

XFD750 Anti-human CD27 Antibody *LT27*

Catalog Number: 102701B0,

102701B1

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	Lot specific (please consult certificate of analysis for given lot)
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	AF750

Biological Properties

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF750 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined. <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate	AF750
Excitation Wavelength	752 nm
Emission Wavelength	776 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 XFD750 (ex/em = 752/776 nm). XFD750 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 750 (Alexa Fluor® is the trademark of Thermo Fisher).