

# Syndecan-2 (human), (recombinant) (Fc-tag)

The key players in the miR-20a-5p/SDC2 axis may be a potential diagnostic biomarker and therapeutic target for osteosarcoma patients. Also, SDC2 methylation as a blood-based DNA test for early detection of colorectal cancer.

## Ordering Information

[Order Online »](#)

ENZ-PRT321-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

kDa

M

116

66.2

45.0

35.0

25.0

18.4

14.4



As a result of glycosylation, the recombinant protein migrates as an approximately 109.5 kDa protein in SDS-PAGE under reducing conditions.

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** CD362, Fibroglycan, Heparan sulfate proteoglycan core protein, HSPG

**Endotoxin Content** <1 EU/μg (LAL method)

**Formulation** Lyophilized from sterile PBS, pH 7.4.

**MW** ~59.5 kDa (SDS-PAGE)

**Purity** ≥95% (SDS-PAGE)

**Reconstitution** Reconstitute with sterile deionized water. Reconstitution instructions are lot specific.

**Source** Produced in HEK293 cells. A DNA sequence encoding the human SDC2 (Glu19-Glu144, H229Y) was expressed with the Fc region of human IgG1 at the C-terminus.

**UniProt ID** P34741



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)