

**mFluor™ Violet 450 Anti-human CD27
Antibody *O323***Catalog number: 102710Z0, 102710Z1
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	mFluor™ Violet 450

Biological Properties

Appearance	Light yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Violet 450 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Violet 450
Excitation Wavelength	406 nm
Emission Wavelength	445 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 mFluor™ Violet 450 (ex/em = 406/445 nm). It is compatible with the 405 nm laser and 450/50 nm bandpass filter (for example, as in the Miltenyi Biotec MACSQuant X).