

Product datasheet

Mouse IL-17C ELISA Kit ab213868

1 Image

Overview

Product name Mouse IL-17C ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
1	16	150pg/ml	11.85	= 7.9%
2	16	1098pg/ml	68.07	= 6.2%
3	16	6281pg/ml	414.54	= 6.6%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	140pg/ml	13.58	= 9.7%
2	24	1153pg/ml	85.32	= 7.4%
3	24	5857pg/ml	447.63	= 7.6%

Sample type Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma

Assay type Sandwich (quantitative)

Sensitivity < 10 pg/ml

Range 125 pg/ml - 8000 pg/ml

Assay time 3h 30m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview The Mouse IL-17C Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213868) is designed for the quantitative measurement of Mouse IL-17C in cell culture supernatants, cell lysates, serum and plasma (heparin, EDTA).

The ELISA kit is based on standard sandwich enzyme-linked immune-sorbent assay technology.

A monoclonal antibody from rat specific for IL-17C has been pre-coated onto 96-well plates. Standards (Expression system for standard: E.coli; Immunogen sequence: D17-Q194) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for IL-17C is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. HRP substrate TMB is used to visualize HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the Mouse IL-17C amount of sample captured in plate.

Notes IL-17C, also known as CX2, is a protein that in humans is encoded by the IL-17C gene. IL-17C is mapped to 16q24.3. The protein encoded by this gene is a T cell-derived cytokine that shares the sequence similarity with IL-17. This cytokine was reported to stimulate the release of tumor necrosis factor alpha and interleukin 1 beta from a monocytic cell line. The expression of this cytokine was found to be restricted to activated T cells. IL-17C is an essential autocrine cytokine regulating innate epithelial immune responses. It also plays an important role in the pathogenesis of inflammatory arthritis.

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

Properties

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

Components	Identifier	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Adhesive Plate Seal		4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-Mouse IL-17C coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti- Mouse IL-17C antibody		1 x 100µl
Lyophilized recombinant Mouse IL-17C standard		2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml

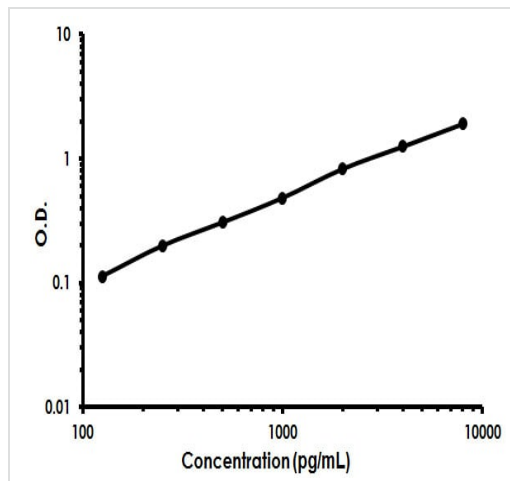
Function Stimulates the release of tumor necrosis factor alpha and IL-1-beta from the monocytic cell line THP-1.

Sequence similarities Belongs to the IL-17 family.

Cellular localization Secreted.

Images

Mouse IL-17C ELISA Kit (ab213868) Standard Curve.



Mouse IL-17C ELISA Kit (ab213868) Standard Curve.

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