

iFluor™ 710 Anti-human CD27 Antibody
LT27Catalog number: 102700K0, 102700K1
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
Isotype	Mouse IgG2a
Immunogen	CD27 (T14, S152, TNFRSF7)
Clone	LT27
Conjugate	iFluor™ 710

Biological Properties

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 710 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 710
Excitation Wavelength	717 nm
Emission Wavelength	739 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 iFluor™ 710 (ex/em = 717/739 nm).