

Mouse TGF beta 1 ELISA Kit ab119557

★★★★★ [5 Abreviews](#)   [49 References](#)   [画像数 3](#)

製品の概要

製品名	Mouse TGF beta 1 ELISA Kit
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	4			7.9%

サンプル	N	平均値	SD	CV%
Overall	4			5.8%

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	7.8 pg/ml
検出範囲	31.3 pg/ml - 2000 pg/ml
添加回収試験	特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant		124% - 141%
Serum		92% - 103%

ステップ	Multiple steps standard assay
種交差性	交差種: Mouse
製品の概要	<p>Abcam's TGF beta 1 Mouse <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for accurate quantitative measurement of Mouse TGF beta 1 concentrations in Cell culture supernatant, serum and plasma (EDTA, citrate).</p> <p>Mouse TGF beta 1 specific antibodies have been precoated onto 96-well plates. Standards and test samples are added to the wells and then incubated at room temperature. After washing, a Biotin-conjugated anti-Mouse TGF beta 1 detection antibody is added then incubated at room temperature. Following washing Streptavidin-HRP conjugate is added to each well, incubated at</p>

room temperature then again washed. TMB is added and then catalyzed by HRP to produce a blue color product that changes into yellow after the addition of an acidic stop solution. The density of yellow coloration is directly proportional to the amount of Mouse TGF beta 1 captured on the plate

#### 試験プラットフォーム

Microplate

#### 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	2 x 5ml
20X Wash Buffer Concentrate	1 x 50ml
Adhesive Films	6 units
Biotinylated mouse anti-TGF beta 1 monoclonal antibody	1 x 120µl
Microplate coated with monoclonal antibody to mouse TGF beta 1 (12 x 8 wells)	1 unit
Mouse TGF beta 1 Standard lyophilized	2 vials
Stop Solution (1M Phosphoric acid)	1 x 15ml
Streptavidin-HRP	1 x 150µl
TMB Substrate Solution	1 x 15ml

#### 機能

Multifunctional protein that controls proliferation, differentiation and other functions in many cell types. Many cells synthesize TGFB1 and have specific receptors for it. It positively and negatively regulates many other growth factors. It plays an important role in bone remodeling as it is a potent stimulator of osteoblastic bone formation, causing chemotaxis, proliferation and differentiation in committed osteoblasts.

#### 組織特異性

Highly expressed in bone. Abundantly expressed in articular cartilage and chondrocytes and is increased in osteoarthritis (OA). Co-localizes with ASPN in chondrocytes within OA lesions of articular cartilage.

#### 関連疾患

Defects in TGFB1 are the cause of Camurati-Engelmann disease (CE) [MIM:131300]; also known as progressive diaphyseal dysplasia 1 (DPD1). CE is an autosomal dominant disorder characterized by hyperostosis and sclerosis of the diaphyses of long bones. The disease typically presents in early childhood with pain, muscular weakness and waddling gait, and in some cases other features such as exophthalmos, facial paralysis, hearing difficulties and loss of vision.

#### 配列類似性

Belongs to the TGF-beta family.

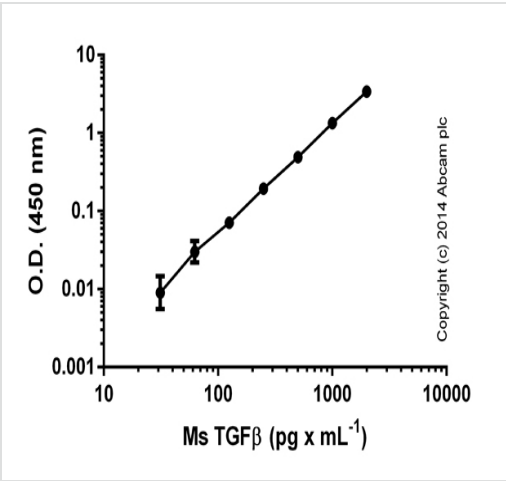
#### 翻訳後修飾

Glycosylated.

The precursor is cleaved into mature TGF-beta-1 and LAP, which remains non-covalently linked to mature TGF-beta-1 rendering it inactive.

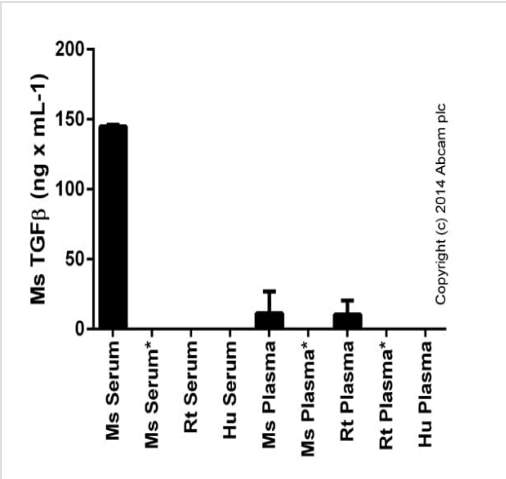
#### 細胞内局在

Secreted > extracellular space > extracellular matrix.



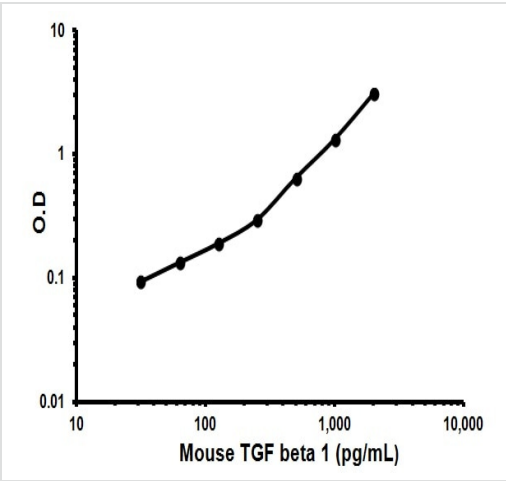
Standard curve of Ms TGFβ, with background signal subtracted (duplicates; +/- SD).

Sandwich ELISA - TGF beta 1 Mouse ELISA Kit (ab119557)



Mouse TGFβ measured in activated biological fluids, background signal subtracted (duplicates +/- SD). \* Samples without activation.

Sandwich ELISA - TGF beta 1 Mouse ELISA Kit (ab119557)



Representative Standard Curve using ab119557.

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