

PE/iFluor™ 647 Anti-human CD305 Antibody
NKTA255Catalog number: 130501P0, 130501P1, 130501P2
Unit size: 25 tests, 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	PE/iFluor™ 647

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 647 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/iFluor™ 647
Excitation Wavelength	569 nm
Emission Wavelength	666 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 PE/iFluor™ 647 (ex/em = 569/666 nm). It is

compatible with the 561 nm laser and 661/20 nm bandpass filter (for example, as in the Miltenyi Biotec MACSQuant VYB).