

PerCP Anti-human CD195 Antibody *T21/8*Catalog number: 119501S0, 119501S1, 119501S2
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
Isotype	Mouse IgG1 kappa
Immunogen	CD195 (CCR5)
Clone	T21/8
Conjugate	PerCP

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PerCP under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PerCP
Excitation Wavelength	477 nm
Emission Wavelength	678 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 PerCP (ex/em = 477/678 nm). It is compatible with the 488 nm laser and 695/40 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).