

Product datasheet

Mouse CCL21α Matched Antibody Pair Kit ab216060

1 Image

Overview

Product name	Mouse CCL21a Matched Antibody Pair Kit
Detection method	Colorimetric
Assay type	ELISA set
Sensitivity	2.9 pg/ml
Range	3.9 pg/ml - 250 pg/ml
Species reactivity	Reacts with: Mouse
Product overview	<p>Mouse CCL21a Matched Antibody Pair Kits include a capture and a biotinylated detector antibody pair, along with a calibrated protein standard, suitable for sandwich ELISA. The Matched Antibody Pair Kit can be used to quantify native and recombinant mouse CCL21a.</p>

Optimization of the kit reagents to sample type, immunoassay format or instrumentation may be required. Guidelines for use of this kit in a standard 96-well microplate sandwich ELISA using HRP/TMB system of colorimetric detection is described in this assay procedure for the purposes of quantification.

Protocol information and tips on the use of the Matched Antibody Pair kits for sandwich ELISA can be found on our [website](#). An accessory pack can be purchased which includes buffer reagents required to perform 10 x 96-well plate sandwich ELISAs ([ab210905](#)).

For additional information on the performance of the antibody pair used in this kit, please see our equivalent SimpleStep ELISA kit [ab208985](#). Please note that while the antibody pair is the same provided in the corresponding SimpleStep ELISA Kit, due to differences in their formulation, this antibody pair cannot be used with the consumables provided with our SimpleStep ELISA Kits.

Tested applications	Suitable for: ELISA
Platform	Reagents

Properties

Storage instructions	Store at -20°C. Please refer to protocols.
-----------------------------	--

Components	10 x 96 tests	5 x 96 tests
Mouse CCL21a Capture Antibody	2 x 50µg	1 x 50µg
Mouse CCL21a Detector Antibody	2 x 12.5µg	1 x 12.5µg
Mouse CCL21a Lyophilized Protein	2 vials	1 vial

Function	Inhibits hemopoiesis and stimulates chemotaxis. Chemotactic in vitro for thymocytes and activated T-cells, but not for B-cells, macrophages, or neutrophils. Potent mesangial cell chemoattractant. Shows preferential activity towards naive T-cells. May play a role in mediating homing of lymphocytes to secondary lymphoid organs.
Tissue specificity	Expressed strongly in lung, spleen, thymus, peripheral and mesenteric lymph nodes. Also expressed in the testis, kidney, liver, and heart.
Involvement in disease	Note=Mice carrying an autosomal recessive mutation designated paucity of lymph node T-cells (plt) show dramatically reduced numbers of T-cells in lymph nodes, Peyer patches, and the white pulp of the spleen. Plt seems to correspond to Scya21b.
Sequence similarities	Belongs to the intercrine beta (chemokine CC) family.
Cellular localization	Secreted.

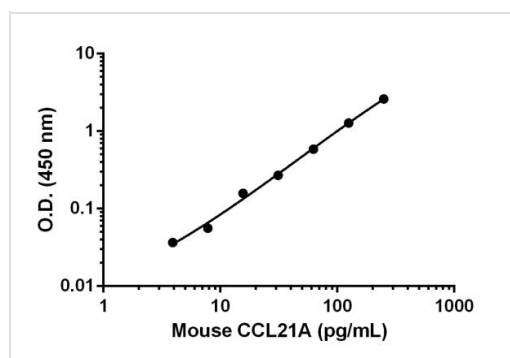
Applications

Our [Abpromise guarantee](#) covers the use of **ab216060** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
ELISA		Use at an assay dependent concentration.

Images



Mouse CCL21a standard curve.

Standard calibration curve. Background subtracted values are graphed.

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