

**XFD647 Anti-human CD205 Antibody \*HD30\***

Catalog Number: 12050170, 12050171

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 kappa |
| Immunogen          | CD205 (DEC205)   |
| Clone              | HD30             |
| Conjugate          | AF647            |

**Biological Properties**

---

|                       |   |
|-----------------------|---|
| Appearance            | Dark blue liquid  |
| Preparation           | Antibody purified by affinity chromatography and then conjugated with AF647 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                    AF647

Excitation Wavelength   650 nm

Emission Wavelength    671 nm

## Applications

---

HD30 is an anti-human monoclonal antibody that targets the CD205 antigen. CD205 (alternatively called DEC205) is a 210 kD glycoprotein that is found on the surface of cells such as B cells. CD205 is associated with a variety of biologically interesting macromolecules/ligands, for example, MIR98. CD205 is a fairly uncommon antibody target, with a little more than 1200 publications in the last decade. Even still, CD205 has a variety of applications in immunology research, often serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to XFD647 (ex/em = 650/671 nm). XFD647 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 647 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 642 nm laser and 702/85 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).