

Human TNF beta ELISA Kit ab119576

1 References 画像数 1

製品の概要

製品名	Human TNF beta ELISA Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma, Other biological fluids, Tissue Extracts
アッセイタイプ	Sandwich (quantitative)
検出感度	< 5 pg/ml
検出範囲	62.5 pg/ml - 4000 pg/ml
ステップ	Multiple steps standard assay
種交差性	交差種: Human
製品の概要	<p>Abcam's Human TNF beta <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Human TNF beta in cell culture supernatants, serum and plasma (heparin, EDTA, citrate).</p> <p>A TNF beta specific mouse monoclonal antibody has been precoated onto 96-well plates. Standards and test samples are added to the wells and incubated. A biotinylated detection polyclonal antibody from goat, specific for TNF beta is then added followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. TMB is then used to visualize the HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the Human TNF Beta amount of sample captured in plate.</p>
試験プラットフォーム	Microplate

製品の特性

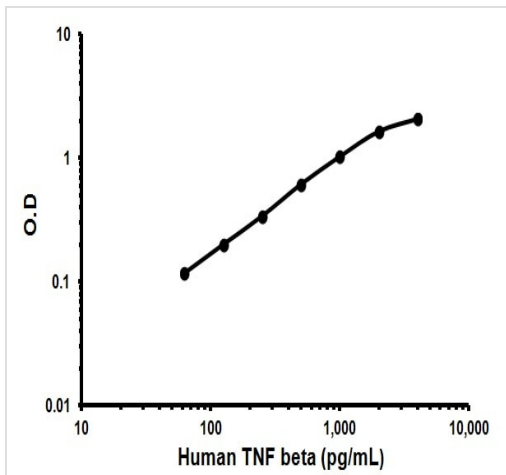
保存方法 Store at -20°C. Please refer to protocols.

内容	ラベル	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-Human TNF beta Antibody Microplate (12 x 8 wells)		1 unit

内容	ラベル	1 x 96 tests
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti-Human TNF beta antibody		1 x 100µl
Lyophilized recombinant Human TNF beta standard		2 x 10ng
Plate Seal		1 x 4 units
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Amber Bottle	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml
Wash Buffer (25X)		1 x 20ml

機能	Cytokine that in its homotrimeric form binds to TNFRSF1A/TNFR1, TNFRSF1B/TNFR2 and TNFRSF14/HVEM. In its heterotrimeric form with LTβ binds to TNFRSF3/LTβR. Lymphotoxin is produced by lymphocytes and cytotoxic for a wide range of tumor cells in vitro and in vivo.
関連疾患	Genetic variations in LTA are a cause of susceptibility psoriatic arthritis (PSORAS) [MIM:607507]. PSORAS is an inflammatory, seronegative arthritis associated with psoriasis. It is a heterogeneous disorder ranging from a mild, non-destructive disease to a severe, progressive, erosive arthropathy. Five types of psoriatic arthritis have been defined: asymmetrical oligoarthritis characterized by primary involvement of the small joints of the fingers or toes; asymmetrical arthritis which involves the joints of the extremities; symmetrical polyarthritis characterized by a rheumatoidlike pattern that can involve hands, wrists, ankles, and feet; arthritis mutilans, which is a rare but deforming and destructive condition; arthritis of the sacroiliac joints and spine (psoriatic spondylitis).
配列類似性	Belongs to the tumor necrosis factor family.
細胞内局在	Secreted. Membrane. The homotrimer is secreted. The heterotrimer is membrane-associated.

画像



Representative Standard Curve using ab119576.

Typical Standard Curve

**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