

TRITC Anti-human CD27 Antibody *O323*Catalog number: 102711H0, 102711H1
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 IgG1, κ
Immunogen	CD27 (T14, S152, TNFRSF7)
Clone	O323
Conjugate	TRITC

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with TRITC under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	TRITC
Excitation Wavelength	544 nm
Emission Wavelength	570 nm

Applications

O323 is an anti-human monoclonal antibody that is specific for the CD27 antigen. CD27 (also known as S152, TNFRSF7 or T14) is a 50 - 55 kD single-pass type I membrane protein that is found on the surface of cells such as T cells, B cells and NK cells. In some organisms, CD27 promotes B cell differentiation, inhibits apoptotic process and positively regulates NIK/NF-kappaB signaling. Additionally, it is a member of vital cellular pathways, for instance, the extrinsic apoptotic signaling pathway, cell surface receptor signaling pathway and tumor necrosis factor-mediated signaling pathway. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands such as CD70. CD27 is a moderately popular antibody target, with over 11000 publications in the last decade. CD27 is vital to 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 TRITC (ex/em = 544/570 nm). It is compatible with the 532 nm laser and 575/25 nm bandpass filter (for example, as in the BD Special Order LSRFortessa™ Cell Analyzer).