

XFD594 Anti-human CD133 Antibody *293C3*

Catalog Number: 11330160, 11330161

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	Lot specific (please consult certificate of analysis for given lot)
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	AF594

Biological Properties

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF594 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined. <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate AF594

Excitation Wavelength 590 nm

Emission Wavelength 618 nm

Applications

The 293C3 monoclonal antibody recognizes human CD133, a 120 kD transmembrane protein typically located on the surface of epithelials 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 XFD594 (ex/em = 590/618 nm). XFD594 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 594 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 592 nm laser and 610/30 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).