

Product datasheet

Human Decorin ELISA Kit (DCN) α b99998

★★★★★ [1 Abreviews](#) [3 References](#) [2 Images](#)

Overview

Product name	Human Decorin ELISA Kit (DCN)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 1.5 pg/ml
Range	0.96 pg/ml - 700 pg/ml
Recovery	> 93 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	93.82	84% - 104%
Serum	122.5	111% - 130%
Plasma	111.8	103% - 119%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's Decorin (DCN) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human Decorin in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human Decorin coated on a 96-well plate. Standards and samples are pipetted into the wells and Decorin present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti Human Decorin antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of Decorin bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Notes Optimisation may be required with urine samples.

The standard is affinity purified.

Platform

Microplate

Properties**Storage instructions**

Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
Biotinylated anti-Human Decorin	2 vials
Decorin Microplate (12 x 8 well strips)	1 unit
Recombinant Human Decorin Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Function

May affect the rate of fibrils formation.

Involvement in disease

Defects in DCN are the cause of congenital stromal corneal dystrophy (CSCD) [MIM:610048]. Corneal dystrophies are inherited, bilateral, primary alterations of the cornea that are not associated with prior inflammation or secondary to systemic disease. Most show autosomal dominant inheritance.

Sequence similarities

Belongs to the small leucine-rich proteoglycan (SLRP) family. SLRP class I subfamily. Contains 12 LRR (leucine-rich) repeats.

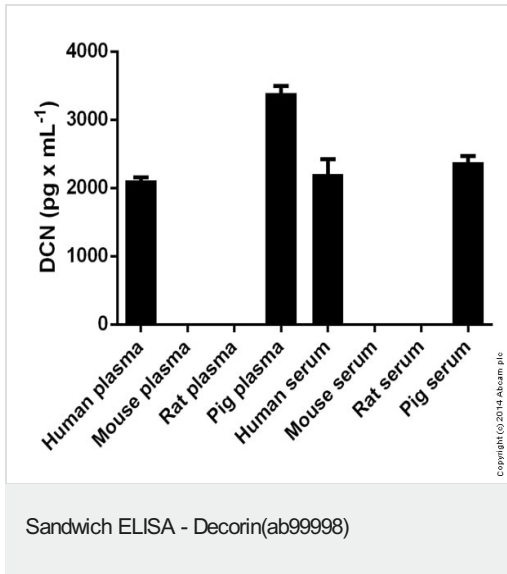
Post-translational modifications

The attached glycosaminoglycan chain can be either chondroitin sulfate or dermatan sulfate depending upon the tissue of origin.

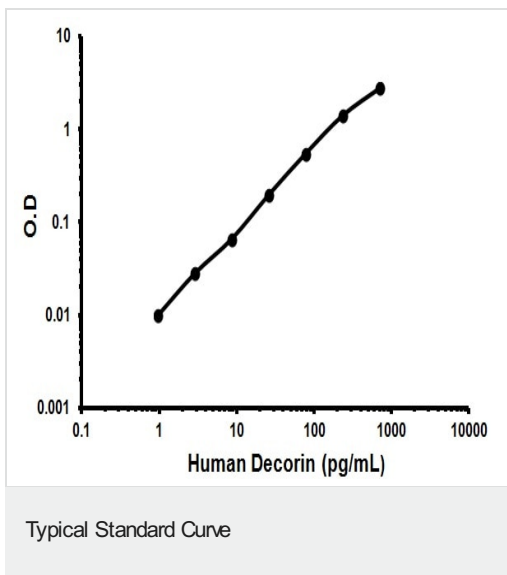
Cellular localization

Secreted > extracellular space > extracellular matrix.

Images



DCN measured in biological fluids showing quantity (pg) per mL of tested sample



Representative Standard Curve using ab99998.

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