

Product datasheet

Mouse OB Cadherin ELISA Kit ab208346

Recombinant SimpleStep ELISA

[5 Images](#)

Overview

Product name Mouse OB Cadherin ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
serum	5			3.5%

Inter-assay

Sample	n	Mean	SD	CV%
serum	3			14.2%

Sample type

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

Assay type

Sandwich (quantitative)

Sensitivity

7 pg/ml

Range

46.9 pg/ml - 3000 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	111	109% - 115%
Cell culture media	87	76% - 94%
Hep Plasma	116	112% - 122%
EDTA Plasma	114	106% - 125%
Cit plasma	103	101% - 109%

Assay time

1h 30m

Assay duration One step assay

Species reactivity **Reacts with:** Mouse
Does not react with: Cow, Pig

Product overview Mouse OB Cadherin ELISA Kit (ab208346) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of OB Cadherin protein in cell culture extracts, cell culture supernatant, serum, tissue extracts, urine, and plasma. It uses our proprietary SimpleStep ELISA® technology. Quantitate Mouse OB Cadherin with 7 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 OB Cadherin (also known as Cadherin-11) is a member of the Cadherin family of calcium-dependent cell adhesion proteins. OB Cadherin is a single-pass transmembrane protein on the plasma membrane. The extracellular portion contains 5 cadherin domains that bind calcium and form homodimeric interactions.

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 OB Cadherin Capture Antibody	1 x 600µl
10X Mouse OB Cadherin 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

Components	1 x 96 tests
Antibody Diluent 5BR	1 x 6ml
Mouse OB Cadherin Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
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

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.

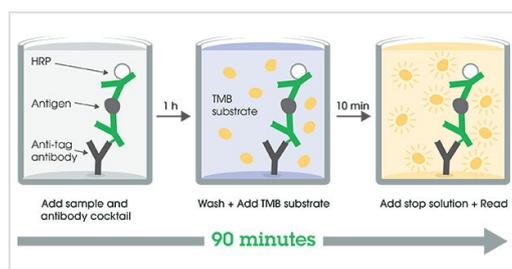
Sequence similarities

Contains 5 cadherin domains.

Cellular localization

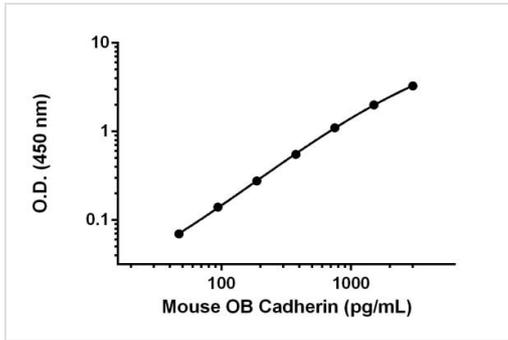
Cell membrane.

Images



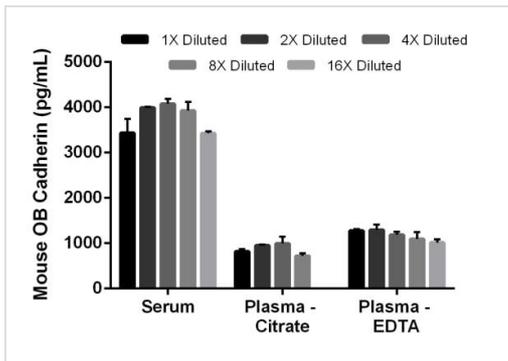
Other - Mouse OB Cadherin ELISA Kit (ab208346)

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.



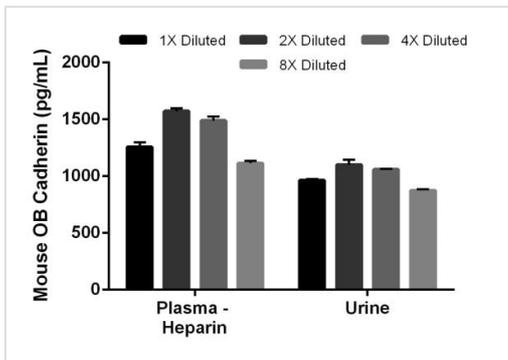
Example of mouse OB Cadherin standard curve in Sample Diluent NS

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



Interpolated concentrations of native OB-Cadherin in mouse serum, plasma (citrate) and plasma (EDTA) samples

The concentrations of OB Cadherin were measured in duplicates, interpolated from the OB Cadherin standard curves and corrected for sample dilution. 1X diluted samples are as follows: serum 25%, plasma (citrate) 25%, and plasma (EDTA) 25%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean OB Cadherin concentration was determined to be 3,771 pg/mL in serum, 827 pg/mL in plasma (citrate) and 1,172 pg/mL in plasma (EDTA).



Interpolated concentrations of native OB-Cadherin in mouse plasma (heparin) and urine samples

The concentrations of OB Cadherin were measured in duplicates, interpolated from the OB Cadherin standard curves and corrected for sample dilution. 1X diluted samples are as follows: plasma (heparin) 25%, and urine 20%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean OB Cadherin concentration was determined to be 1,343 pg/mL in plasma (heparin) and 1,000 pg/mL in urine.

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 - Mouse OB Cadherin ELISA Kit
(ab208346)

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