

**PE/Texas Red® Anti-human CD133 Antibody
*293C3***Catalog number: 113301R0, 113301R1, 113301R2
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 IgG2b
Immunogen	CD133 (AC133, Prominin-1, PROM1)
Clone	293C3
Conjugate	PE/Texas Red®

Biological Properties

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

Spectral Properties

Conjugate	PE/Texas Red®
Excitation Wavelength	567 nm
Emission Wavelength	615 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

PE/Texas Red® (ex/em = 567/615 nm). It is compatible with the 561 nm laser and 610/20 nm bandpass filter (for example, as in the BD FACS Aria™ Fusion).