

TRITC Anti-human CD28 Antibody *CD28.2*Catalog number: 102801I0, 102801I1
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
Immunogen	CD28 (Tp44, T44)
Clone	CD28.2
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

CD28.2 is an anti-human monoclonal antibody that forms an immune complex with the CD28 antigen. CD28 (sometimes called T44) is a 44 kD member of the Ig superfamily that is located on the surface of cells such as T cells. CD28 acts in essential cellular pathways, in particular, the cell surface receptor signaling pathway, T cell receptor signaling pathway and apoptotic signaling pathway. Furthermore, in some organisms, it positively regulates isotype switching to IgG isotypes, is an enhancer of interleukin-4 production and upregulates translation. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands such as PI3-kinase, CD80 and CD86. CD28 is a very popular antibody target, with over 30000 publications in the last decade. CD28 is essential for costimulatory molecules and immunology 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 561 nm laser and 572/28 nm

bandpass filter (for example, as in the Agilent Technologies NovoCyte Advanteon).