

iFluor™ 800 Anti-human CD123 Antibody
12H7Catalog number: 112310N0, 112310N1
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, κ
Immunogen	CD123 (IL-3R α)
Clone	12H7
Conjugate	iFluor™ 800

Biological Properties

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 800 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 800
Excitation Wavelength	801 nm
Emission Wavelength	820 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 iFluor™ 800 (ex/em = 801/820 nm).