

**PE/iFluor™ 700 Anti-human CD231 Antibody**  
**\*B2D\***Catalog number: 123101X0, 123101X1, 123101X2  
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          | CD231 (Tspan-7, TALLA-1, TM4SF2) |
| Clone              | B2D                              |
| Conjugate          | PE/iFluor™ 700                   |

**Biological Properties**

|             |   |
|-------------|---|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 700 under optimal conditions |
| Application | Flow Cytometry (FACS)   |

**Spectral Properties**

|                       |                |
|-----------------------|----------------|
| Conjugate             | PE/iFluor™ 700 |
| Excitation Wavelength | 566 nm         |
| Emission Wavelength   | 708 nm         |

**Applications**

The B2D monoclonal antibody binds with human CD231, a transmembrane protein frequently found on the surface of skeletal muscles, neurons, spleens and leukemias. CD231 is associated with a variety of biologically interesting macromolecules/ligands. CD231 is a relatively rare antibody target, with fewer than 30 publications in the last decade. Even still, CD231 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of immunology. This antibody was purified through affinity chromatography and conjugated to PE/iFluor™ 700 (ex/em = 566/708 nm). It is compatible with the 561 nm laser and 695/40 nm bandpass filter

(for example, as in the Agilent Technologies NovoCyte Advanteon).