

Mouse Decorin ELISA Kit (DCN) ab155454

3 References 1 Image

Overview

Product name Mouse Decorin ELISA Kit (DCN)

Detection method Colorimetric

Precision Intra-assay

Sample	n	Mean	SD	CV%
All				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
All				< 12%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity < 5 pg/ml

Range 4.1 pg/ml - 1000 pg/ml

Recovery > 68 %

Sample specific recovery

Sample type	Average %	Range
Serum	104.7	83% - 126%
Plasma	93.91	74% - 119%
Cell culture media	111.1	68% - 130%

Assay time 5h 00m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview Abcam's Decorin (DCN) Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay designed for the quantitative measurement of mouse Decorin in serum, plasma (collect plasma using EDTA and heparin as an anticoagulant. Citrate is

not recommended as anticoagulants for use in this assay).

This assay employs an antibody specific for mouse 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-mouse 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.

Get better reproducibility in only 90 minutes with Mouse Decorin ELISA Kit ([ab207618](#)) from our SimpleStep ELISA® range.

Notes Optimization may be required with urine samples.

Platform Microplate

Properties

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

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent	1 x 15ml
Biotinylated anti-mouse Decorin	2 vials
Decorin Microplate (12 x 8 wells)	1 unit
Recombinant Mouse Decorin Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

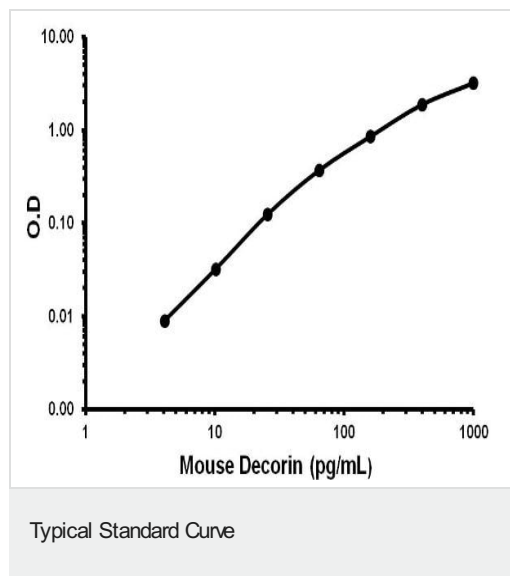
Function May affect the rate of fibrils formation.

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.



Representative standard curve using ab155454.

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