

Product datasheet

Mouse TARC ELISA Kit (CCL17) ab213474

Recombinant SimpleStep ELISA

[9 Images](#)

Overview

Product name Mouse TARC ELISA Kit (CCL17)

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall	5			9.6%

Inter-assay

Sample	n	Mean	SD	CV%
Overall	3			6.8%

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

Assay type Sandwich (quantitative)

Sensitivity 2.5 pg/ml

Range 7.8 pg/ml - 500 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	86	81% - 92%
Cell culture media	98	93% - 101%
Hep Plasma	81	76% - 84%
EDTA Plasma	107	103% - 112%
Cit plasma	101	95% - 107%

Assay time 1h 30m

Assay duration One step assay

Species reactivity **Reacts with:** Mouse

Product overview Mouse TARC ELISA Kit (CCL17) (ab213474) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of TARC (CCL17) protein in cell culture supernatant, cit plasma, hep plasma, serum, and edta plasma. It uses our proprietary SimpleStep ELISA® technology. Quantitate Mouse TARC (CCL17) with 2.5 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 TARC (also known as C-C motif chemokine 17) is a chemokine that functions through chemokine receptors CCR4 and CCR8. TARC is expressed in the thymus as well as by dendritic cells, keratinocytes and bronchial epithelial cells.

Platform Pre-coated microplate (12 x 8 well strips)

Properties

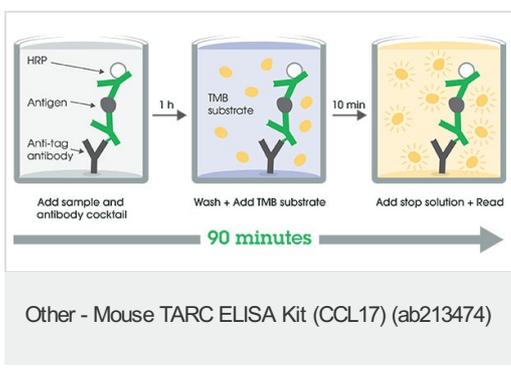
Storage instructions Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X Mouse TARC Capture Antibody	1 x 600µl
10X Mouse TARC Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 4BR	1 x 6ml
Mouse TARC Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Sample Diluent NBS	1 x 20ml

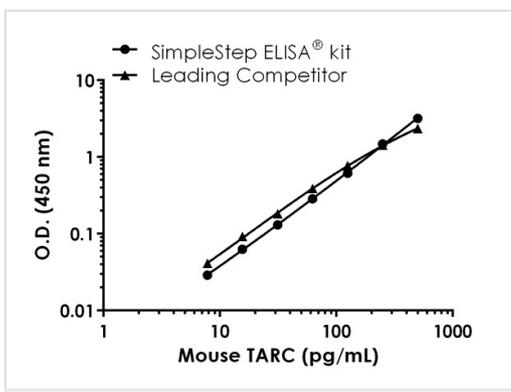
Components	1 x 96 tests
Sample Diluent NS (ab193972)	1 x 50ml
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

Function	Chemotactic factor for T-lymphocytes but not monocytes or granulocytes. May play a role in T-cell development in thymus and in trafficking and activation of mature T-cells. Binds to CCR4.
Tissue specificity	Expressed at high levels in thymus and at low levels in the lung, colon and small intestine.
Sequence similarities	Belongs to the intercrine beta (chemokine CC) family.
Cellular localization	Secreted.

Images

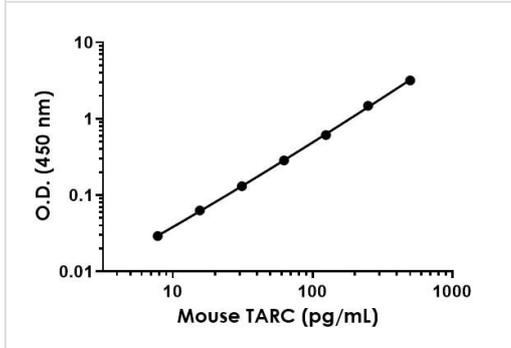


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



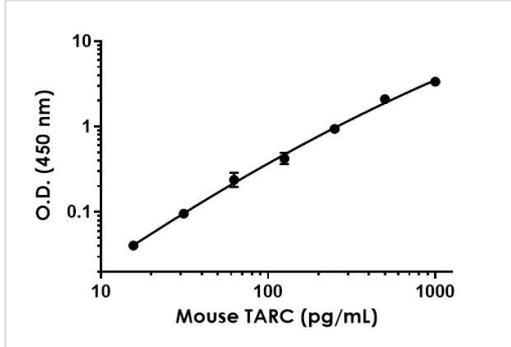
Standard curve comparison between mouse TARC SimpleStep ELISA[®] kit and traditional ELISA kit from leading competitor. SimpleStep ELISA kit shows comparable sensitivity.

