Perlecan monoclonal antibody (A76)

Perlecan is an extracellular matrix proteoglycan. It has a large core protein of 400-450kDa and is often produced with heparan sulfate side chains. Perlecan is found in basement membranes where it contributes to permeability charateristics, serves as a substrate for vascular cells, and binds growth factors involved in vascular remodelling.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BPD-CSI-001-76-1	1mg
BPD-CSI-001-76-04	400µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Basement membrane-specific heparan sulfate

proteoglycan core protein, HSPG, PLC

Application Affinity chromatography, ELISA, IHC (FS), IHC (PS), IP,

WB

Clone A76

CrossreactivityThere is no evidence for cross-reactivity with other

connective tissue proteins (vitronectin, fibronectin, elastin,

collagen, laminin).

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M

sodium chloride and 15mM sodium azide.

GenBank ID 444872

Host Mouse

Immunogen Lysed bovine corneal endothelial cells and extracellular

matrix.

lgG1κ

Purity Detail Protein A/G-purified.

Source From culture supernatant.

Species Reactivity Bovine, Human

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

-

Last modified: May 29, 2024



info-

eu@enzolifesciences.com