

**mFluor™ Green 620 Anti-human CD314
Antibody *1D11***Catalog number: 131400U0, 131400U1
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	CD314 (NKG2D)
Clone	1D11
Conjugate	mFluor™ Green 620

Biological Properties

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Green 620 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Green 620
Excitation Wavelength	525 nm
Emission Wavelength	623 nm

Applications

The 1D11 monoclonal antibody binds to human CD314, a transmembrane protein commonly found on the surface of natural killer cells and T cells. CD314 is associated with a variety of biologically interesting macromolecules/ligands, namely, ULBP2, MICB, ULBP1 and MICA. CD314 is a relatively rare antibody target, with fewer than 300 publications in the last decade. Even still, CD314 is essential for costimulatory molecules

research, typically serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to mFluor™ Green 620 (ex/em = 525/623 nm). It is compatible with the 532 nm laser and 609/30 nm bandpass filter (for example, as in the Luminex Guava easyCyte).