

Instruction Manual

Catalog Number	PK-AB577-3644
Description	Syndecans are type I integral membrane proteoglycans that are involved in cell-extracellular matrix adhesion and growth factor binding. Syndecan-1 is an matrix receptor which binds to Collagens, Fibronectin and Thrombospondin. Syndecan-1 and Syndecan-3 interact with monkey (midkine), a growth/differentiation factor involved in embryogenesis. Syndecan-2 is highly expressed at areas of high morphogenetic activity, such as epithelial-mesenchymal interfaces and the prechondrogenic and preosteogenic mesenchymal condensations. Syndecan-4 functions cooperatively with integrins in the processes of cell spreading and focal adhesion assembly.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Synthetic peptide surrounding amino acid 184 of human Syndecan-4.
Clone / IgG Subtype	Rabbit IgG
Species Reactivity	human, mouse, rat
Specificity	See Applications or Species Reactivity.
Formulation	100 µg (0.5 mg/ml) Protein A purified rabbit anti-Syndecan-4 polyclonal antibody in phosphate buffered saline (PBS), pH 7.2, containing 30% glycerol, 0.5% BSA, 0.01% thimerosal.
Reconstitution	Can be diluted in other aqueous buffers at the concentrations determined for the respective application.
Storage & Stability	Store at -20°C. For long-term storage, aliquot and refreeze at -70°C. Avoid repeated freeze/thaw cycles.
Applications	Western blotting (0.5-4 µg/ml), immunoprecipitation (5-10 µg/ml) and immunohistochemistry (frozen and paraffin sections; 15-25 µg/ml). The optimal conditions should be determined individually. Antibody might be suitable for other applications not tested so far.

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.