

**Purified Anti-human CD334 Antibody
*4FR6D3***

Catalog number: 13340000

Unit size: 100 ug

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	0.5 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1 kappa
Immunogen	CD334 (FGFR4)
Clone	4FR6D3

Biological Properties

Appearance	Liquid
Preparation	Antibody purified by affinity chromatography
Application	Flow Cytometry (FACS), ELISA, HC, Western Blot

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.