

APC/XFD750 Anti-human CD164 Antibody *67D2*

Catalog Number: 116401D0,
116401D1, 116401D2
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	Lot specific (please consult certificate of analysis for given lot)
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/AF750

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/AF750 under optimal conditions
Application	Flow Cytometry (FACS)
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined. <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate	APC/AF750
-----------	-----------

Excitation Wavelength 651 nm

Emission Wavelength 785 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/XFD750 (ex/em = 756/785 nm). XFD750 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 750 (Alexa Fluor® is the trademark of Thermo Fisher).