

Mouse Decorin ELISA Kit ab207618

Recombinant SimpleStep ELISA[®]

[6 Images](#)

Overview

Product name Mouse Decorin ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Serum	5			2.52%

Inter-assay

Sample	n	Mean	SD	CV%
Serum	3			6%

Sample type

Cell culture supernatant, Urine, Serum, Hep Plasma, EDTA Plasma, Cit plasma

Assay type

Sandwich (quantitative)

Sensitivity

0.13 ng/ml

Range

0.63 ng/ml - 40 ng/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Urine	79	71% - 86%
Serum	102	97% - 106%
Cell culture media	98	97% - 98%
Hep Plasma	106	97% - 110%
EDTA Plasma	90	89% - 91%
Cit plasma	86	84% - 90%

Assay time	1h 30m
Assay duration	One step assay
Species reactivity	Reacts with: Mouse Does not react with: Goat, Cow, Pig

Product overview Mouse Decorin ELISA Kit (ab207618) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Decorin protein in cell culture supernatant, cit plasma, edta plasma, hep plasma, serum, and urine. It uses our proprietary SimpleStep ELISA® technology. Quantitate Mouse Decorin with NULL NULL sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate (**ab203359**) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

Notes Decorin is a secreted proteoglycan of the small leucine-rich proteoglycan (SLRP) family. Decorin binds to type I and type II collagen fibrils as well as to fibronectin, TGF-beta and epidermal growth factor receptor.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

Platform Microplate (12 x 8 well strips)

Properties

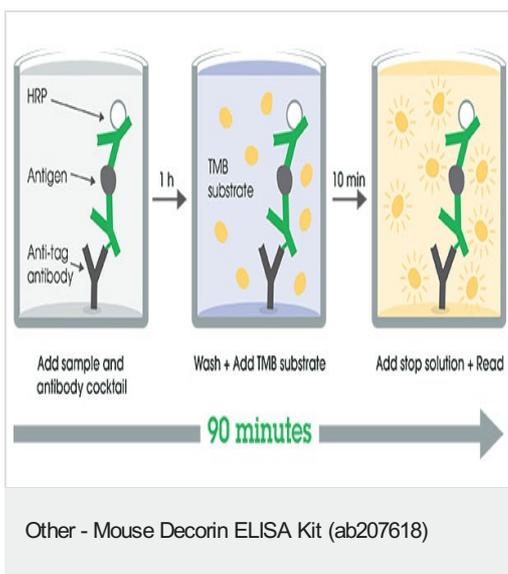
Storage instructions Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X Mouse Decorin Capture Antibody	1 x 600µl
10X Mouse Decorin Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
50X Cell Extraction Enhancer Solution (ab193971)	1 x 1ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml

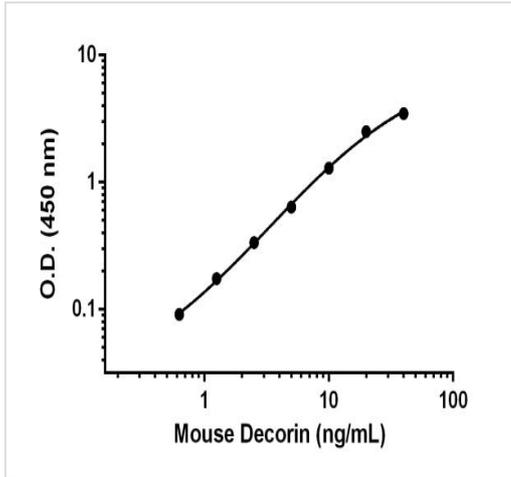
Components	1 x 96 tests
Antibody Diluent 5BR	1 x 6ml
Mouse Decorin Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NBS	1 x 20ml
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 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

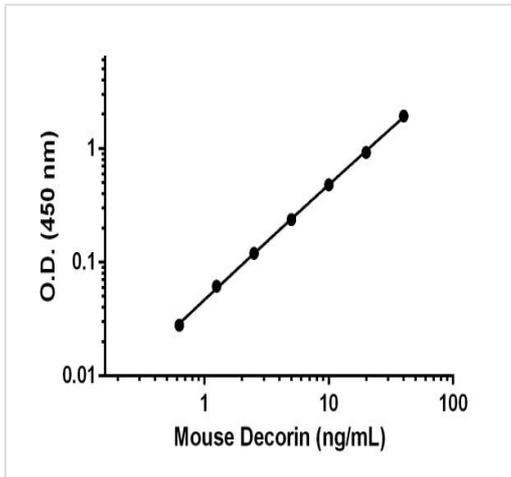


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



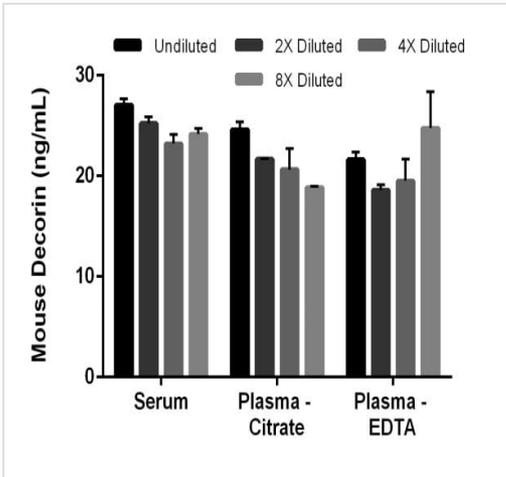
Background-subtracted data values (mean +/- SD) are graphed.

Example of mouse Decorin standard curve in Sample Diluent NS with enhancer.



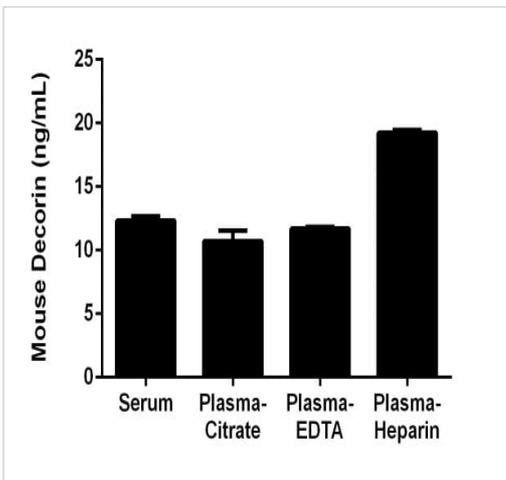
Background-subtracted data values (mean +/- SD) are graphed.

Example of mouse Decorin standard curve in Sample Diluent 15BS with enhancer.



Interpolated concentrations of spiked Decorin in mouse serum, and plasma samples.

The concentrations of Decorin were measured in duplicates, interpolated from the Decorin standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 25%, plasma (citrate) 50%, and plasma (EDTA) 25%.



Interpolated concentrations of Decorin in mouse serum, and plasma samples.

The concentrations of Decorin were measured in duplicates, interpolated from the Decorin standard curves and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean Decorin concentration was determined to be 12.3 ng/mL in serum, 10.7 ng/mL in plasma (citrate), 11.7 ng/mL in plasma (EDTA), and 19.2 ng/mL in plasma (heparin).

Powered by recombinant antibodies

- Research with confidence**
Consistent and reproducible results
- Long-term and scalable supply**
Recombinant technology
- Success from the first experiment**
Confirmed specificity
- Ethical standards compliant**
Animal-free production

To learn more about the advantages of recombinant antibodies see [here](#).

Sandwich ELISA - Mouse Decorin ELISA Kit
(ab207618)

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