

PE/Cy7 Anti-human CD5 Antibody *UCHT2*Catalog number: 100521N0, 100521N1, 100521N2
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/Cy7

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/Cy7 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/Cy7
Excitation Wavelength	566 nm
Emission Wavelength	778 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/Cy7 (ex/em = 566/778 nm). It is compatible with the 561 nm laser and 780/60 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).