

**PerCP/Cy5.5 Anti-human CD305 Antibody**  
**\*NKTA255\***Catalog number: 130501T0, 130501T1, 130501T2  
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          | PerCP/Cy5.5   |

**Biological Properties**

|             |                                                                                                            |
|-------------|------------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PerCP/Cy5.5 under optimal conditions |
| Application | Flow Cytometry (FACS)                                                                                      |

**Spectral Properties**

|                       |             |
|-----------------------|-------------|
| Conjugate             | PerCP/Cy5.5 |
| Excitation Wavelength | 489 nm      |
| Emission Wavelength   | 679 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 PerCP/Cy5.5 (ex/em = 489/679 nm). It is

compatible with the 488 nm laser and 670/30 nm bandpass filter (for example, as in the BD LSRFortessa™ X-20).