

PE/iFluor™ 647 Anti-human CD152 Antibody
BN13Catalog number: 115201Q0, 115201Q1, 115201Q2
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 igg2a, κ
Immunogen	CD152 (CTLA-4)
Clone	BN13
Conjugate	PE/iFluor™ 647

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 647 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/iFluor™ 647
Excitation Wavelength	569 nm
Emission Wavelength	666 nm

Applications

The BN13 monoclonal antibody binds to human CD152, a 33 kD transmembrane protein frequently expressed on the surface of B cells and T cells. CD152 is associated with a variety of biologically interesting macromolecules/ligands, namely, PTP1D and PI3-kinase. CD152 is a relatively rare antibody target, with fewer than 1000 publications in the last decade. Even still, CD152 has a variety of applications in immunology research, often serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to PE/iFluor™ 647 (ex/em = 569/666 nm). It is compatible with the 561 nm laser and 661/20 nm

bandpass filter (for example, as in the Miltenyi Biotec MACSQuant VYB).