

iFluor™ 560 Anti-human CD268 Antibody
11C1Catalog number: 126800A0, 126800A1
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	CD268 (BAFFR, TNFRSF13C)
Clone	11C1
Conjugate	iFluor™ 560

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 560 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 560
Excitation Wavelength	560 nm
Emission Wavelength	571 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 iFluor™ 560 (ex/em =

560/571 nm). It is compatible with the 561 nm laser and 582/15 nm bandpass filter (for example, as in the BD FACSAria™ III).