

Mouse DLL1 ELISA Kit ab208342

Recombinant SimpleStep ELISA[®]

[5 Images](#)

Overview

Product name Mouse DLL1 ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Mouse serum	5			2.8%

Inter-assay

Sample	n	Mean	SD	CV%
Mouse serum	3			7.2%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity 24 pg/ml

Range 62.5 pg/ml - 4000 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	99	94% - 104%
Serum	100	99% - 101%
Hep Plasma	101	98% - 104%
EDTA Plasma	103	102% - 105%
Cit plasma	104	103% - 106%

Assay time 1h 30m

Assay duration One step assay

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

Mouse DLL1 ELISA Kit (ab208342) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of DLL1 protein in cell culture supernatant, serum, and plasma. It uses our proprietary SimpleStep ELISA® technology. Quantitate Mouse DLL1 with 24 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

DLL1 (Delta-like protein 1) is a transmembrane glycoprotein that is needed during development for cell-to-cell communication in embryos and in the nervous system development. In adult tissues, DLL1 is involved in morphogenesis and remodeling and is overexpressed in several types of cancers.

Platform

Pre-coated microplate (12 x 8 well strips)

Properties**Storage instructions**

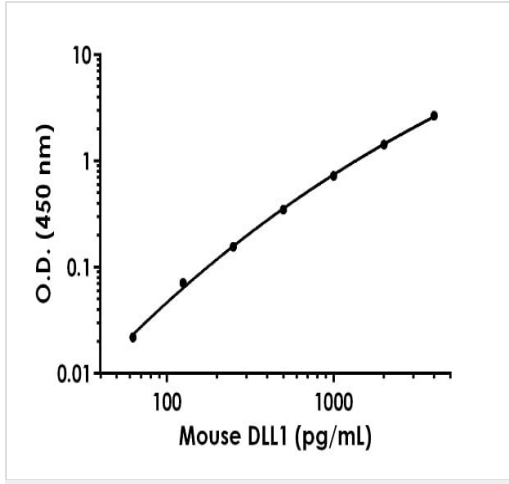
Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X Mouse DLL1 Capture Antibody	1 x 600µl
10X Mouse DLL1 Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 5BR	1 x 6ml
Mouse DLL1 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

Components	1 x 96 tests
TMB Development Solution	1 x 12ml

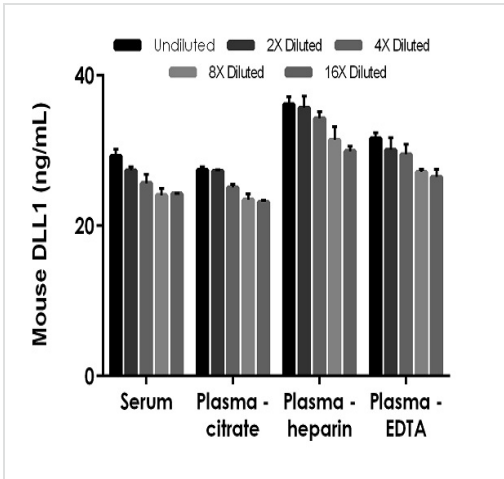
Function	Acts as a ligand for Notch receptors. Blocks the differentiation of progenitor cells into the B-cell lineage while promoting the emergence of a population of cells with the characteristics of a T-cell/NK-cell precursor.
Tissue specificity	Expressed in heart and pancreas, with lower expression in brain and muscle and almost no expression in placenta, lung, liver and kidney.
Sequence similarities	Contains 1 DSL domain. Contains 8 EGF-like domains.
Post-translational modifications	Ubiquitinated by MIB (MIB1 or MIB2), leading to its endocytosis and subsequent degradation.
Cellular localization	Membrane.

Images



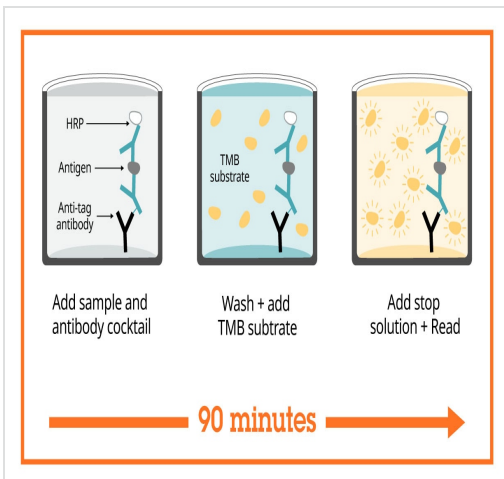
Background-subtracted data values (mean +/- SD) are graphed.

Example of mouse DLL1 standard



Interpolated concentrations of native DLL1 in (mouse) serum and plasma samples.

The concentrations of DLL1 were measured in duplicates, interpolated from the DLL1 standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 5%, plasma (citrate) 5%, plasma (heparin) 2.5% and spleen supernatant 25%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean DLL1 concentration was determined to be 26 ng/mL in serum, 26 ng/mL in plasma (citrate), 34 ng/mL in plasma (heparin) and 29 ng/mL in plasma (EDTA).



Sandwich ELISA - Mouse DLL1 ELISA Kit (ab208342)

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.

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 DLL1 ELISA Kit (ab208342)

To learn more about the advantages of recombinant antibodies see [here](#).

Get more done with SimpleStep ELISA



Easy to use
Single-wash 90-minute protocol



Flexible
Matched antibody pairs available



Precision antibodies
High sensitivity, specificity and reproducibility



Scalable
Now in 10-pack and 384-well formats

Sandwich ELISA - Mouse DLL1 ELISA Kit
(ab208342)

To learn more about the advantages of SimpleStep ELISA[®] kits 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