

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

Catalog number: V1031890
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	Phosphate-buffered saline (PBS, pH 7.2), 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
Conjugate	PE

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE under optimal conditions
Application	FC (QC TESTED)

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.