

Mouse IL-31 ELISA Kit ab213872

[1 Image](#)

Overview

Product name Mouse IL-31 ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
1	16	387pg/ml	21.3	5.5%
2	16	1228pg/ml	78.6	6.4%
3	16	2155pg/ml	99.13	4.6%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	453pg/ml	31.26	6.9%
2	24	1396pg/ml	100.5	7.2%
3	24	2370pg/ml	128	5.4%

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

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-31 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213872) is designed for the quantitative measurement of Mouse IL-31 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-31 has been pre-coated onto 96-well plates. Standards (Expression system for standard: E.coli; Immunogen sequence: T24-C163) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for IL-31 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-31 amount of sample captured in plate.

Notes Interleukin-31, also known as IL-31 is a protein that in humans is encoded by the IL31 gene. The IL-31 gene is mapped to chromosome 12q24.31 and the mouse gene maps to a syntenic region on chromosome 5. IL-31, which is made principally by activated Th2-type T cells, interacts with a heterodimeric receptor consisting of IL-31RA and OSMR that is constitutively expressed on epithelial cells and keratinocytes. IL-31 may be involved in the promotion of allergic skin disorders and in regulating other allergic diseases, such as asthma.

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

Properties

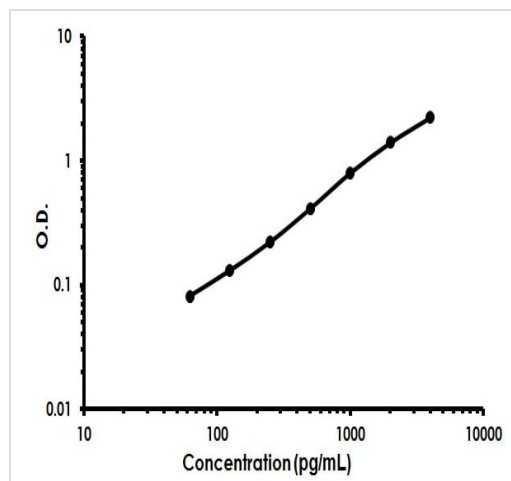
Storage instructions Store at -20°C. Please refer to protocols.

Components	Identifier	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Adhesive Plate Seal		4 units	4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-Mouse IL-31 coated Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti- Mouse IL-31 antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant Mouse IL-31 standard		2 vials	2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

Function Activates STAT3 and possibly STAT1 and STAT5 through the IL31 heterodimeric receptor composed of IL31RA and OSMR. IL31 may function in skin immunity.

Tissue specificity Detected at low levels in testis, bone marrow, skeletal muscle, kidney, colon, thymus, small intestine and trachea.

Cellular localization Secreted.



Mouse IL-31 ELISA Kit (ab213872) Standard Curve.

Mouse IL-31 ELISA Kit (ab213872) 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