

PE/Cy7 Anti-human CD268 Antibody *11C1*Catalog number: 12680100, 12680101, 12680102
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 IgG1 kappa
Immunogen	CD268 (BAFFR, TNFRSF13C)
Clone	11C1
Conjugate	PE/Cy7

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/Cy7 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/Cy7
Excitation Wavelength	566 nm
Emission Wavelength	778 nm

Applications

11C1 is an anti-human monoclonal antibody that forms an immune complex with the CD268 antigen. CD268 (alternatively called TNFRSF13C) is a 19 kD glycoprotein that is found on the surface of cells such as macrophages and granulocytes. CD268 is associated with a variety of biologically interesting macromolecules/ligands, for example, BAFF. CD268 is a relatively rare antibody target, with fewer than publications in the last decade. Even still, CD268 is typically used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of immunology. This antibody was purified through affinity chromatography and conjugated to PE/Cy7 (ex/em = 566/778 nm). It is compatible with the 561 nm laser and 780/60 nm bandpass filter (for example, as in the BD LSRFortessa™ X-20).