

PE/iFluor™ 750 Anti-human CD193 Antibody
5E8Catalog number: 119301Q0, 119301Q1, 119301Q2
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 IgG2b kappa
Immunogen	CD193 (CCR3, CC CKR3)
Clone	5E8
Conjugate	PE/iFluor™ 750

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 750 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/iFluor™ 750
Excitation Wavelength	566 nm
Emission Wavelength	778 nm

Applications

The 5E8 monoclonal antibody reacts with human CD193, a transmembrane protein typically found on the surface of mononuclear phagocytes and platelets. CD193 is associated with a variety of biologically interesting macromolecules/ligands, for instance, RANTES(CCL5) and MCP-3 (CCL7). CD193 is a relatively rare antibody target, with fewer than 5 publications in the last decade. Even still, CD193 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of immunology. This antibody was purified through affinity chromatography and conjugated to PE/iFluor™ 750 (ex/em = 566/778 nm). It is compatible with the 561 nm laser and

780/60 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte).