

iFluor™ 350 Anti-human CD160 Antibody
BY55Catalog number: 11600010, 11600011
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 IgM kappa
Immunogen	CD160 (BY55, NK1, NK28)
Clone	BY55
Conjugate	iFluor™ 350

Biological Properties

Appearance	Off-white liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 350 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 350
Excitation Wavelength	345 nm
Emission Wavelength	450 nm

Applications

The BY55 monoclonal antibody recognizes human CD160, a 27 kD transmembrane protein typically located on the surface of natural killer cells and T cells. CD160 is associated with a variety of biologically interesting macromolecules/ligands, for instance, MHC class I. CD160 is a fairly uncommon antibody target, with a little more than 1200 publications in the last decade. Even still, CD160 is commonly used in flow cytometry

applications as a phenotypic marker for differentiation of cell types, especially in the study of costimulatory molecules and immunology. This antibody was purified through affinity chromatography and conjugated to iFluor™ 350 (ex/em = 345/450 nm). It is compatible with the 355 nm laser and 447/60 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell Analyzer).