

**XFD700 Anti-human CD305 Antibody \*NKTA255\***

Catalog Number: 13050190, 13050191

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      | Lot specific (please consult certificate of analysis for given lot)         |
| 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          | AF700         |

**Biological Properties**

---

|                       |   |
|-----------------------|---|
| Appearance            | Dark blue liquid  |
| Preparation           | Antibody purified by affinity chromatography and then conjugated with AF700 under optimal conditions  |
| Application           | Flow Cytometry (FACS), Fluorescence Imaging   |
| Recommended Dilutions | For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined.<br><i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i> |

**Spectral Properties**

---

|           |       |
|-----------|-------|
| Conjugate | AF700 |
|-----------|-------|

Excitation Wavelength 696 nm

Emission Wavelength 719 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 XFD700 (ex/em = 696/719 nm). XFD700 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 700 (Alexa Fluor® is the trademark of Thermo Fisher).