

Mouse E-Cadherin ELISA Kit (CDH1) ab100678

1 References 2 Images

Overview

Product name	Mouse E-Cadherin ELISA Kit (CDH1)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 6 pg/ml
Range	4.12 pg/ml - 3000 pg/ml
Recovery	94 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	114.7	103% - 123%
Serum	78.78	69% - 88%
Plasma	88.93	76% - 95%

Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Mouse
Product overview	

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

This assay employs an antibody specific for mouse E-Cadherin coated on a 96- well plate. Standards and samples are pipetted into the wells and E-Cadherin present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse E-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 E-Cadherin bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Get higher sensitivity in only 90 minutes with Mouse E-Cadherin ELISA Kit ([ab197751](#)) from our

SimpleStep ELISA® range.

Platform

Microplate

Properties

Storage instructions

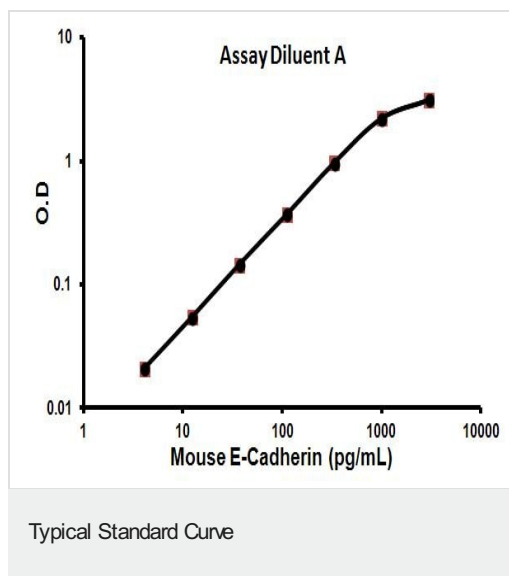
Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Mouse E-Cadherin	2 vials
E Cadherin Microplate (12 x 8 wells)	1 unit
Recombinant Mouse E-Cadherin Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

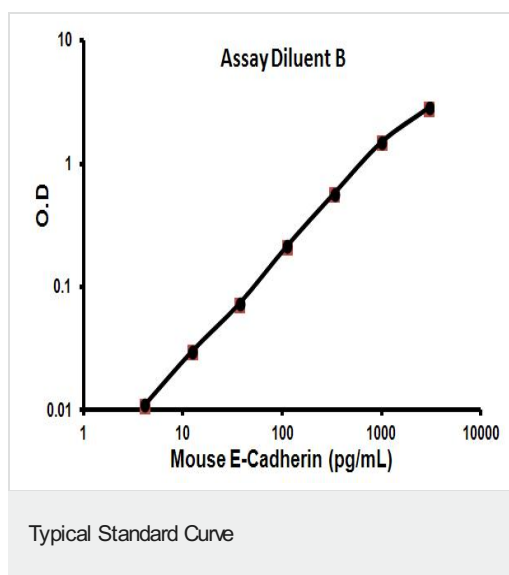
Relevance

One of the epithelial cell adhesion molecules, E Cadherin, plays an important role in the formation of cell-cell contacts in epithelia irrespective their origin from ecto-, meso- or endodermal tissue. This early adhesion event between epithelial cells is a prerequisite for the assembly of all elements of the junctional complex. Furthermore, E Cadherin plays a crucial role in the maintenance of the epithelial junctional complex and is as such an important molecule in maintaining epithelial integrity. Over 90% of the malignant tumors are carcinomas. One of the prerequisites for the release of carcinoma cells from the primary site might be a defect in intercellular adhesion mediated by the absence of E Cadherin expression. Therefore, the expression of E Cadherin might be an important parameter for the determination of the invasive potential of epithelial neoplasms, and for the transition of a benign to a malignant neoplasm.

Images



Representative Standard Curve using [ab200678](#)



Representative Standard Curve using [ab200678](#)

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