

iFluor™ 514 Anti-human CD305 Antibody
NKTA255Catalog number: 13050060, 13050061
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	CD305 (LAIR1)
Clone	NKTA255
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 NKTA255 monoclonal antibody binds to human CD305, a 40 kD single-pass type I membrane protein commonly expressed on the surface of macrophages and natural killer cells. CD305 is associated with a variety of biologically interesting macromolecules/ligands, for example, PTPN11 and PTPN6. CD305 is a relatively rare antibody target, with fewer than 100 publications in the last decade. Even still, CD305 is typically used in

flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of inhibitory molecules and 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 Luminex Guava easyCyte).