

**mFluor™ Violet 540 Anti-human CD160
Antibody *BY55***Catalog number: 11600120, 11600121
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	mFluor™ Violet 540

Biological Properties

Appearance	Yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Violet 540 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Violet 540
Excitation Wavelength	394 nm
Emission Wavelength	537 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 mFluor™ Violet 540 (ex/em = 394/537 nm). It is compatible with the 405 nm laser and 525/50 nm bandpass filter (for example, as in the Miltenyi Biotec MACSQuant Analyzer 16).