

Product datasheet

Rat IL-17A ELISA Kit ab119536

[1 References](#) [1 Image](#)

Overview

Product name Rat IL-17A ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall	7			8.5%

Inter-assay

Sample	n	Mean	SD	CV%
Overall	7			7.6%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity 1 pg/ml

Range 1.6 pg/ml - 100 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant		92% - 112%
Serum		34% - 44%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Rat

Product overview

Abcam's IL-17A Rat *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for accurate quantitative measurement of Rat IL-17 (Interleukin-17) concentrations in Cell culture supernatant, serum and plasma (EDTA, citrate, heparin).

IL-17A specific antibodies have been precoated onto 96-well plates. Standards and test samples are added to the wells along with a biotin-conjugated IL-17A detection antibody then incubated at

room temperature. Following washing, a Streptavidin-HRP conjugate is added to each well, incubated at room temperature and washed. TMB is added and then catalyzed by HRP to produce a blue color product that changes into yellow after addition of acidic stop solution. The density of yellow coloration is directly proportional to the amount of IL-17A captured on the plate.

Tested applications

Suitable for: Sandwich ELISA

Platform

Microplate

Properties

Storage instructions

Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
20X Assay Buffer Concentrate	1 x 5ml
20X Wash Buffer Concentrate	1 x 50ml
Adhesive Films	4 units
Biotin-Conjugate anti-rat IL-17A monoclonal antibody	1 x 70µl
Microplate coated with monoclonal antibody to rat IL-17A (12 x 8 wells)	1 unit
Rat IL-17A Standard Lyophilized (200 pg/mL on reconstitution)	2 vials
Sample Diluent	1 x 12ml
Stop Solution (1M Phosphoric acid)	1 x 15ml
Streptavidin-HRP	1 x 150µl
TMB Substrate Solution	1 x 15ml

Function

Induces stromal cells to produce proinflammatory and hematopoietic cytokines. Enhances the surface expression of the intracellular adhesion molecule-1 (ICAM-1) in fibroblasts.

Tissue specificity

Restricted to activated memory T-cells.

Sequence similarities

Belongs to the IL-17 family.

Post-translational modifications

Found both in glycosylated and nonglycosylated forms.

Cellular localization

Secreted.

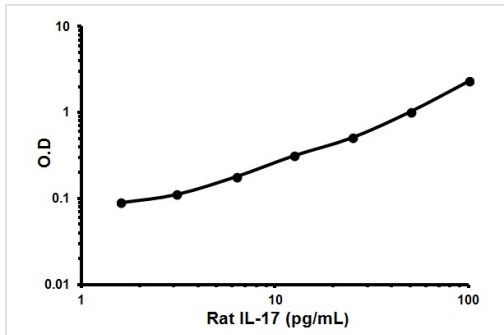
Applications

Our [Abpromise guarantee](#) covers the use of **ab119536** 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
Sandwich ELISA		Use at an assay dependent concentration.

Images



Representative Standard Curve using ab119536.

Typical 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