

Human Syndecan-1 ELISA Kit (CD138) ab46506

[26 References](#) [画像数 4](#)

製品の概要

製品名 Human Syndecan-1 ELISA Kit (CD138)

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	6			6.2%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Overall	6			10.2%

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 = 4.94 ng/ml

検出範囲 8 ng/ml - 256 ng/ml

添加回収試験 101 %

全工程の試験時間 1h 45m

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 Human Syndecan-1 ELISA (CD138) kit is designed for the quantitative measurement of Syndecan-1 in Human serum, plasma, buffered solutions and supernatants.

A monoclonal antibody specific for Syndecan-1 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known Syndecan-1 concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for Syndecan-1 are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of Syndecan-1 present in the samples.

This kit will recognize both endogenous and recombinant Human Syndecan-1.

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	ラベル	1 x 96 tests	2 x 96 tests	1 x 96 tests	2 x 96 tests
10X Standard Diluent Buffer	Black	1 x 15ml	1 x 25ml	1 x 15ml	1 x 25ml
200X Wash Buffer	White	1 x 10ml	2 x 10ml	1 x 10ml	2 x 10ml
Biotinylated Antibody Diluent	Red	1 x 7ml	1 x 13ml	1 x 7ml	1 x 13ml
Biotinylated anti-Syndecan-1	Red	1 x 400µl	2 x 400µl	1 x 400µl	2 x 400µl
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml	1 x 11ml	1 x 24ml
Control	Yellow	2 vials	4 vials	2 vials	4 vials
HRP Diluent	Red	1 x 12ml	1 x 23ml	1 x 12ml	1 x 23ml
Plastic Plate Covers		2 units	4 units	2 units	4 units
Stop Reagent	Black	1 x 11ml	2 x 11ml	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl	2 x 5µl	4 x 5µl
Syndecan-1 Microplate (12 x 8 well strips)		1 unit	2 units	1 unit	2 units
Syndecan-1 Standard (lyophilized)	Orange	2 vials	4 vials	2 vials	4 vials

機能

Cell surface proteoglycan that bears both heparan sulfate and chondroitin sulfate and that links the cytoskeleton to the interstitial matrix.

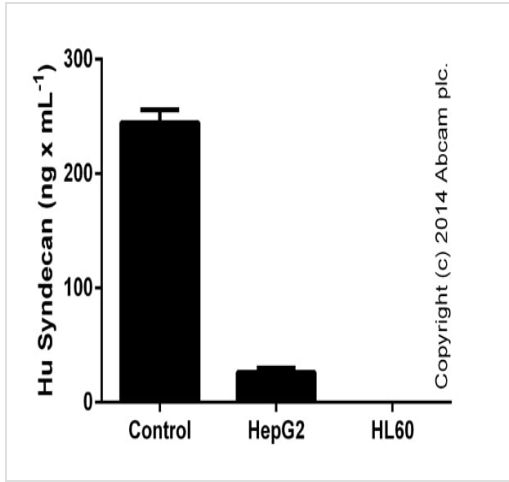
配列類似性

Belongs to the syndecan proteoglycan family.

細胞内局在

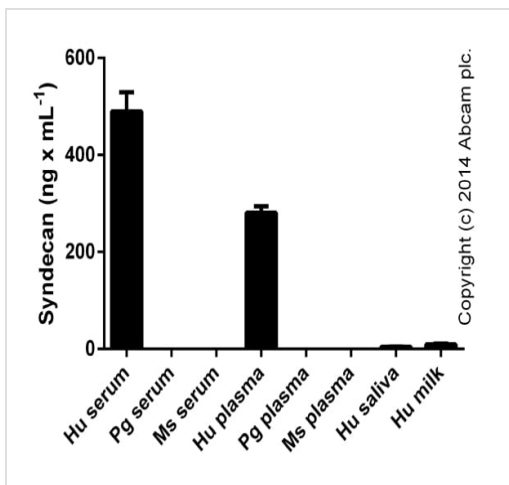
Membrane.

画像



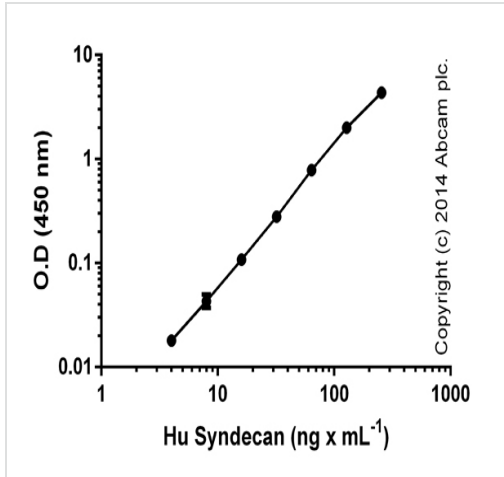
Syndecan measured in undiluted cell culture supernatants, including the positive control from kit (duplicates +/- SD).

ELISA: Syndecan (CD138) Human ELISA Kit (ab46506)



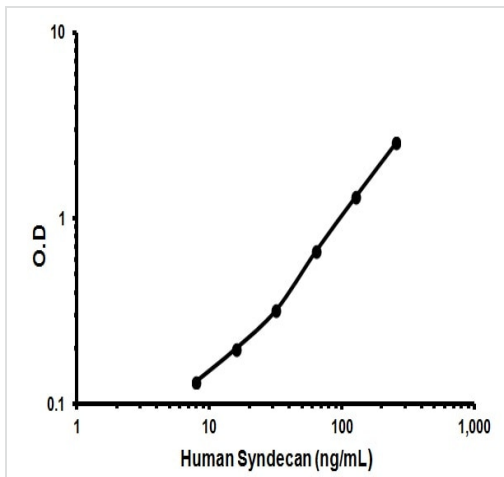
Syndecan measured in human biological fluids (dilution range 1/1-1/10); mouse and pig samples (plasma and serum) were below limit of detection (duplicates +/- SD).

ELISA: Syndecan (CD138) Human ELISA Kit (ab46506)



Standard curve of human syndecan with background signal subtracted (duplicates; +/- SD).

ELISA: Syndecan (CD138) Human ELISA Kit
(ab46506)



Representative Standard Curve using ab46506

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