

APC/Cy7 Anti-human CD101 Antibody
BB27Catalog number: 110101C0, 110101C1, 110101C2
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	Mouse IgG1
Immunogen	CD101 (IGSF2)
Clone	BB27
Conjugate	APC/Cy7

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/Cy7 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	APC/Cy7
Excitation Wavelength	754 nm
Emission Wavelength	779 nm

Applications

The BB27 monoclonal antibody binds to human CD101, a 120 kD member of the Ig superfamily commonly found on the surface of T cells, granulocytes and dendritic cells. CD101 acts in important cellular pathways, for instance, the cell surface receptor signaling pathway. Additionally, in some organisms, it is an enhancer of myeloid leukocyte differentiation. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands. CD101 is a relatively rare antibody target, with fewer than 200 publications in the last decade. Even still, CD101 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of immunology. This antibody was purified through affinity chromatography and conjugated to APC/Cy7 (ex/em = 754/779 nm).