

PE Anti-human CD5 Antibody *UCHT2*

Catalog number: 100521L0, 100521L1, 100521L2

Unit size: 25 tests, 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	CD5 (Leu1, Ly-1, T1, Tp67)
Clone	UCHT2
Conjugate	PE

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE
Excitation Wavelength	566 nm
Emission Wavelength	574 nm

Applications

UCHT2 is an anti-human monoclonal antibody that is specific for the CD5 antigen. CD5 (sometimes referred to as Tp67) is a 67 kD transmembrane protein that is found on the surface of cells like T cells and B cells. CD5 is involved with essential cellular pathways, for instance, the apoptotic signaling pathway. From a research standpoint, it is of biological interest due to its association with important macromolecules/ligands such as CD72, gp35-37, ZAP-70 and TCR. CD5 is a moderately popular antibody target, with over 11000 publications in the last decade. CD5 has been widely used in immunology and costimulatory molecules research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to PE (ex/em = 566/574 nm). It is compatible with the 561 nm laser and 586/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte).