


Mouse Fibronectin ELISA Kit ab108849

6 References 画像数 2

製品の概要

製品名	Mouse Fibronectin 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			CV%
				4.5%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				9.6%
サンプルの種類	Cell culture supernatant, Urine, Serum, Plasma, Tissue, Cell Lysate			
アッセイタイプ	Sandwich (quantitative)			
検出感度	= 0.48 ng/ml			
検出範囲	3.125 ng/ml - 200 ng/ml			
添加回収試験	101 %			
全工程の試験時間	3h 30m			
ステップ	Multiple steps standard assay			
種交差性	交差種: Mouse			

交差が予測される動物種: Rat 

製品の概要

Mouse Fibronectin *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit (ab108849) is designed for the quantitative measurement of Fibronectin levels in plasma, serum, urine and cell culture supernatants.

A Fibronectin specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Fibronectin specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Complex 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 Fibronectin captured in plate.

**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 N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Mouse Fibronectin Antibody	1 x 120µl
Chromogen Substrate	1 x 7ml
Fibronectin Microplate (12 x 8 well strips)	1 unit
Fibronectin Standard (lyophilized)	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

#### 機能

Fibronectins bind cell surfaces and various compounds including collagen, fibrin, heparin, DNA, and actin. Fibronectins are involved in cell adhesion, cell motility, opsonization, wound healing, and maintenance of cell shape. Involved in osteoblast compaction through the fibronectin fibrillogenesis cell-mediated matrix assembly process, essential for osteoblast mineralization. Participates in the regulation of type I collagen deposition by osteoblasts.

Anastellin binds fibronectin and induces fibril formation. This fibronectin polymer, named superfibronectin, exhibits enhanced adhesive properties. Both anastellin and superfibronectin inhibit tumor growth, angiogenesis and metastasis. Anastellin activates p38 MAPK and inhibits lysophospholipid signaling.

#### 組織特異性

Plasma FN (soluble dimeric form) is secreted by hepatocytes. Cellular FN (dimeric or cross-linked multimeric forms), made by fibroblasts, epithelial and other cell types, is deposited as fibrils in the extracellular matrix. Ugl-Y1, Ugl-Y2 and Ugl-Y3 are found in urine.

#### 関連疾患

Glomerulopathy with fibronectin deposits 2

#### 配列類似性

Contains 12 fibronectin type-I domains.  
Contains 2 fibronectin type-II domains.  
Contains 16 fibronectin type-III domains.

#### 発生段階

Ugl-Y1, Ugl-Y2 and Ugl-Y3 are present in the urine from 0 to 17 years of age.

#### 翻訳後修飾

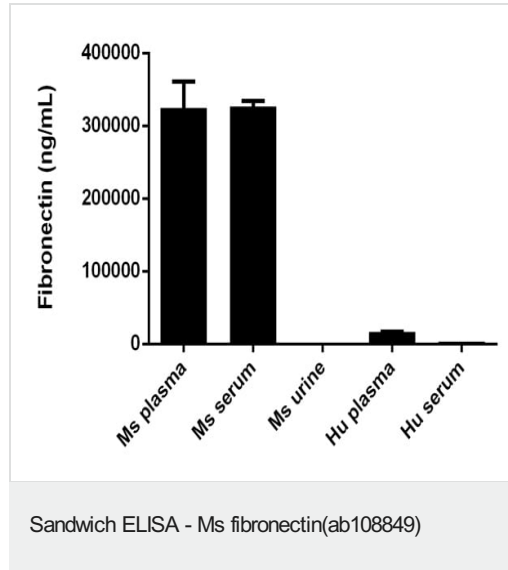
Sulfated.

It is not known whether both or only one of Thr-2064 and Thr-2065 are/is glycosylated.  
 Forms covalent cross-links mediated by a transglutaminase, such as F13A or TGM2, between a glutamine and the epsilon-amino group of a lysine residue, forming homopolymers and heteropolymers (e.g. fibrinogen-fibronectin, collagen-fibronectin heteropolymers).  
 Phosphorylated by FAM20C in the extracellular medium.  
 Proteolytic processing produces the C-terminal NC1 peptide, anastellin.

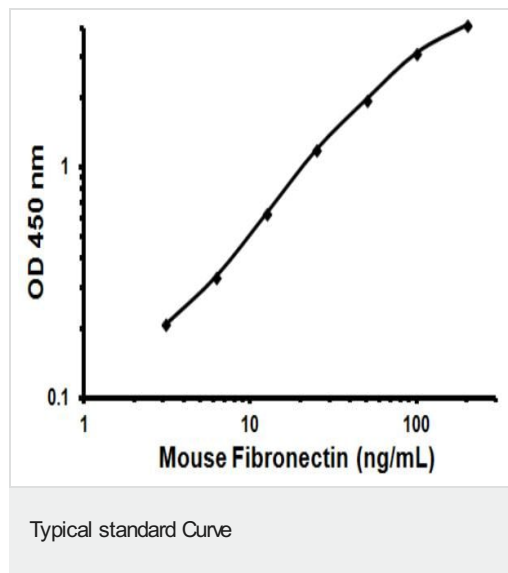
細胞内局在

Secreted, extracellular space, extracellular matrix.

## 画像



Fibronectin measured in biological fluids showing quantity (ng) per mL of sample tested



Representative Standard Curve using ab108849

**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