

## Product datasheet

# Monkey IL-6 ELISA Kit ab119549

1 Image

### Overview

**Product name** Monkey IL-6 ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Overall	6			< 5%

Inter-assay

Sample	n	Mean	SD	CV%
Overall	6			< 10%

**Sample type** Cell culture supernatant, Serum

**Assay type** Sandwich (quantitative)

**Sensitivity** 1.3 U/ml

**Range** 1.9 U/ml - 120 U/ml

**Recovery** 90 %

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Monkey

**Product overview** Abcam's IL-6 Monkey ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Monkey IL-6 in culture supernatant and serum (baboon, cynomolgus).

IL-6 specific antibodies have been precoated onto 96-well plates. Standards and test samples are added to the wells along with a biotin-conjugated IL-6 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-6 captured on the plate.

**Notes** Standard is recombinant protein produced in *E.coli*.

**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-monkey IL-6 monoclonal antibody	1 x 70µl
IL-6 Standard lyophilized (240 U/mL upon reconstitution)	2 vials
Microplate coated with monoclonal antibody to Monkey IL-6 (12 x 8 wells)	1 unit
Stop Solution (1M Phosphoric acid)	1 x 15ml
Streptavidin-HRP	1 x 150µl
TMB Substrate Solution	1 x 15ml

**Function** Cytokine with a wide variety of biological functions. It is a potent inducer of the acute phase response. Plays an essential role in the final differentiation of B-cells into Ig-secreting cells. Involved in lymphocyte and monocyte differentiation. It induces myeloma and plasmacytoma growth and induces nerve cells differentiation. Acts on B-cells, T-cells, hepatocytes, hematopoietic progenitor cells and cells of the CNS. Also acts as a myokine. It is discharged into the bloodstream after muscle contraction and acts to increase the breakdown of fats and to improve insulin resistance.

**Involvement in disease** Genetic variations in IL6 are associated with susceptibility to rheumatoid arthritis systemic juvenile (RASJ) [MIM:604302]. An inflammatory articular disorder with systemic-onset beginning before the age of 16. It represents a subgroup of juvenile arthritis associated with severe extraarticular features and occasionally fatal complications. During active phases of the disorder, patients display a typical daily spiking fever, an evanescent macular rash, lymphadenopathy, hepatosplenomegaly, serositis, myalgia and arthritis.  
Note=A IL6 promoter polymorphism is associated with a lifetime risk of development of Kaposi sarcoma in HIV-infected men.

**Sequence similarities** Belongs to the IL-6 superfamily.

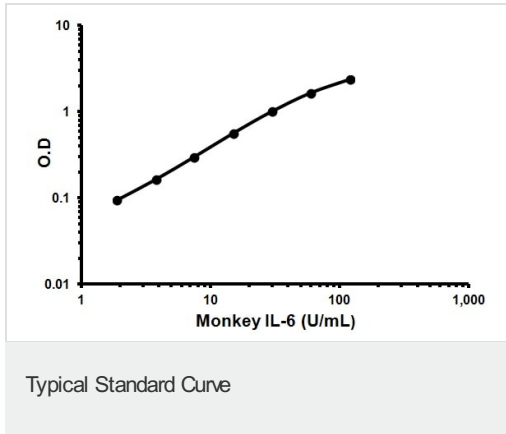
**Post-translational modifications** N- and O-glycosylated.

**Cellular localization** Secreted.

---

## Images

---



Representative Standard Curve using ab119549.

**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