

Product datasheet

Human 4-1BBL ELISA Kit ab246536

Recombinant SimpleStep ELISA

[4 Images](#)

Overview

Product name Human 4-1BBL ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
supernatant	8			7%

Inter-assay

Sample	n	Mean	SD	CV%
supernatant	3			3.8%

Sample type Cell culture supernatant, Serum, Cit plasma

Assay type Sandwich (quantitative)

Sensitivity 2.2 pg/ml

Range 6.25 pg/ml - 400 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	101	99% - 104%
Serum	88	86% - 89%
Cit plasma	92	90% - 94%

Assay time 1h 30m

Assay duration One step assay

Species reactivity **Reacts with:** Human

Product overview Human 4-1BBL ELISA Kit (ab246536) is a single-wash 90 min sandwich ELISA designed for the

quantitative measurement of 4-1BBL protein in cell culture supernatant, cit plasma, and serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human 4-1BBL with 2.2 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 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 4-1BBL Capture Antibody	1 x 600µl
10X Human 4-1BBL Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 5BI	1 x 6ml
Human 4-1BBL 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 Cytokine that binds to TNFRSF9. Induces the proliferation of activated peripheral blood T-cells. May have a role in activation-induced cell death (AICD). May play a role in cognate interactions between T-cells and B-cells/macrophages.

Tissue specificity Expressed in brain, placenta, lung, skeletal muscle and kidney.

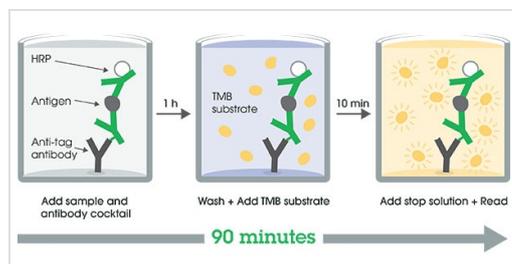
Sequence similarities

Belongs to the tumor necrosis factor family.

Cellular localization

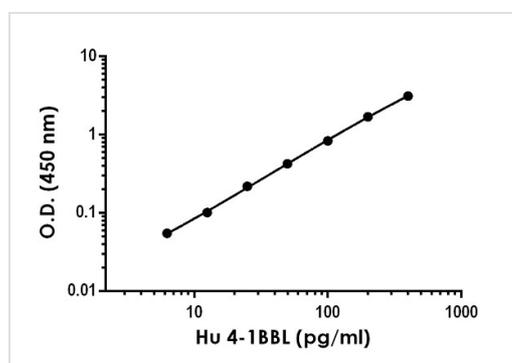
Membrane.

Images



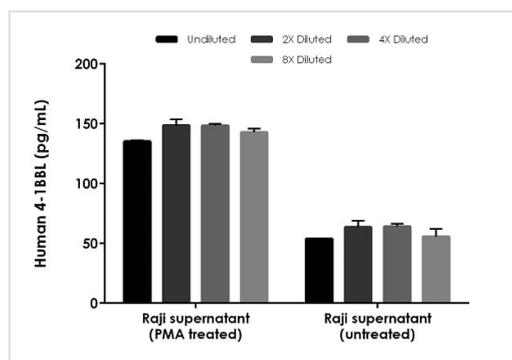
Other - Human 4-1BBL ELISA Kit (ab246536)

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 4-1BBL standard curve in Sample Diluent NS.

The 4-1BBL standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean \pm SD) are graphed.



Interpolated concentrations of native 4-1BBL in human cell culture supernatant samples.

The concentrations of 4-1BBL were measured in duplicates, interpolated from the 4-1BBL standard curves and corrected for sample dilution. Undiluted samples are as follows: Raji (PMA treated) cell culture supernatant 50% and Raji (untreated) cell culture supernatant 100%. The interpolated dilution factor corrected values are plotted (mean \pm SD, $n=2$). The mean 4-1BBL concentration was determined to be 143.7 pg/mL in Raji (PMA treated) cell culture supernatant, and 59.2 pg/mL in Raji (untreated) cell culture 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 - Human 4-1BBL ELISA Kit
(ab246536)

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