

**XFD555 Anti-human CD180 Antibody \*G28-8\***

Catalog Number: 11800150, 11800151

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	Lot specific (please consult certificate of analysis for given lot)
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	AF555

**Biological Properties**

---

Appearance	Red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF555 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined.  <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

**Spectral Properties**

---

Conjugate AF555

Excitation Wavelength 553 nm

Emission Wavelength 568 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 XFD555 (ex/em = 553/568 nm). XFD555 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 555 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 561 nm laser and 577/15 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell Analyzer).