Mouse TARC standard curve comparison data.



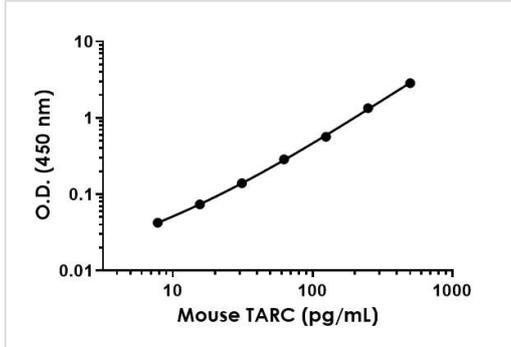
Background-subtracted data values (mean +/- SD) are graphed.

Example of mouse TARC standard curve in Sample Diluent NS.



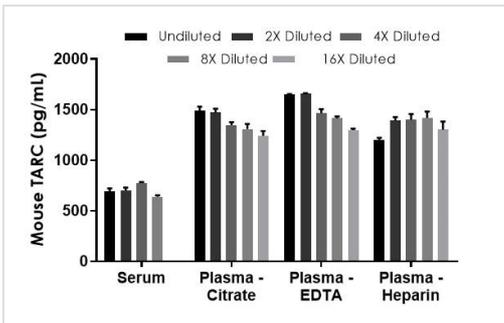
Background-subtracted data values (mean +/- SD) are graphed.

Example of mouse TARC standard curve in Sample Diluent 50BS.



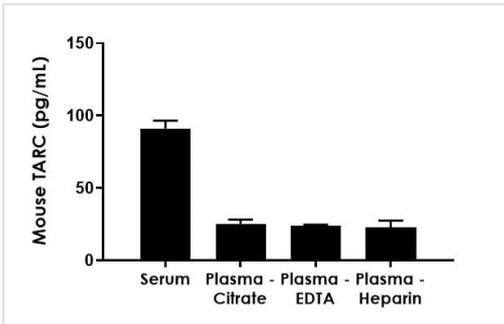
Background-subtracted data values (mean +/- SD) are graphed.

Example of mouse TARC standard curve in Sample Diluent 5BS.



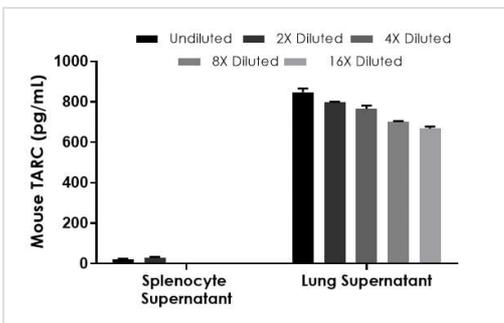
Interpolated concentrations of spiked TARC in mouse serum and plasma samples.

The concentrations of TARC were measured in duplicates, interpolated from the TARC standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 50%, plasma (citrate) 50%, plasma (EDTA) 50%, and plasma (heparin) 25%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2).



Interpolated concentrations of native TARC in mouse serum and plasma samples.

The concentrations of TARC were measured in duplicates, interpolated from the TARC standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 50%, plasma (citrate) 50%, plasma (EDTA) 50%, and plasma (heparin) 25%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean TARC concentration was determined to be 91 pg/mL in serum, 25.4 pg/mL in plasma (citrate), 23.0 pg/mL in plasma (EDTA) and 25.4 pg/mL in plasma (heparin).



Interpolated concentrations of native TARC in mouse splenocyte and lung supernatant samples.

The concentrations of TARC were measured in duplicates, interpolated from the TARC standard curves and corrected for sample dilution. Undiluted samples are as follows: splenocyte supernatant 50%, and lung supernatant 50%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean TARC concentration was determined to be 25 pg/mL in splenocyte supernatant and 756 pg/mL in lung 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 TARC ELISA Kit (CCL17)
(ab213474)

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