

**iFluor™ 647 Anti-human CD334 Antibody
*4FR6D3***Catalog number: 133400F0, 133400F1
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	CD334 (FGFR4)
Clone	4FR6D3
Conjugate	iFluor™ 647

Biological Properties

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 647 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 647
Excitation Wavelength	656 nm
Emission Wavelength	670 nm

Applications

4FR6D3 is an anti-human monoclonal antibody that recognizes the CD334 antigen. CD334 (sometimes called FGFR4) is a 88 kD glycoprotein that is located on the surface of cells such as epithelial cells. CD334 is associated with a variety of biologically interesting macromolecules/ligands, for instance, acidic FGF. CD334 is a relatively rare antibody target, with fewer than 10 publications in the last decade. Even still, CD334 has a

variety of applications in stem cells research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to iFluor™ 647 (ex/em = 656/670 nm). It is compatible with the 638 nm laser and 660/20 nm bandpass filter (for example, as in the Beckman Coulter Navios EX).