

★★★★★ **5 Abreviews** **49 References** 3 Images

Product name	Mouse TGF beta 1 ELISA Kit
Detection method	Colorimetric
Precision	Intra-assay

				Inter-assay
Sample	n	Mean	SD	CV%
Overall	4			5.8%

Range 31.3 pg/ml - 2000 pg/ml

Recovery	Sample specific recovery
----------	--------------------------

Sample type	Average %	Range
Cell culture supernatant		124% - 141%
Serum		92% - 103%

Product overview	<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

Platform Microplate

Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	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

Function 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.

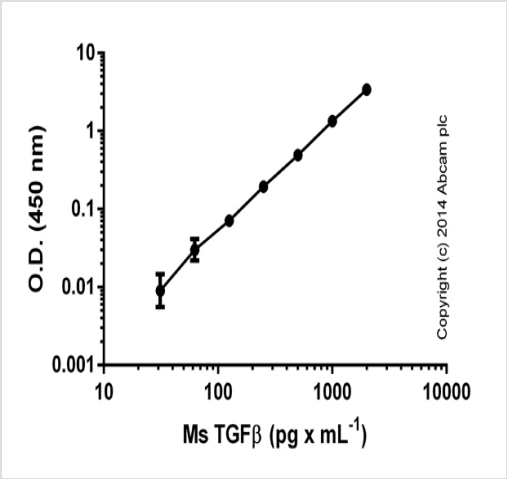
Tissue specificity Highly expressed in bone. Abundantly expressed in articular cartilage and chondrocytes and is increased in osteoarthritis (OA). Co-localizes with ASPN in chondrocytes within OA lesions of articular cartilage.

Involvement in disease 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.

Sequence similarities Belongs to the TGF-beta family.

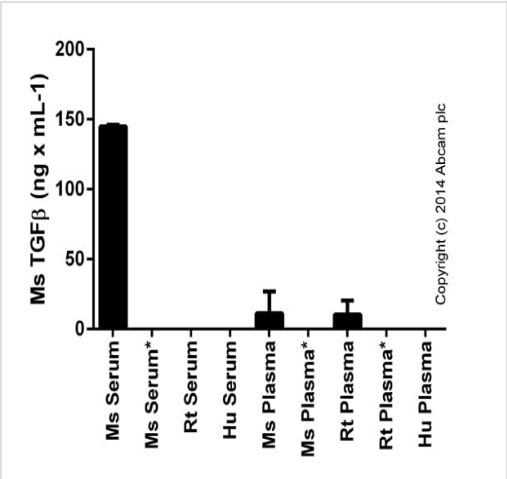
Post-translational modifications Glycosylated.
The precursor is cleaved into mature TGF-beta-1 and LAP, which remains non-covalently linked to mature TGF-beta-1 rendering it inactive.

Cellular localization Secreted > extracellular space > extracellular matrix.



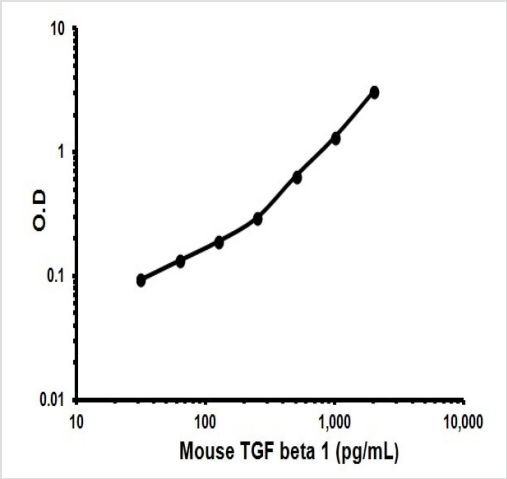
Standard curve of Ms TGFβ, with background signal subtracted (duplicates; +/- SD).

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.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors