

**mFluor™ Violet 540 Anti-human CD5
Antibody *UCHT2***Catalog number: 10052120, 10052121
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 | UCHT2 |
| Conjugate | mFluor™ Violet 540 |

Biological Properties

| | |
|-------------|---|
| Appearance | Yellow liquid |
| Preparation | Antibody purified by affinity chromatography and then conjugated with mFluor™ Violet 540 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging |

Spectral Properties

| | |
|-----------------------|--------------------|
| Conjugate | mFluor™ Violet 540 |
| Excitation Wavelength | 394 nm |
| Emission Wavelength | 537 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 mFluor™ Violet 540 (ex/em = 394/537 nm). It is compatible with the 405 nm laser and 528/46 nm bandpass filter (for example, as in the Luminex Amnis CellStream).