

## Product datasheet

# Mouse CD27 ELISA Kit ab233619

Recombinant **SimpleStep ELISA**

[5 Images](#)

### Overview

**Product name** Mouse CD27 ELISA Kit

**Detection method** Colorimetric

#### Precision

Intra-assay

Sample	n	Mean	SD	CV%
Serum	5			5.3%

Inter-assay

Sample	n	Mean	SD	CV%
Serum	3			2%

**Sample type** Cell culture supernatant, Serum, Hep Plasma, EDTA Plasma, Cit plasma

**Assay type** Sandwich (quantitative)

**Sensitivity** 10.9 pg/ml

**Range** 31.3 pg/ml - 2000 pg/ml

#### Recovery

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	104	98% - 108%
Serum	100	93% - 106%
Hep Plasma	105	99% - 109%
EDTA Plasma	90	79% - 99%
Cit plasma	94	88% - 98%

**Assay time** 1h 30m

**Assay duration** One step assay

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

**Product overview** Mouse CD27 ELISA Kit (ab233619) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of CD27 protein in cell culture supernatant, cit plasma, edta plasma, hep plasma, and serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Mouse CD27 with 10.9 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.

**Platform** Microplate (12 x 8 well strips)

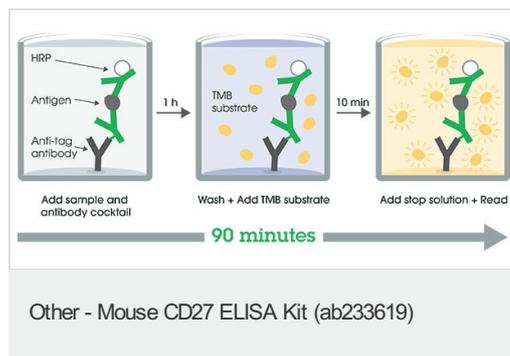
## Properties

**Storage instructions** Store at +4°C. Please refer to protocols.

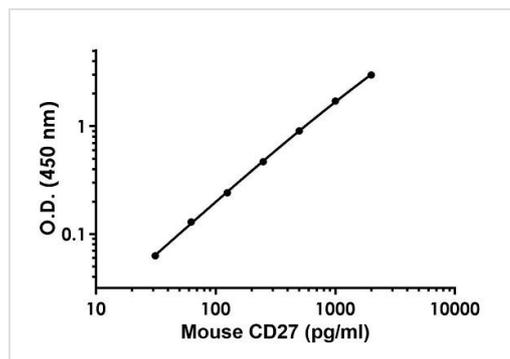
Components	1 x 96 tests
10X Mouse CD27 Capture Antibody	1 x 600µl
10X Mouse CD27 Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 4BR	1 x 6ml
Mouse CD27 Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

<b>Function</b>	Receptor for CD70/CD27L. May play a role in survival of activated T-cells. May play a role in apoptosis through association with SIVA1.
<b>Tissue specificity</b>	Found in most T-lymphocytes.
<b>Sequence similarities</b>	Contains 3 TNFR-Cys repeats.
<b>Post-translational modifications</b>	Phosphorylated and O-glycosylated.
<b>Cellular localization</b>	Membrane.

## Images

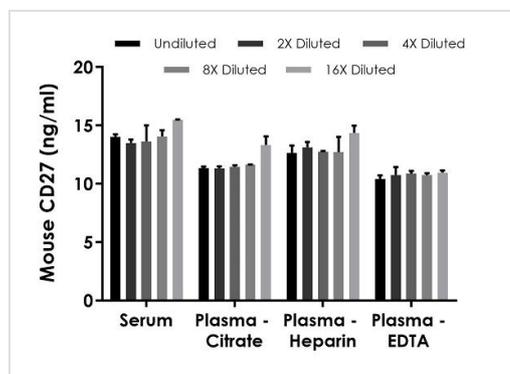


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



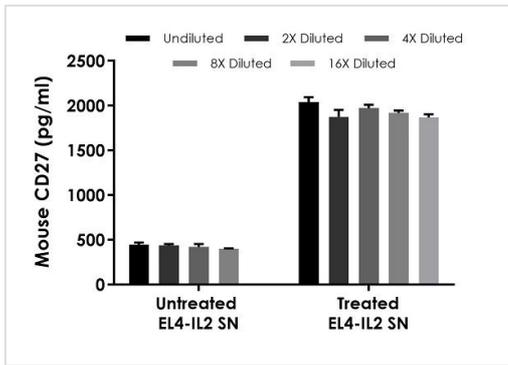
Background-subtracted data values (mean +/- SD) are graphed.

Example of mouse CD27 standard curve in Sample Diluent NS



The concentrations of CD27 were measured in duplicates, interpolated from the CD27 standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 12.5%, plasma (citrate) 12.5%, plasma (heparin) 12.5%, and plasma (EDTA) 12.5%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean CD27 concentration was determined to be 14.3 ng/mL in serum, 11.9 ng/mL in plasma (citrate), 13.4 g/ml in plasma (heparin), and 10.8 ng/ml in plasma (EDTA).

Interpolated concentrations of native CD27 in mouse serum and plasma



Interpolated concentrations of native CD27 in mouse EL4-IL2 cell culture supernatant samples, both untreated and treated with 10 µg/ml PHA and 10 ng/ml PMA for two days

The concentrations of CD27 were measured in duplicates, interpolated from the CD27 standard curves and corrected for sample dilution. Undiluted samples are as follows: untreated EL4-IL2 supernatant 100% and treated EL4-IL2 supernatant 50%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean CD27 concentration was determined to be 426 pg/mL in untreated EL4-IL2 supernatant and 1,978 pg/ml in treated EL4-IL2 supernatant.

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 CD27 ELISA Kit  
(ab233619)

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