

## Product datasheet

# Human CD85d ELISA Kit (LILRB2) ab269551

Recombinant SimpleStep ELISA

[4 Images](#)

### Overview

**Product name** Human CD85d ELISA Kit (LILRB2)

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Serum	8			2.1%

Inter-assay

Sample	n	Mean	SD	CV%
Serum	3			5.6%

**Sample type** Serum, Hep Plasma, EDTA Plasma, Cit plasma

**Assay type** Sandwich (quantitative)

**Sensitivity** 15.9 pg/ml

**Range** 62.5 pg/ml - 4000 pg/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
Serum	91	100% - 95%
Hep Plasma	84	93% - 89%
EDTA Plasma	102	94% - 107%
Cit plasma	95	100% - 94%

**Assay time** 1h 30m

**Assay duration** One step assay

**Species reactivity** **Reacts with:** Human

## Product overview

Human CD85d ELISA Kit (LILRB2) (ab269551) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of CD85d (LILRB2) protein in edta plasma, hep plasma, serum, and cit plasma. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human CD85d (LILRB2) with 15.9 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.

## Notes

CD85d (LILRB2), Leukocyte immunoglobulin-like receptor subfamily B member 2, is a key protein in the regulation of the immune response where it plays a pivotal as a negative regulator. It is primarily found on immune cells where it interacts with MHC domains on antigen presenting cells. Given its important role in immunosuppression, it is currently under investigation as a drug-targetable protein for a host of disease states. The antibodies for this kit were raised against a recombinant protein fragment targeting the extracellular domain of this receptor. There is little homology between mouse and rat for this fragment, so reactivity is not likely.

## Platform

Pre-coated microplate (12 x 8 well strips)

## Properties

### Storage instructions

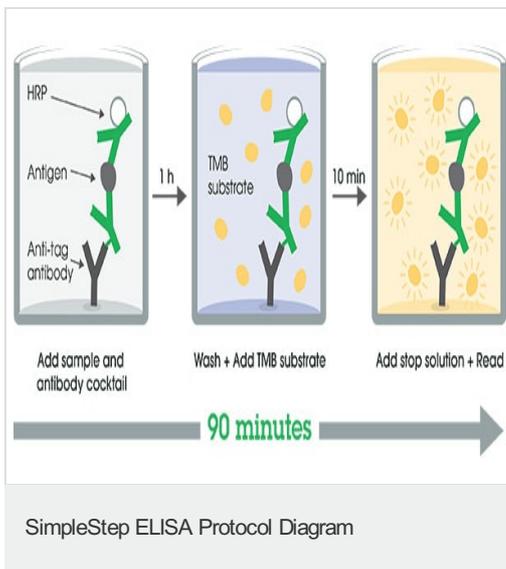
Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X Human CD85d (LILRB2) Capture Antibody	1 x 600µl
10X Human CD85d (LILRB2) Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent CPI - HAMA Blocker (ab193969)	1 x 6ml
Human CD85d (LILRB2) Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 12ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit

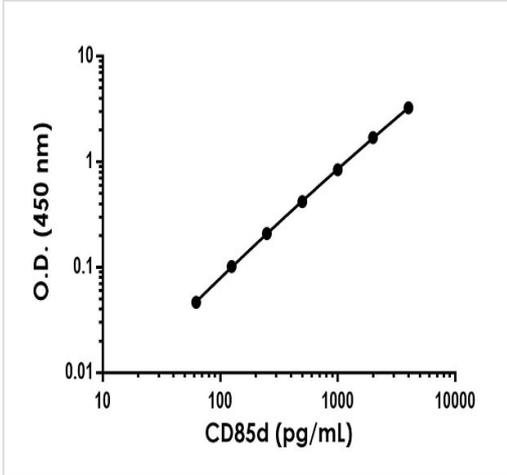
Components	1 x 96 tests
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

<b>Function</b>	Receptor for class I MHC antigens. Recognizes a broad spectrum of HLA-A, HLA-B, HLA-C and HLA-G alleles. Involved in the down-regulation of the immune response and the development of tolerance. Competes with CD8A for binding to class I MHC antigens. Inhibits FCGR1A-mediated phosphorylation of cellular proteins and mobilization of intracellular calcium ions.
<b>Tissue specificity</b>	Expressed on monocytes and B-cells, and at lower levels on dendritic cells. Detected at low levels in natural killer (NK) cells.
<b>Sequence similarities</b>	Contains 4 Ig-like C2-type (immunoglobulin-like) domains.
<b>Domain</b>	Contains 3 copies of a cytoplasmic motif that is referred to as the immunoreceptor tyrosine-based inhibitor motif (ITIM). This motif is involved in modulation of cellular responses. The phosphorylated ITIM motif can bind the SH2 domain of several SH2-containing phosphatases.
<b>Post-translational modifications</b>	Phosphorylated on tyrosine residues. Dephosphorylated by PTPN6.
<b>Cellular localization</b>	Membrane.

## Images

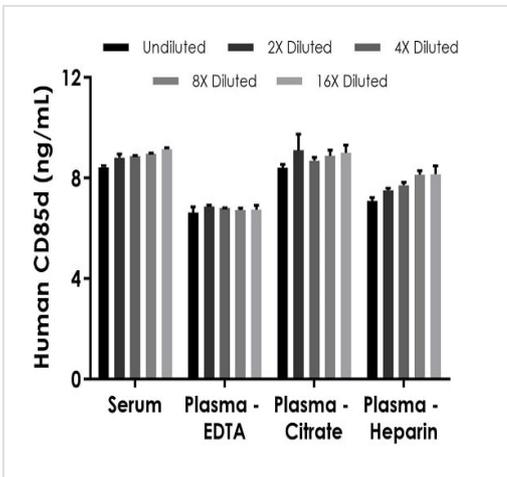


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



Example of human CD85d standard curve in Sample Diluent NS.

The CD85d standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



Interpolated concentrations of native CD85d in human serum and plasma samples.

The concentrations of CD85d were measured in duplicates, interpolated from the CD85d standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 30%, plasma (EDTA) 30%, plasma (citrate) 30%, and plasma (heparin) 30%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean CD85d concentration was determined to be 8.85 ng/mL in serum, 6.75 ng/mL in plasma (EDTA), 8.82 ng/mL in plasma (citrate), and 7.72 ng/mL in plasma (Heparin).

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 - Human CD85d ELISA Kit  
(LILRB2) (ab269551)

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