

mFluor™ Red 700 Anti-human CD5 Antibody
HISM2Catalog number: 100500V0, 100500V1
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
Isotype	Mouse IgG1
Immunogen	CD5 (Leu1, Ly-1, T1, Tp67)
Clone	HISM2
Conjugate	mFluor™ Red 700

Biological Properties

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Red 700 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Red 700
Excitation Wavelength	680 nm
Emission Wavelength	695 nm

Applications

HISM2 is an anti-human monoclonal antibody that forms an immune complex with the CD5 antigen. CD5 (sometimes referred to as LEU1, Tp67 or Lyt-1) is a 67 kD single-pass type I membrane protein that is located on the surface of cells like B cells and T cells. CD5 is a component of vital 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 like ZAP-70, TCR and BCR. CD5 is a moderately popular antibody target, with over 11000 publications in the last decade. CD5 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of costimulatory molecules. This antibody was purified through affinity chromatography and conjugated to mFluor™ Red 700 (ex/em = 680/695 nm). It is compatible with the 642 nm laser and 702/85 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).