

Product datasheet

Human Tenascin C ELISA Kit ab277081

Recombinant SimpleStep ELISA

[5 Images](#)

Overview

Product name Human Tenascin C ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

| Sample | n | Mean | SD | CV% |
|--------|---|------|----|------|
| Serum | 8 | | | 3.3% |

Inter-assay

| Sample | n | Mean | SD | CV% |
|--------|---|------|----|-------|
| Serum | 3 | | | 12.3% |

Sample type Serum, Hep Plasma, EDTA Plasma, Cit plasma

Assay type Sandwich (quantitative)

Sensitivity 27.672 pg/ml

Range 312.5 pg/ml - 20000 pg/ml

Recovery

Sample specific recovery

| Sample type | Average % | Range |
|-------------|-----------|------------|
| Serum | 97 | 88% - 106% |
| Hep Plasma | 90 | 84% - 94% |
| EDTA Plasma | 98 | 96% - 99% |
| Cit plasma | 93 | 83% - 106% |

Assay time 1h 30m

Assay duration One step assay

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

Human Tenascin C ELISA kit (ab277081) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Human Tenascin C protein in human serum and plasma. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human Tenascin C with 27.67 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

Tenascin C (TNC), is an extracellular matrix glycoprotein involved in tissue morphogenesis like embryogenesis, wound healing and tumorigenesis. Tenascin C is implicated in guidance of migrating neurons and axons during development, synaptic plasticity and neuronal regeneration. Tenascin C also as a role in neurite outgrowth from cortical neurons grown on a monolayer of astrocytes. In tumors, Tenascin C stimulates angiogenesis by elongation, migration and sprouting of endothelial cell. Tenascin C shares 69.0%, 70.6%, 72.0%, and 97.3% sequence homology with mouse, rat, bovine, and monkey, respectively.

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 Human Tenascin C Capture Antibody | 1 x 600µl |
| 10X Human Tenascin C Detector Antibody | 1 x 600µl |
| 10X Wash Buffer PT (ab206977) | 1 x 20ml |
| Antibody Diluent 4BI | 1 x 6ml |
| Human Tenascin C Lyophilized Recombinant Protein | 2 vials |
| Plate Seals | 1 unit |
| Sample Diluent NS (ab193972) | 1 x 50ml |

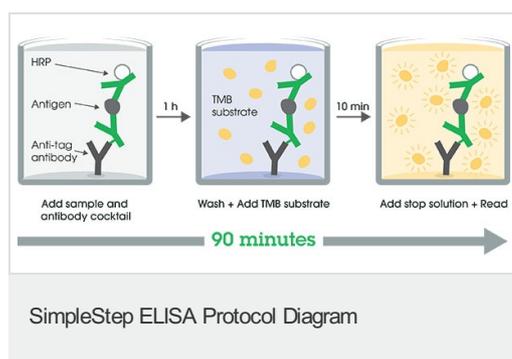
| Components | 1 x 96 tests |
|---|--------------|
| SimpleStep Pre-Coated 96-Well Microplate (ab206978) | 1 unit |
| Stop Solution | 1 x 12ml |
| TMB Development Solution | 1 x 12ml |

Function Extracellular matrix protein implicated in guidance of migrating neurons as well as axons during development, synaptic plasticity as well as neuronal regeneration. Promotes neurite outgrowth from cortical neurons grown on a monolayer of astrocytes. Ligand for integrins alpha-8/beta-1, alpha-9/beta-1, alpha-V/beta-3 and alpha-V/beta-6.

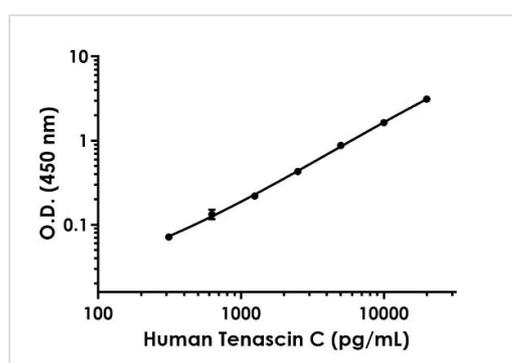
Sequence similarities Belongs to the tenascin family.
 Contains 15 EGF-like domains.
 Contains 1 fibrinogen C-terminal domain.
 Contains 15 fibronectin type-III domains.

Cellular localization Secreted > extracellular space > extracellular matrix.

Images

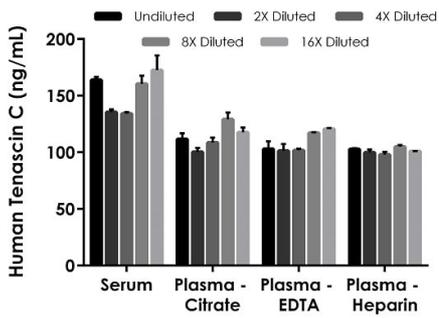


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



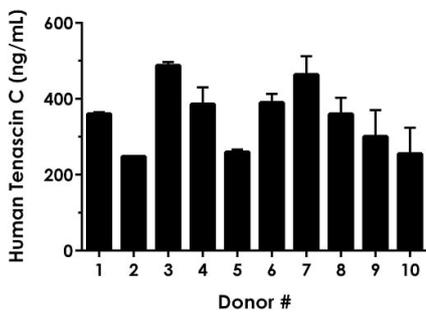
The Tenascin C standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.

Example of human Tenascin C standard curve in Sample Diluent NS. The Tenascin C standard curve was prepared as described in Section 10.



Interpolated concentrations of native Tenascin C in human serum and plasma samples.

The concentrations of Tenascin C were measured in duplicates, interpolated from the Tenascin C standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 10%, plasma (citrate) 10%, plasma (EDTA) 10%, and plasma (heparin) 10%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Tenascin C concentration was determined to be 153.25 ng/mL in serum, 113.37 ng/mL in plasma (citrate), 108.74 ng/mL in plasma (EDTA), and 101.20 ng/mL in plasma (heparin).



5% serum from ten individual healthy human female donors was measured in duplicate.

Interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Tenascin C concentration was determined to be 351.07 ng/mL with a range of 247.61 – 487.97 ng/mL.

Powered by recombinant antibodies



Research with confidence
Consistent and reproducible results



Long-term and scalable supply
Recombinant technology



Success from the first experiment
Confirmed specificity



Ethical standards compliant
Animal-free production

Sandwich ELISA - Human Tenascin C ELISA Kit

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