

Human Entactin ELISA Kit (NID1) ab113338

[1 Image](#)

Overview

Product name	Human Entactin ELISA Kit (NID1)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 0.1 ng/ml
Range	0.123 ng/ml - 30 ng/ml
Recovery	> 100 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	125.7	111% - 138%
Serum	132.3	123% - 145%
Plasma	131.4	122% - 147%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

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

This assay employs an antibody specific for Human Entactin coated on a 96-well plate. Standards and samples are pipetted into the wells and Entactin present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Entactin 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 Entactin 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.

Platform Microplate

Properties

Storage instructions

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

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

Function

Sulfated glycoprotein widely distributed in basement membranes and tightly associated with laminin. Also binds to collagen IV and perlecan. It probably has a role in cell-extracellular matrix interactions.

Sequence similarities

Contains 6 EGF-like domains.
Contains 4 LDL-receptor class B repeats.
Contains 1 NIDO domain.
Contains 1 nidogen G2 beta-barrel domain.
Contains 1 thyroglobulin type-1 domain.

Post-translational modifications

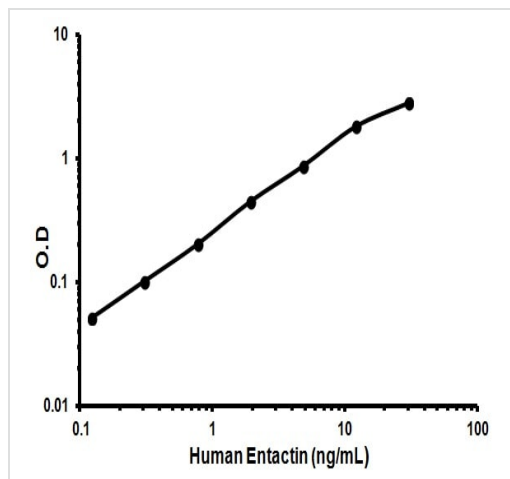
N- and O-glycosylated.

Cellular localization

Secreted > extracellular space > extracellular matrix > basement membrane.

Images

Representative Standard Curve using ab113338



Representative 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.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