

iFluor™ 568 Anti-human CD5 Antibody
L17F12Catalog number: 100510B0, 100510B1
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 | CD5 (Leu1, Ly-1, T1, Tp67) |
| Clone | L17F12 |
| Conjugate | iFluor™ 568 |

Biological Properties

| | |
|-------------|--|
| Appearance | Purple liquid |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 568 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging |

Spectral Properties

| | |
|-----------------------|-------------|
| Conjugate | iFluor™ 568 |
| Excitation Wavelength | 568 nm |
| Emission Wavelength | 587 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 iFluor™ 568 (ex/em = 568/587 nm). It is compatible with the 561 nm laser and 586/15 nm bandpass filter (for example, as in the Miltenyi Biotec MACSQuant VYB).