

**mFluor™ Violet 540 Anti-human CD193  
Antibody \*5E8\***Catalog number: 11930120, 11930121  
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 IgG2b kappa
Immunogen	CD193 (CCR3, CC CKR3)
Clone	5E8
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**

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 mFluor™ Violet 540 (ex/em = 394/537 nm). It is compatible with the 405 nm laser and 528/45 nm bandpass filter (for example, as in the BD FACSMelody™).