

**mFluor™ Green 620 Anti-human CD2
Antibody *RPA-2.10***Catalog number: 100210U0, 100210U1
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
Immunogen	CD2 (LFA-2, Erythrocyte R, T11)
Clone	RPA-2.10
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

RPA-2.10 is an anti-human monoclonal antibody that is specific for the CD2 antigen. CD2 (sometimes referred to as T11, Rosette receptor, Erythrocyte receptor or LFA-2) is a 45 kD single-pass type I membrane protein that is located on the surface of cells such as T cells, B cells and NK cells. In certain organisms, CD2 positively regulates interleukin-8 secretion, enhances myeloid dendritic cell activation and is an enhancer of interferon-gamma secretion. Also, it has been thought to be involved with important biological processes like cell-cell adhesion, specifically heterotypic cell-cell adhesion. CD2 is involved with key cellular pathways, in particular, the cell surface receptor signaling pathway. From a

research standpoint, it is of biological interest due to its association with key macromolecules/ligands like CD48, LFA-3 and CD58. CD2 is a moderately popular antibody target, with over 16000 publications in the last decade. CD2 is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of costimulatory molecules and immunology. 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 620/15 nm bandpass filter (for example, as in the Thermo Fisher Attune NxT).