

Purified Mouse Anti-species independent N-sulfated heparan sulfate Antibody
HepSS-1, monoclonal

Catalog number: V1031880

Unit size: 0.1 mg

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Tris buffered saline (TBS) solution, 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Species independent
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	N-sulfated heparan sulfate
Clone	HepSS-1

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	WB, ICC, FC (QC TESTED), IHC, IP

Applications

Heparan sulfate has demonstrated use in a wide range of applications, especially in the field of biotechnology. It has been widely used in the application of key cellular activities such as anti-inflammatory activities and antiresorption activities, specifically in cell-lines like escort cells, Purkinje cells and 293-EBNA cells. It, moreover, has a variety of applications in the examination of key biological diseases like neurodegenerative disease, respiratory disease and aging-associated diseases, and it has been widely used in the study of important cell-based receptors such as phagocytic receptors and Dpp receptors.