

**mFluor™ Green 620 Anti-human CD111
Antibody *R1.302***Catalog number: 111100U0, 111100U1
Unit size: 100 tests, 500 tests**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	0.1 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1 kappa
Immunogen	CD111 (Nectin-1, HVEC, PVRL1, PRR1)
Clone	R1.302
Conjugate	mFluor™ Green 620

Biological Properties

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Green 620 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Green 620
Excitation Wavelength	525 nm
Emission Wavelength	623 nm

Applications

R1.302 is an anti-human monoclonal antibody that forms an immune complex with the CD111 antigen. CD111 (also known as PVRL1 or Nectin-1) is a 75 kD single-pass type I membrane protein that is found on the surface of cells like erythrocytes, endothelial cells, epithelial cells, stem cells and macrophages. CD111 has been thought to be involved with critical biological processes like cell-cell adhesion, particularly

homophilic cell adhesion via plasma membrane adhesion molecules, and is associated with a variety of biologically interesting macromolecules/ligands, for instance, nectin3 and afadin gd. CD111 is a relatively rare antibody target, with fewer than 90 publications in the last decade. Even still, CD111 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of neuroscience. This antibody was purified through affinity chromatography and conjugated to mFluor™ Green 620 (ex/em = 525/623 nm). It is compatible with the 532 nm laser and 610/20 nm bandpass filter (for example, as in the BD FACSymphony™ A5).