

**mFluor™ UV375 Anti-human CD195
Antibody *T21/8***Catalog number: 119500X0, 119500X1
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 kappa
Immunogen	CD195 (CCR5)
Clone	T21/8
Conjugate	mFluor™ UV375

Biological Properties

Appearance	Yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ UV375 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ UV375
Excitation Wavelength	351 nm
Emission Wavelength	387 nm

Applications

The T21/8 monoclonal antibody binds to human CD195, a 45 kD glycoprotein frequently located on the surface of T cells and monocytes. CD195 is associated with a variety of biologically interesting macromolecules/ligands, for example, RANTES. CD195 is a relatively rare antibody target, with fewer than 300 publications in the last decade. Even still, CD195 is frequently used in flow cytometry applications as a phenotypic marker

for differentiation of cell types, especially in the study of immunology, neuroinflammation and neuroscience. This antibody was purified through affinity chromatography and conjugated to mFluor™ UV375 (ex/em = 351/387 nm). It is compatible with the 355 nm laser and 379/28 nm bandpass filter (for example, as in the BD FACSCelesta™).