

**TRITC Anti-human CD2 Antibody**  
**\*RPA-2.10\***Catalog number: 10021110, 10021111  
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	TRITC

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with TRITC under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

---

Conjugate	TRITC
Excitation Wavelength	544 nm
Emission Wavelength	570 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 TRITC (ex/em = 544/570 nm). It is compatible with the 561 nm laser and 586/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte).