

Product datasheet

Human Syndecan-1 ELISA Set (without plates) ab47352

[3 References](#) [1 Image](#)

Overview

Product name	Human Syndecan-1 ELISA Set (without plates)
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 4.94 ng/ml
Range	8 ng/ml - 256 ng/ml
Assay time	1h 45m
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human
Product overview	<p>Abcam's Syndecan-1 Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of Syndecan-1 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>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.</p> <p>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.</p> <p>Recognizes both natural and recombinant human Syndecan-1.</p>
Tested applications	Suitable for: Sandwich ELISA
Platform	Reagents

Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	Identifier	15 x 96 tests	10 x 96 tests	5 x 96 tests
Chromogen TMB Substrate Solution		6 vials	4 vials	2 vials

Components	Identifier	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

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

Sequence similarities Belongs to the syndecan proteoglycan family.

Cellular localization Membrane.

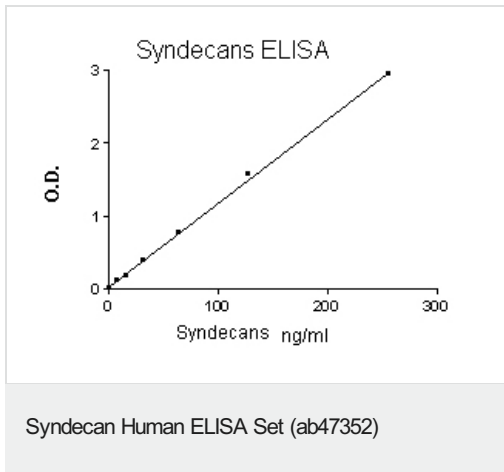
Applications

The Abpromise guarantee Our **Abpromise guarantee** covers the use of ab47352 in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

Images



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.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors