

TRITC Anti-human CD268 Antibody *11C1*Catalog number: 126801J0, 126801J1
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	TRITC

Biological Properties

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

Spectral Properties

Conjugate	TRITC
Excitation Wavelength	544 nm
Emission Wavelength	570 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 TRITC (ex/em = 544/570 nm). It is compatible with the 561 nm laser and 582/15 nm bandpass filter (for example, as in the BD FACSMelody™).