

APC Anti-human CD2 Antibody *AHC0031*Catalog number: 100221C0, 100221C1, 100221C2
Unit size: 25 tests, 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	IgG1
Immunogen	CD2 (LFA-2, Erythrocyte R, T11)
Clone	AHC0031
Conjugate	APC

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	APC
Excitation Wavelength	651 nm
Emission Wavelength	660 nm

Applications

AHC0031 is an anti-human monoclonal antibody that recognizes the CD2 antigen. CD2 (also known as T11, SRBC or LFA-3 receptor) is a 45 kD single-pass type I membrane protein that is found on the surface of cells like NK cells and T cells. CD2 has been closely linked to vital biological processes like cell-cell adhesion, especially heterotypic cell-cell adhesion. Also, in certain organisms, it promotes interleukin-8 secretion, plays a role in the upregulation of myeloid dendritic cell activation and is an enhancer of tumor necrosis factor production. CD2 is a member of critical cellular pathways, namely, the cell surface receptor signaling pathway. From a research standpoint, it is of biological interest due to its association with important macromolecules/ligands like CD58. CD2 is a moderately popular antibody target, with over 16000 publications in the last decade. CD2 is essential for costimulatory molecules and immunology research, frequently serving as a phenotypic marker for

differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to APC (ex/em = 651/660 nm). It is compatible with the 642 nm laser and 702/85 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).