

**PacOrange Anti-human CD334 Antibody
*4FR6D3***Catalog number: 133401L0, 133401L1
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 | PacOrange |

Biological Properties

| | |
|-------------|--|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PacOrange under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging |

Spectral Properties

| | |
|-----------------------|-----------|
| Conjugate | PacOrange |
| Excitation Wavelength | 400 nm |
| Emission Wavelength | 551 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 PacOrange (ex/em = 400/551 nm). It is compatible

with the 405 nm laser and 525/50 nm bandpass filter (for example, as in the Miltenyi Biotec MACSQuant VYB).