

Mouse IL-17D ELISA Kit ab213869

1 Image

Overview

Product name	Mouse IL-17D ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	1	16	715pg/ml	30.03	4.2%
	2	16	1410pg/ml	64.86	4.6%
	3	16	2703ng/ml	137.9	5.1%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	1	24	694pg/ml	38.17	5.5%
	2	24	1582pg/ml	99.67	6.3%
	3	24	2691pg/ml	180.3	6.7%
Sample type	Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma, Tissue Homogenate				
Assay type	Sandwich (quantitative)				
Sensitivity	< 10 pg/ml				
Range	62.5 pg/ml - 4000 pg/ml				
Assay time	3h 30m				
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Mouse				
Product overview	The Mouse IL-17D Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213869) is designed for the quantitative measurement of Mouse IL-17D in cell culture supernatants, cell lysates, tissue homogenates, 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-17D has been pre-coated onto 96-well plates. Standards (Expression system for standard: E.coli; Immunogen sequence: A25-R205) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for IL-17D 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-17D amount of sample captured in plate.

Notes IL-17D is a member of the unique IL-17 family, which share a conserved cysteine motif. It is mapped to 13q12.11. IL-17D regulates cytokine production in endothelial cells and inhibits hemopoiesis *in vitro*. Purified recombinant IL-17D inhibited myeloid progenitor cell colony formation. IL-17D can enhance endothelial cell production of IL6, IL8, and GM-CSF, but not other cytokines, and it is preferentially expressed in skeletal muscle, brain, adipose tissue, heart, lung, and pancreas.

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

Properties

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

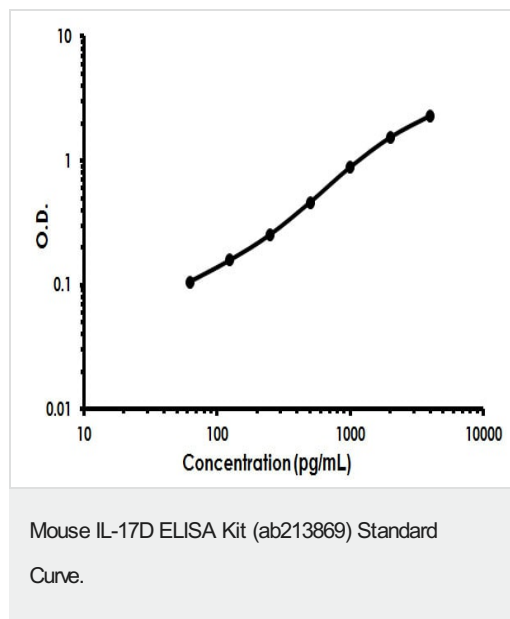
Components	Identifier	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Adhesive Plate Seal		4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-Mouse IL-17D coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti- Mouse IL-17D antibody		1 x 100µl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

Function Induces expression of IL-6, IL-8, and GM-CSF from endothelial cells.

Tissue specificity Expressed preferentially in adipose, skeletal muscle and CNS.

Sequence similarities Belongs to the IL-17 family.

Cellular localization Secreted.



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