

Human Syndecan-1 ELISA Set (without plates) ab47352

3 References [画像数 1](#)

製品の概要

製品名 Human Syndecan-1 ELISA Set (without plates)

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

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 Abcam's Syndecan-1 Human *in vitro* ELISA Set is designed for the quantitative measurement of Syndecan-1 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.

This ELISA Set is intended to measure human Syndecan-1 in cell -culture supernatants, as well as in serum and plasma samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.

The Syndecan-1 ELISA Set includes both the capture and the detection antibodies as well as Horse-Radish peroxidase conjugated Streptavidin (HRP-Strep) and TMB which are the reagents used in the colour development step.

Recognizes both natural and recombinant human Syndecan-1.

アプリケーション **適用あり:** Sandwich ELISA

試験プラットフォーム Reagents

製品の特性

保存方法 Store at +4°C. Please refer to protocols.

内容	ラベル	15 x 96 tests	10 x 96 tests	5 x 96 tests
Chromogen TMB Substrate Solution		6 vials	4 vials	2 vials

内容	ラベル	15 x 96 tests	10 x 96 tests	5 x 96 tests
Detection biotinylated anti-Syndecan-1 antibody		3 x 0.55ml	2 x 0.55ml	1 x 0.55ml
Streptavidin-HRP		3 vials	2 vials	1 vial
Syndecan-1 Capture Antibody		3 x 0.5ml	2 x 0.5ml	1 x 0.5ml
Syndecan-1 Standard (lyophilized)	Yellow	15 vials	10 vials	5 vials

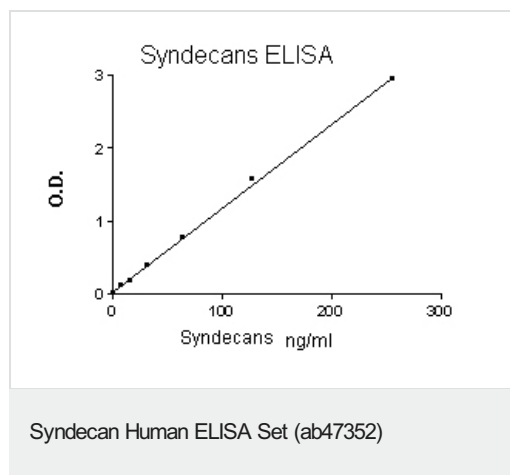
機能	Cell surface proteoglycan that bears both heparan sulfate and chondroitin sulfate and that links the cytoskeleton to the interstitial matrix.
配列類似性	Belongs to the syndecan proteoglycan family.
細胞内局在	Membrane.

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ab47352の使用に適用されず**
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
Sandwich ELISA		Use at an assay dependent concentration.

画像



Syndecans standard curve ranging from 8 to 256 ng/ml

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