

Human Fibronectin Matched Antibody Pair Kit ab222262

リコンビナント

画像数 2

製品の概要

製品名	Human Fibronectin Matched Antibody Pair Kit
検出方法	Colorimetric
アッセイタイプ	ELISA set
検出感度	17.7 pg/ml
検出範囲	62.5 pg/ml - 4000 pg/ml
種交差性	交差種: Human
製品の概要	<p>Human Fibronectin Matched Antibody Pair Kits include a capture and a biotinylated detector antibody pair, along with a calibrated protein standard, suitable for sandwich ELISA. The Matched Antibody Pair Kit can be used to quantify native and recombinant human Fibronectin.</p> <p>Optimization of the kit reagents to sample type, immunoassay format or instrumentation may be required. Guidelines for use of this kit in a standard 96-well microplate sandwich ELISA using HRP/TMB system of colorimetric detection is described in this assay procedure for the purposes of quantification.</p> <p>Protocol information and tips on the use of the Matched Antibody Pair kits for sandwich ELISA can be found on our website. An accessory pack can be purchased which includes buffer reagents required to perform 10 x 96-well plate sandwich ELISAs (ab210905).</p> <p>For additional information on the performance of the antibody pair used in this kit, please see our equivalent SimpleStep ELISA kit ab219046. Please note that while the antibody pair is the same provided in the corresponding SimpleStep ELISA Kit, due to differences in their formulation, this antibody pair cannot be used with the consumables provided with our SimpleStep ELISA Kits.</p>
アプリケーション	適用あり: ELISA
試験プラットフォーム	Reagents

製品の特性

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

内容	5 x 96 tests
Human Fibronectin Capture Antibody	1 x 50µg
Human Fibronectin Detector Antibody	1 x 12.5µg
Human Fibronectin Lyophilized Protein	1 vial

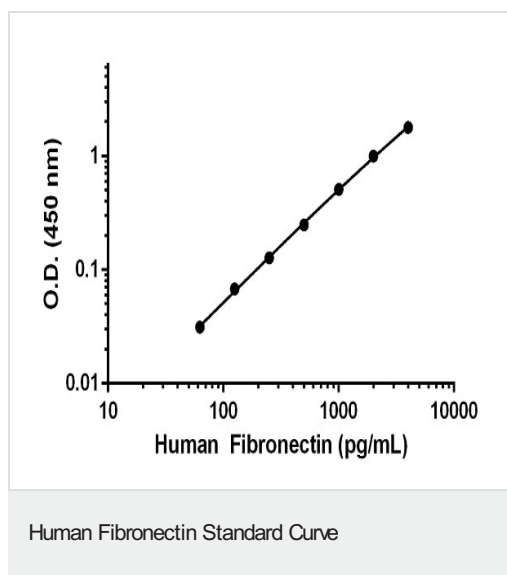
機能	<p>Fibronectins bind cell surfaces and various compounds including collagen, fibrin, heparin, DNA, and actin. Fibronectins are involved in cell adhesion, cell motility, opsonization, wound healing, and maintenance of cell shape. Involved in osteoblast compaction through the fibronectin fibrillogenesis cell-mediated matrix assembly process, essential for osteoblast mineralization. Participates in the regulation of type I collagen deposition by osteoblasts.</p> <p>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.</p>
組織特異性	<p>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.</p>
関連疾患	<p>Glomerulopathy with fibronectin deposits 2</p>
配列類似性	<p>Contains 12 fibronectin type-I domains.</p> <p>Contains 2 fibronectin type-II domains.</p> <p>Contains 16 fibronectin type-III domains.</p>
発生段階	<p>Ugl-Y1, Ugl-Y2 and Ugl-Y3 are present in the urine from 0 to 17 years of age.</p>
翻訳後修飾	<p>Sulfated.</p> <p>It is not known whether both or only one of Thr-2064 and Thr-2065 are/is glycosylated.</p> <p>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).</p> <p>Phosphorylated by FAM20C in the extracellular medium.</p> <p>Proteolytic processing produces the C-terminal NC1 peptide, anastellin.</p>
細胞内局在	<p>Secreted, extracellular space, extracellular matrix.</p>

アプリケーション

The Abpromise guarantee **Abpromise保証は、次のテスト済みアプリケーションにおけるab222262の使用に適用されます**
 アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

アプリケーション	Abreviews	特記事項
ELISA		Use at an assay dependent concentration.

画像



Standard calibration curve. Background subtracted values are graphed.

Powered by
recombinant antibodies

Research with confidence
Consistent and reproducible results

Long-term and scalable supply
Recombinant technology

Success from the first experiment
Confirmed specificity

Ethical standards compliant
Animal-free production

Sandwich ELISA - Human Fibronectin Matched
Antibody Pair Kit (ab222262)

To learn more about the advantages of recombinant antibodies see [here](#).

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