

画像数 1

製品名	Mouse IL-6 ELISA Set (without plates)
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 10 pg/ml
検出範囲	15.6 pg/ml - 500 pg/ml
ステップ	Multiple steps standard assay
種交差性	<b>交差種:</b> Mouse
製品の概要	<p>Abcam's IL-6 Murine in vitro ELISA Set is designed for the quantitative measurement of Murine IL-6 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>This ELISA Set is intended to measure murine IL-6 in cell-culture supernatants, as well as in serum samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.</p> <p>The murine IL-6 ELISA Set includes both the capture and the detection antibodies as well as Horse-Radish peroxidase conjugated Streptavidin (HRP-Strep) and TMB which are the reagents used in the colour development step.</p> <p>Recognizes both natural and recombinant Murine IL6</p>
アプリケーション	<b>適用あり:</b> Sandwich ELISA
試験プラットフォーム	Reagents

## 製品の特長

**保存方法** Store at +4°C. Please refer to protocols.

内容	10 x 96 tests	15 x 96 tests	5 x 96 tests
Chromogen TMB Substrate Solution	4 vials	6 vials	2 vials
Detection biotinylated anti-murine IL-6 antibody	2 x 0.55ml	3 x 0.55ml	1 x 0.55ml

内容	10 x 96 tests	15 x 96 tests	5 x 96 tests
Murine IL-6 Capture antibody	2 x 0.1ml	3 x 0.1ml	1 x 0.1ml
Murine IL-6 standard	10 vials	15 vials	5 vials
Streptavidin-HRP	2 vials	3 vials	1 vial

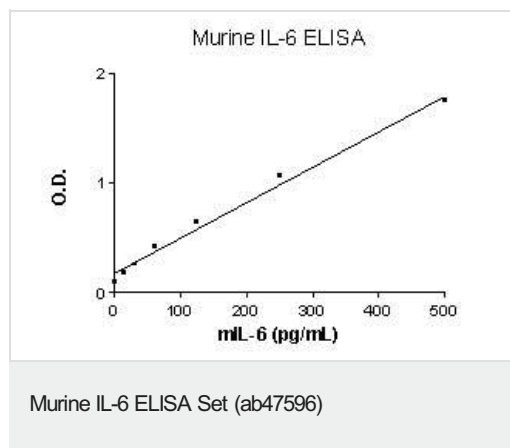
機能	Cytokine with a wide variety of biological functions. It is a potent inducer of the acute phase response. Plays an essential role in the final differentiation of B-cells into Ig-secreting cells. Involved in lymphocyte and monocyte differentiation. It induces myeloma and plasmacytoma growth and induces 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

**The Abpromise guarantee**      **Abpromise保証は、次のテスト済みアプリケーションにおけるab47596の使用に適用されます**  
 アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

アプリケーション	Abreviews	特記事項
Sandwich ELISA		Use at an assay dependent concentration.

## 画像



Mouse IL-6 Standard curve ranging from 15.6 to 500 pg/mL

**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.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