

**mFluor™ Violet 450 Anti-human CD180
Antibody *G28-8***Catalog number: 118000Z0, 118000Z1
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 kappa
Immunogen	CD180 (RP105)
Clone	G28-8
Conjugate	mFluor™ Violet 450

Biological Properties

Appearance	Light yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Violet 450 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Violet 450
Excitation Wavelength	406 nm
Emission Wavelength	445 nm

Applications

The G28-8 monoclonal antibody binds with human CD180, a 105 kD transmembrane protein typically found on the surface of B cells, monocytes and dendritic cells. CD180 is associated with a variety of biologically interesting macromolecules/ligands, for example, MD-1. CD180 is a relatively rare antibody target, with fewer than 400 publications in the last decade. Even still, CD180 is commonly used in flow cytometry

applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology. This antibody was purified through affinity chromatography and conjugated to mFluor™ Violet 450 (ex/em = 406/445 nm). It is compatible with the 405 nm laser and 452/45 nm bandpass filter (for example, as in the Miltenyi Biotec MACSQuant VYB).