

Product datasheet

Human P cadherin ELISA Kit ab100621

[2 Images](#)

Overview

Product name	Human P cadherin ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 20 pg/ml
Range	2.47 pg/ml - 18000 pg/ml
Recovery	> 100 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	129.1	117% - 138%
Serum	121.3	111% - 131%
Plasma	114.9	105% - 125%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

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

This assay employs an antibody specific for Human P cadherin coated on a 96-well plate. Standards and samples are pipetted into the wells and P cadherin present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human P cadherin 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 P cadherin 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
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
600X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Human P cadherin	2 vials
P cadherin Microplate (12 x 8 wells)	1 unit
Recombinant Human P cadherin Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Function Cadherins are calcium dependent cell adhesion proteins. They preferentially interact with themselves in a homophilic manner in connecting cells; cadherins may thus contribute to the sorting of heterogeneous cell types.

Tissue specificity Expressed in some normal epithelial tissues and in some carcinoma cell lines.

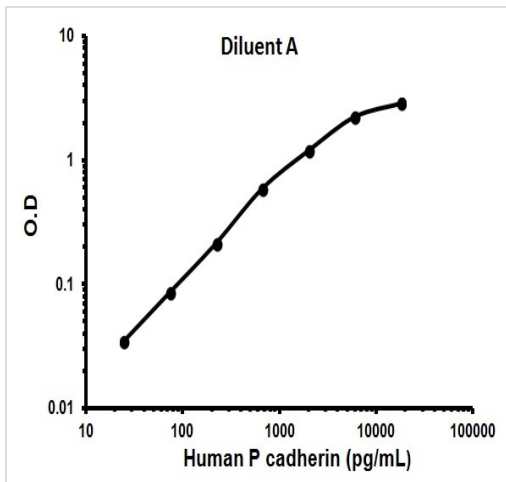
Involvement in disease Defects in CDH3 are the cause of hypotrichosis with juvenile macular dystrophy (HJMD) [MIM:601553]. HJMD is a rare autosomal recessive disorder characterized by early hair loss heralding severe degenerative changes of the retinal macula and culminating in blindness during the second to third decade of life.

Defects in CDH3 are the cause of ectodermal dysplasia with ectrodactyly and macular dystrophy (EEM) [MIM:225280]; also known as EEM syndrome, Albrechtsen-Svendsen syndrome or Ohdo-Hirayama-Terawaki syndrome. Ectodermal dysplasia defines a heterogeneous group of disorders due to abnormal development of two or more ectodermal structures. EEM is an autosomal recessive condition characterized by features of ectodermal dysplasia such as sparse eyebrows and scalp hair, and selective tooth agenesis associated with macular dystrophy and ectrodactyly.

Sequence similarities Contains 5 cadherin domains.

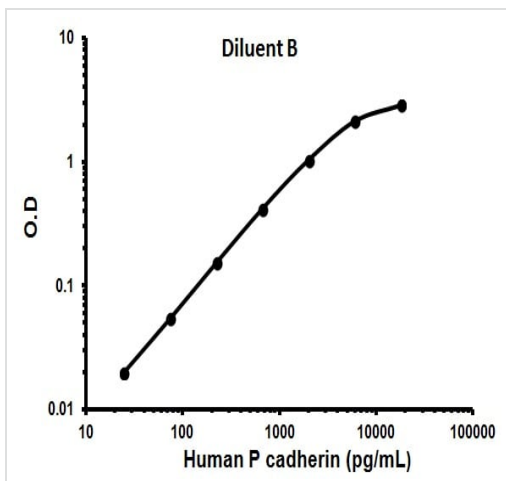
Cellular localization Cell membrane.

Images



Representative Standard Curve using ab100621.

Typical Standard Curve



Representative Standard Curve using ab100621.

Typical 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