

Human OB Cadherin ELISA Kit ab267594

1 Image

Overview

Product name Human OB Cadherin ELISA Kit

Detection method Colorimetric

Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 10%

	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				< 12%

Sample type Serum, Plasma, Cell culture media

Assay type Sandwich (quantitative)

Sensitivity 0.8 ng/ml

Range 0.819 ng/ml - 200 ng/ml

Recovery	Sample specific recovery		
	Sample type	Average %	Range
	Serum	93.61	87% - 100%
	Plasma	87	79% - 91%
	Cell culture media	114	102% - 128%

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Human

Product overview Human OB Cadherin ELISA Kit is designed for the quantitative determination of OB Cadherin in cell culture supernatants, plasma and serum samples.

This assay employs an antibody specific for human OB Cadherin coated on a 96-well plate. Standards and samples are pipetted into the wells and OB Cadherin present in a sample is

bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-human OB 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 OB Cadherin bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

This ELISA kit detects human OB Cadherin, no cross-reactivity with Cadherin-4, 6, 8, 12, 13, 17, E-Cadherin, N-Cadherin, P-Cadherin, or VE-Cadherin is observed.

Platform Pre-coated microplate (12 x 8 well strips)

Properties

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

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
5X Assay Diluent D	1 x 15ml
Anti-Human OB Cadherin coated Microplate (12 x 8 wells)	1 unit
Biotinylated Anti-Human OB Cadherin Detection Antibody	2 vials
Human OB Cadherin Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	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.

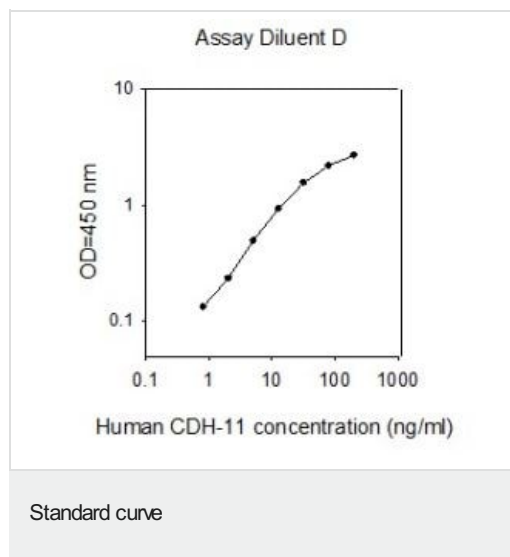
Tissue specificity Expressed mainly in brain but also found in other tissues. Expressed in neuroblasts.

Involvement in disease Note=A chromosomal aberration involving CDH11 is a common genetic feature of aneurysmal bone cyst, a benign osseous neoplasm. Translocation t(16;17)(q22;p13) with USP6. The translocation generates a fusion gene in which the strong CDH11 promoter is fused to the entire USP6 coding sequence, resulting in USP6 transcriptional up-regulation.

Sequence similarities Contains 5 cadherin domains.

Cellular localization Cell membrane.

Images



Human OB Cadherin ELISA kit (ab267594) Standard curve.

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.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