

**APC/iFluor™ 700 Anti-human CD27
Antibody *LT27***Catalog number: 102701F0, 102701F1, 102701F2
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 IgG2a
Immunogen	CD27 (T14, S152, TNFRSF7)
Clone	LT27
Conjugate	APC/iFluor™ 700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/iFluor™ 700 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	APC/iFluor™ 700
Excitation Wavelength	685 nm
Emission Wavelength	710 nm

Applications

LT27 is an anti-human monoclonal antibody that recognizes the CD27 antigen. CD27 (sometimes referred to as T14, S152 or TNFRSF7) is a 50 - 55 kD member of the TNF-R superfamily that is expressed on the surface of cells such as NK cells, B cells and T cells. In some organisms, CD27 is a promoter of T cell differentiation, acts to positively regulate NIK/NF-kappaB signaling and is involved in the negative regulation of apoptotic process. Furthermore, it plays a role in key cellular pathways, in particular, the tumor necrosis factor-mediated signaling pathway, cell surface receptor signaling pathway and extrinsic apoptotic signaling pathway. From a research standpoint, it is of biological interest due to its

association with key macromolecules/ligands like TRAF5, TRAF2 and CD70. CD27 is a moderately popular antibody target, with over 11000 publications in the last decade. CD27 is essential for immunology and costimulatory molecules research, typically serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to APC/iFluor™ 700 (ex/em = 685/710 nm).