

**APC/iFluor™ 750 Anti-human CD164
Antibody *67D2***Catalog number: 116401F0, 116401F1, 116401F2
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 IgG1
Immunogen	CD164 (MUC-24, MGC-24)
Clone	67D2
Conjugate	APC/iFluor™ 750

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/iFluor™ 750 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	APC/iFluor™ 750
Excitation Wavelength	754 nm
Emission Wavelength	793 nm

Applications

The 67D2 monoclonal antibody binds with human CD164, a transmembrane glycoprotein typically found on the surface of leukemias, Is and basophils. In certain organisms, CD164 is an inhibitor of cell population proliferation, and is associated with a variety of biologically interesting macromolecules/ligands. CD164 is a relatively rare antibody target, with fewer than 400 publications in the last decade. Even still, CD164 has a variety of applications in immunology 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 APC/iFluor™ 750 (ex/em = 754/793 nm).