

iFluor™ 488 Anti-human CD164 Antibody
67D2Catalog number: 11640050, 11640051
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
Immunogen	CD164 (MUC-24, MGC-24)
Clone	67D2
Conjugate	iFluor™ 488

Biological Properties

Appearance	Orange-red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 488 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 488
Excitation Wavelength	491 nm
Emission Wavelength	516 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 iFluor™ 488 (ex/em = 491/516 nm). It is compatible with the 488 nm laser and 530/40 nm bandpass filter (for example, as in the BD FACSJazz™).