

**Purified Mouse Anti-human NG2 Antibody**  
**\*7.1, monoclonal\***

Catalog number: V1031840

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	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	NG2
Clone	7.1

**Biological Properties**

---

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

**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.