

PE/Texas Red® Anti-human CD5 Antibody
L17F12Catalog number: 100511S0, 100511S1, 100511S2
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	L17F12
Conjugate	PE/Texas Red®

Biological Properties

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

Spectral Properties

Conjugate	PE/Texas Red®
Excitation Wavelength	567 nm
Emission Wavelength	615 nm

Applications

L17F12 is an anti-human monoclonal antibody that is specific for the CD5 antigen. CD5 (also known as Lyt-1 or Ly-1) is a 67 kD single-pass type I membrane protein that is found on the surface of cells such as B cells and T cells. CD5 is a component of important cellular pathways, namely, the apoptotic signaling pathway. From a research standpoint, it is of biological interest due to its association with essential macromolecules/ligands such as gp35-37, TCR and CD21. CD5 is a moderately popular antibody target, with over 11000 publications in the last decade. CD5 is typically used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of immunology. This antibody was purified through affinity chromatography and conjugated to PE/Texas Red® (ex/em = 567/615 nm). It is compatible with the 561 nm laser and 610/20 nm bandpass filter (for example, as in the BD LSRFortessa™ X-20).