

## Product datasheet

# Monkey IL-6 ELISA Kit (Interleukin-6) ab119549

### 画像数 1

#### 製品の概要

**製品名** Monkey IL-6 ELISA Kit (Interleukin-6)

**検出方法** Colorimetric

**再現性** Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
Overall	6			< 5%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Overall	6			< 10%

**サンプルの種類** Cell culture supernatant, Serum

**アッセイタイプ** Sandwich (quantitative)

**検出感度** 1.3 U/ml

**検出範囲** 1.9 U/ml - 120 U/ml

**添加回収試験** 90 %

**ステップ** Multiple steps standard assay

**種交差性** 交差種: Monkey

**製品の概要** Abcam's IL-6 Monkey (Interleukin-6) 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.

**特記事項** Standard is recombinant protein produced in *E.coli*.

**アプリケーション** 適用あり: Sandwich ELISA

## 製品の特性

## 保存方法

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

内容	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 IL6 monoclonal antibody	1 x 70µl
IL-6 Standard lyophilized (4 ng/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

## 機能

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.

## 関連疾患

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.

## 配列類似性

Belongs to the IL-6 superfamily.

## 翻訳後修飾

N- and O-glycosylated.

## 細胞内局在

Secreted.

## アプリケーション

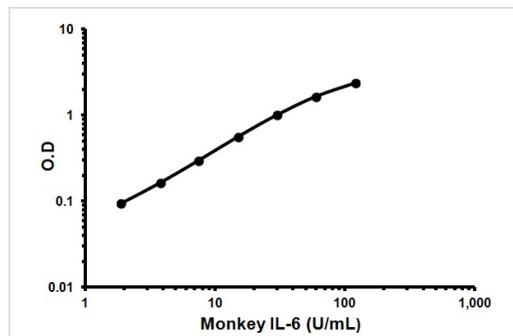
Our [Abpromise guarantee](#) covers the use of **ab119549** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Sandwich ELISA

Use at an assay dependent concentration.

## 画像



Representative Standard Curve using  
ab119549.

Typical Standard Curve

**Please note:** All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

### 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 <http://www.abcam.co.jp/abpromise> or contact our technical team.

### Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors