

iFluor™ 514 Anti-human CD165 Antibody
SN2Catalog number: 11650060, 11650061
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 kappa
Immunogen	CD165 (AD2, gp37)
Clone	SN2
Conjugate	iFluor™ 514

Biological Properties

Appearance	Orange-red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 514 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 514
Excitation Wavelength	528 nm
Emission Wavelength	555 nm

Applications

The SN2 monoclonal antibody recognizes human CD165, a glycoprotein typically expressed on the surface of monocytes, fibroblasts, T cells and thymocytes. CD165 is associated with a variety of biologically interesting macromolecules/ligands. CD165 is a relatively rare antibody target, with fewer than 10 publications in the last decade. Even still, CD165 is commonly used in flow cytometry applications as a phenotypic marker

for differentiation of cell types, specifically in the study of immunology. This antibody was purified through affinity chromatography and conjugated to iFluor™ 514 (ex/em = 528/555 nm). It is compatible with the 532 nm laser and 575/25 nm bandpass filter (for example, as in the BD FACSymphony™ A5).