

**iFluor™ 750 Anti-human CD231 Antibody**  
**\*B2D\***Catalog number: 123100L0, 123100L1  
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      | 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          | iFluor™ 750                      |

**Biological Properties**

|             |                                                                                                            |
|-------------|------------------------------------------------------------------------------------------------------------|
| Appearance  | Dark blue liquid                                                                                           |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 750 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                |

**Spectral Properties**

|                       |             |
|-----------------------|-------------|
| Conjugate             | iFluor™ 750 |
| Excitation Wavelength | 757 nm      |
| Emission Wavelength   | 779 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 iFluor™ 750 (ex/em = 757/779 nm).