

**mFluor™ Green 620 Anti-human CD162
Antibody *TC2***Catalog number: 116200U0, 116200U1
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
Immunogen	CD162 (PSGL-1)
Clone	TC2
Conjugate	mFluor™ Green 620

Biological Properties

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Green 620 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Green 620
Excitation Wavelength	525 nm
Emission Wavelength	623 nm

Applications

TC2 is an anti-human monoclonal antibody that forms an immune complex with the CD162 antigen. CD162 (alternatively called PSGL-1) is a transmembrane glycoprotein that is located on the surface of cells such as macrophages, granulocytes and stem cells. CD162 is associated with a variety of biologically interesting macromolecules/ligands, namely, selectins. CD162 is a relatively rare antibody target, with fewer than 400

publications in the last decade. Even still, CD162 is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of cell adhesion and immunology. This antibody was purified through affinity chromatography and conjugated to mFluor™ Green 620 (ex/em = 525/623 nm). It is compatible with the 532 nm laser and 610/20 nm bandpass filter (for example, as in the BD FACSymphony™ A5).