

XFD700 Anti-human CD123 Antibody *12H7*

Catalog Number: 11231190, 11231191

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	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, κ
Immunogen	CD123 (IL-3R α)
Clone	12H7
Conjugate	AF700

Biological Properties

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF700 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 μ L/million cells in 100 μ L 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 AF700

Excitation Wavelength 696 nm

Emission Wavelength 719 nm

Applications

The 12H7 monoclonal antibody binds to human CD123, a 70 kD single-pass type I membrane protein often located on the surface of macrophages, hematopoietic progenitors and dendritic cells. CD123 acts in vital cellular pathways, for instance, the cytokine-mediated signaling pathway. From a research standpoint, it is of biological interest due to its association with essential macromolecules/ligands like CD131. CD123 is a fairly uncommon antibody target, with a little more than 3900 publications in the last decade. Even still, CD123 is essential for immunology research, typically serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to XFD700 (ex/em = 696/719 nm). XFD700 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 700 (Alexa Fluor® is the trademark of Thermo Fisher).