

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

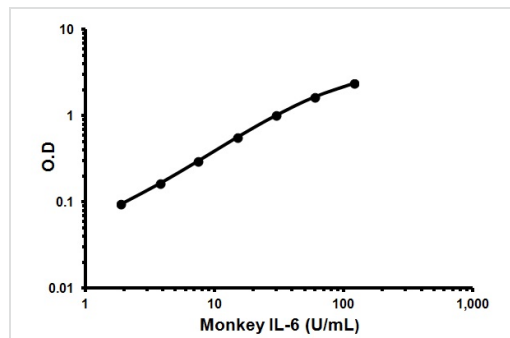
Our [Abpromise guarantee](#) covers the use of **ab119549** in the following tested applications.

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