

Mouse IL-17RC ELISA Kit ab213870

1 Image

Overview

Product name Mouse IL-17RC ELISA Kit

Detection method Colorimetric

Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	1	16	184pg/ml	11.77	= 6.4%
	2	16	1503pg/ml	60.12	= 4%
	3	16	5668pg/ml	345.74	= 6.1%

	Inter-assay				
	Sample	n	Mean	SD	CV%
	1	24	190pg/ml	13.68	= 7.2%
	2	24	1560pg/ml	65.52	= 4.2%
	3	24	5820pg/ml	413.22	= 7.1%

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

Assay type Sandwich (quantitative)

Sensitivity < 10 pg/ml

Range 156 pg/ml - 10000 pg/ml

Assay time 3h 30m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview The Mouse IL-17RC Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213870) is designed for the quantitative measurement of Mouse IL-17RC 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-17RC has been pre-coated onto 96-well plates. Standards (Expression system for standard: NSO; Immunogen sequence: E22-R440) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for IL-17RC 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-17RC amount of sample captured in plate.

**Notes** Interleukin-17 receptor C is a protein that in humans is encoded by the IL-17RC gene. By genomic sequence analysis, IL-17RC gene was mapped to chromosome 3p25.3-p24.1. This gene encodes a single-pass transmembrane protein that shares limited similarity with the interleukin-17 receptor. Multiple alternatively spliced transcript variants encoding different isoforms have been detected for this gene, but the full-length nature of only three have been determined to date. The biologic activity of IL-17 is dependent on a complex composed of IL-17RA and IL-17RC, suggesting a new paradigm for understanding the interactions between the expanded family of IL-17 ligands and their receptors.

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

## Properties

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

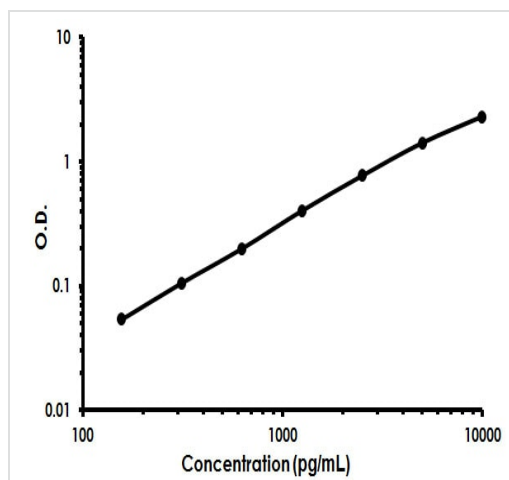
Components	Identifier	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Adhesive Plate Seal		4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-Mouse IL-17RC coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti- Mouse IL-17RC antibody		1 x 100µl
Lyophilized recombinant Mouse IL-17RC 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
Wash Buffer (25X)		1 x 20ml

**Relevance** IL-17RC is a single pass transmembrane protein that shares limited similarity with the interleukin 17 receptor. It is widely expressed in non leukocyte tissues and forms a heteromeric receptor for IL-17 with IL-17RA. At least four different isoforms are known to exist.

**Cellular localization** Cell membrane; Single pass type I membrane protein. Note: soluble isoforms may be produced.

## Images

---



Mouse IL-17RC ELISA Kit (ab213870) Standard Curve.

Mouse IL-17RC ELISA Kit (ab213870) 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