

iFluor™ 405 Anti-human CD111 Antibody
R1.302Catalog number: 11110020, 11110021
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	iFluor™ 405

Biological Properties

Appearance	Light yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 405 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 405
Excitation Wavelength	403 nm
Emission Wavelength	427 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 iFluor™ 405 (ex/em = 403/427 nm). It is compatible with the 405 nm laser and 445/45 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte Quanteon).