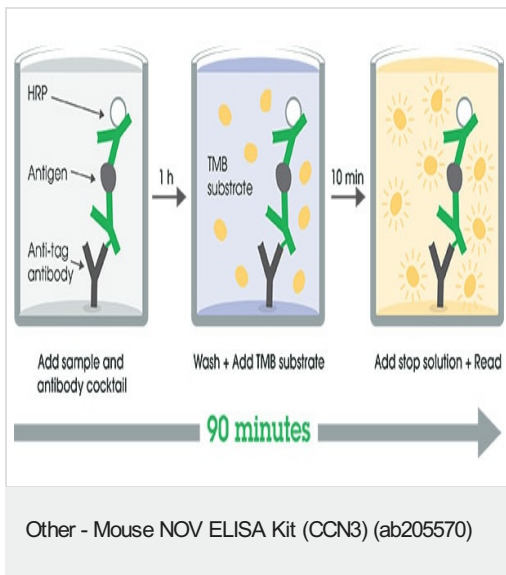
Sequence similarities

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.

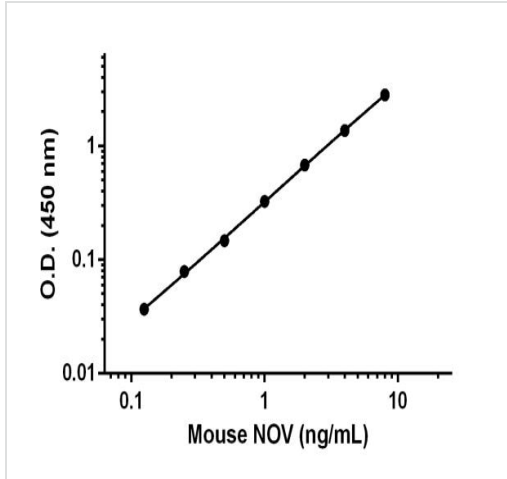
Cellular localization

Secreted.

Images

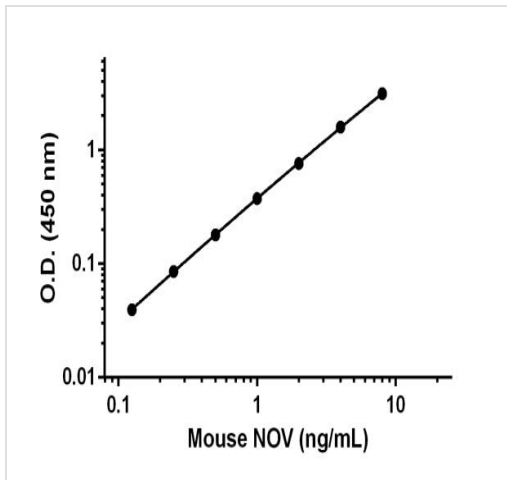


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



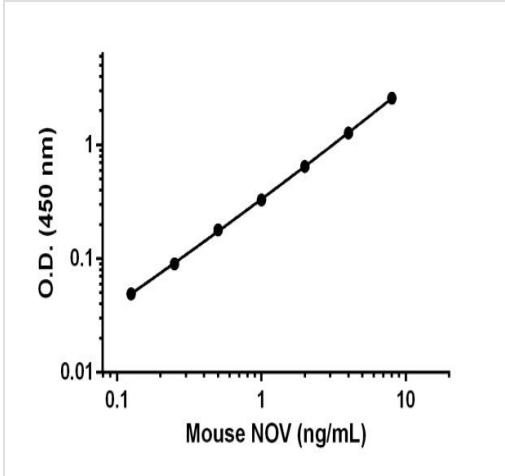
Background-subtracted data values (mean +/- SD) are graphed.

Example of mouse NOV standard curve in Sample Diluent NS.



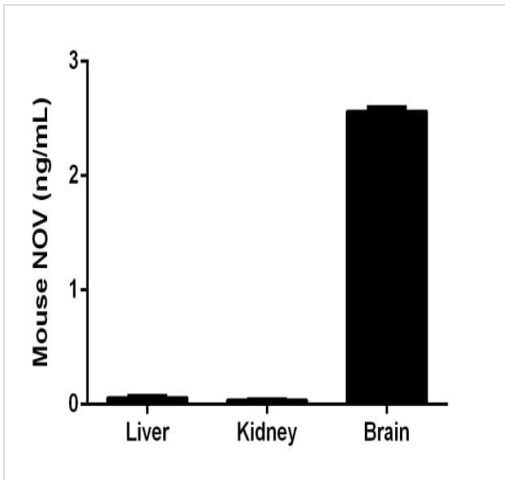
Background-subtracted data values (mean +/- SD) are graphed.

Example of mouse NOV standard curve in 1X Cell Extraction Buffer PTR.



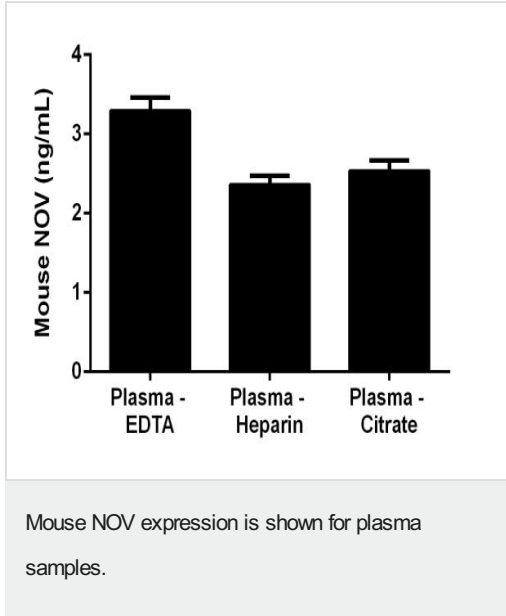
Background-subtracted data values (mean +/- SD) are graphed.

Example of mouse NOV standard curve in Sample Diluent 75BS.



The concentrations of mouse NOV were interpolated from the mouse NOV standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are plotted. Interpolated values of NOV are plotted based on an extract load of 250 μ g/mL.

Linearity of dilution of native mouse NOV protein in mouse tissue extracts.



Native NOV was measured in duplicate in 50% EDTA, heparin and citrate plasma diluted in Sample Diluent 75BS and concentrations interpolated from a standard curve diluted in Sample Diluent 75BS. The interpolated dilution factor corrected values are graphed (mean +/- SD).

Powered by recombinant antibodies

| | |
|--|--|
| <p>Research with confidence Consistent and reproducible results</p> | <p>Long-term and scalable supply Recombinant technology</p> |
| <p>Success from the first experiment Confirmed specificity</p> | <p>Ethical standards compliant Animal-free production</p> |

Sandwich ELISA - Mouse NOV ELISA Kit (CCN3)
(ab205570)

To learn more about the advantages of recombinant antibodies see [here](#).

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