

TRITC Anti-human CD133 Antibody *293C3*Catalog number: 113301H0, 113301H1
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 IgG2b
Immunogen	CD133 (AC133, Prominin-1, PROM1)
Clone	293C3
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

The 293C3 monoclonal antibody recognizes human CD133, a 120 kD transmembrane protein typically located on the surface of epithelial and fetal liver cells. In many organisms, CD133 is a promoter of nephron tubule epithelial cell differentiation, and is associated with a variety of biologically interesting macromolecules/ligands. CD133 is a moderately popular antibody target, with over 18000 publications in the last decade. CD133 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of neuroscience cell markers, cell biology and 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 586/20 nm bandpass filter (for example, as in the Agilent Technologies

NovoCyte).