

PerCP Anti-human CD180 Antibody *G28-8*Catalog number: 11800150, 11800151, 11800152
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 IgG1 kappa
Immunogen	CD180 (RP105)
Clone	G28-8
Conjugate	PerCP

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PerCP under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PerCP
Excitation Wavelength	477 nm
Emission Wavelength	678 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 PerCP (ex/em = 477/678 nm). It is compatible with the 488 nm laser and 690/50 nm bandpass filter (for example, as in the Beckman Coulter DxFLEx).