

### Human IL-1 beta ELISA Set (without plates) ab47351

#### 画像数 1

#### 製品の概要

**製品名** Human IL-1 beta 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** < 5 pg/ml

**検出範囲** 15.6 pg/ml - 500 pg/ml

**全工程の試験時間** 3h 45m

**ステップ** Multiple steps standard assay

**種交差性** 交差種: Human

**製品の概要** Abcam's IL-1 $\beta$  Human *in vitro* ELISA Set is designed for the quantitative measurement of IL-1 $\beta$  in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.

This ELISA Set is intended to measure human IL-1b in cell -culture supernatants, as well as in serum and plasma samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.

The IL-1b 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.

Recognizes both natural and recombinant human IL-1b.

**アプリケーション** 適用あり: Sandwich ELISA

**試験プラットフォーム** Reagents

#### 製品の特性

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

内容	ラベル	15 x 96 tests	10 x 96 tests	5 x 96 tests
Chromogen TMB Substrate Solution		6 vials	4 vials	2 vials

内容	ラベル	15 x 96 tests	10 x 96 tests	5 x 96 tests
Detection biotinylated anti-IL-1 beta antibody		3 x 0.55ml	2 x 0.55ml	1 x 0.55ml
IL-1 beta Capture Antibody		3 x 0.5ml	2 x 0.5ml	1 x 0.5ml
IL1 beta standard	Yellow	15 vials	10 vials	5 vials
Streptavidin-HRP		3 vials	2 vials	1 vial

<b>機能</b>	Potent proinflammatory cytokine. Initially discovered as the major endogenous pyrogen, induces prostaglandin synthesis, neutrophil influx and activation, T-cell activation and cytokine production, B-cell activation and antibody production, and fibroblast proliferation and collagen production. Promotes Th17 differentiation of T-cells.
<b>組織特異性</b>	Expressed in activated monocytes/macrophages (at protein level).
<b>配列類似性</b>	Belongs to the IL-1 family.
<b>翻訳後修飾</b>	Activation of the IL1B precursor involves a CASP1-catalyzed proteolytic cleavage. Processing and secretion are temporarily associated.
<b>細胞内局在</b>	Cytoplasm, cytosol. Lysosome. Secreted, exosome. Cytoplasmic vesicle, autophagosome. Secreted. The precursor is cytosolic. In response to inflammasome-activating signals, such as ATP for NLRP3 inflammasome or bacterial flagellin for NLRC4 inflammasome, cleaved and secreted. IL1B lacks any known signal sequence and the pathway(s) of its secretion is(are) not yet fully understood (PubMed:24201029). On the basis of experimental results, several unconventional secretion mechanisms have been proposed. 1. Secretion via secretory lysosomes: a fraction of CASP1 and IL1B precursor may be incorporated, by a yet undefined mechanism, into secretory lysosomes that undergo Ca(2+)-dependent exocytosis with release of mature IL1B (PubMed:15192144). 2. Secretory autophagy: IL1B-containing autophagosomes may fuse with endosomes or multivesicular bodies (MVBs) and then merge with the plasma membrane releasing soluble IL1B or IL1B-containing exosomes (PubMed:24201029). However, autophagy impacts IL1B production at several levels and its role in secretion is still controversial. 3. Secretion via exosomes: ATP-activation of P2RX7 leads to the formation of MVBs containing exosomes with entrapped IL1B, CASP1 and other inflammasome components. These MVBs undergo exocytosis with the release of exosomes. The release of soluble IL1B occurs after the lysis of exosome membranes (By similarity). 4. Secretion by microvesicle shedding: activation of the ATP receptor P2RX7 may induce an immediate shedding of membrane-derived microvesicles containing IL1B and possibly inflammasome components. The cytokine is then released in the extracellular compartment after microvesicle lysis (PubMed:11728343). 5. Release by translocation through permeabilized plasma membrane. This may occur in cells undergoing pyroptosis due to sustained activation of the inflammasome (By similarity). These mechanisms may not be mutually exclusive.

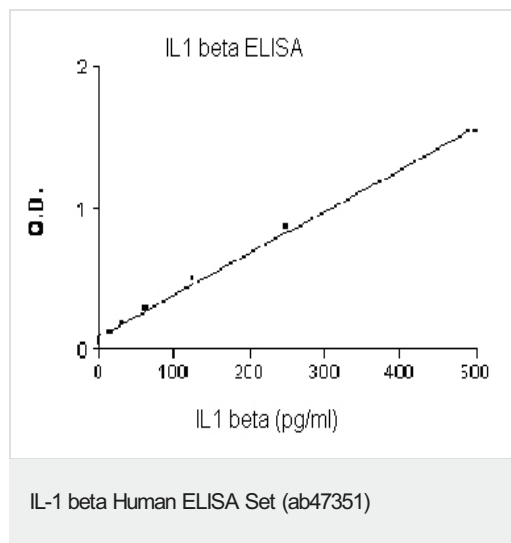
## アプリケーション

**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ab47351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
<b>Sandwich ELISA</b>		Use at an assay dependent concentration.

## 画像

---



IL-1 beta 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