

iFluor™ A7 Anti-human CD127 Antibody
A019D5Catalog number: 112700S0, 112700S1
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	CD127 (IL-7R α , IL-7R)
Clone	A019D5
Conjugate	iFluor™ A7

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ A7 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ A7
-----------	------------

Applications

A019D5 is an anti-human monoclonal antibody that targets the CD127 antigen. CD127 (sometimes called IL7R or IL-7R α) is a 60 - 90 kD single-pass type I membrane protein that is located on the surface of cells such as stem cells and T cells. In some organisms, CD127 is an inhibitor of T cell mediated cytotoxicity, acts to positively regulate cell population proliferation and is an enhancer of gene expression. Also, it is involved with important cellular pathways, for instance, the cell surface receptor signaling pathway, interleukin-7-mediated signaling pathway and positive regulation of receptor signaling pathway via STAT. From a research standpoint, it is of biological interest due to its association with essential macromolecules/ligands such as IL-7, PI3-kinase and CD132. CD127 is a fairly uncommon antibody target, with a little more than 7000 publications in the last decade. Even still, CD127 is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of immunology. This antibody was purified through affinity chromatography and conjugated to iFluor™ A7.