

Human TGF beta 1 ELISA Kit ab108912

4 References 画像数 1

製品の概要

製品名	Human 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
	Overall	60		CV%
				6.5%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Overall	60		CV%
				10.9%
サンプルの種類	Cell culture supernatant			
アッセイタイプ	Sandwich (quantitative)			
検出感度	= 31 pg/ml			
検出範囲	0.063 ng/ml - 4 ng/ml			
添加回収試験	108 %			
全工程の試験時間	5h 00m			
ステップ	Multiple steps standard assay			
種交差性	交差種: Human			
製品の概要	Abcam's TGF beta 1 Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of TGF beta 1 concentrations in cell culture supernatants.			

A TGF beta 1 specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a TGF beta 1 specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase 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 amount of TGF beta 1 captured in plate. This assay recognizes both natural and recombinant Human TGF beta 1.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

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	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent M Concentrate	1 x 30ml
1X Standard Diluent	1 x 2ml
20X Wash Buffer Concentrate	2 x 30ml
30X Biotinylated Human TGF beta 1 Antibody	1 x 180µl
Chromogen Substrate	1 x 7ml
Sealing Tapes	3 units
Stop Solution	1 x 11ml
TGF beta 1 Microplate (12 x 8 well strips)	1 unit
TGF beta 1 Standard	1 vial

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.

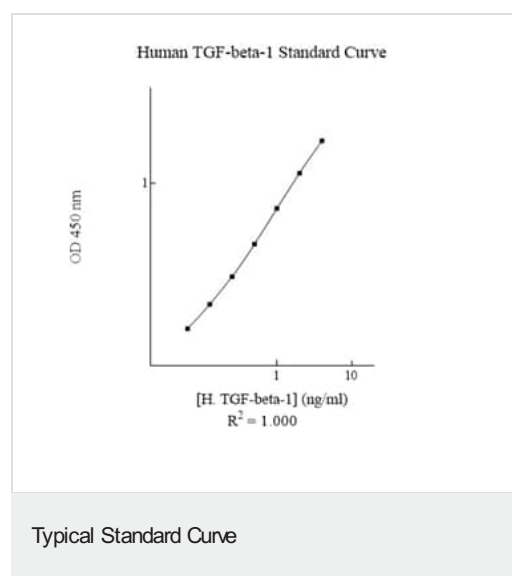
翻訳後修飾

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.



Representative Standard Curve using ab108912.

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