

**PE Mouse Anti-human NG2 Antibody *7.1,
monoclonal***Catalog number: V1031845
Unit size: 100 tests**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	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	NG2
Clone	7.1
Conjugate	PE

Biological Properties

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

Applications

Chondroitin sulfate proteoglycan 4 (sometimes called Melanoma-associated chondroitin sulfate proteoglycan, Melanoma chondroitin sulfate proteoglycan or Chondroitin sulfate proteoglycan NG2) is a transmembrane protein with a molecular weight of 251 kDa, located in the collagen-containing extracellular matrix, cell surface and apical plasma membrane of cells. In Homo sapiens, chondroitin sulfate proteoglycan 4 reacts with protein kinase. It plays a role in the upregulation of peptidyl-tyrosine phosphorylation. Chondroitin sulfate proteoglycan 4 is the subject of intensive study due to the fact that it acts in the platelet-derived growth factor receptor signaling pathway. It aids in organismal processes, namely, angiogenesis, glial cell migration and chondroitin sulfate biosynthetic process. Sequencing of chondroitin sulfate proteoglycan 4 has supported it contains 4 types of conserved structural units: Laminin G-like 1, Laminin G-like 2, cytoplasmic, and extracellular domain.