

TRITC Anti-human CD193 Antibody *5E8*Catalog number: 119301H0, 119301H1
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	TRITC

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with TRITC under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	TRITC
Excitation Wavelength	544 nm
Emission Wavelength	570 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 TRITC (ex/em = 544/570 nm). It is compatible with the 532 nm laser and 575/25 nm bandpass filter (for example, as in the BD FACSymphony™ A5).