

**APC/XFD750 Anti-human CD231 Antibody**  
**\*B2D, XFD750 Same Structure to Alexa**  
**Fluor™ 750\***

Catalog number: 123101E0, 123101E1, 123101E2  
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          | APC/AF750                        |

### Biological Properties

|             |                                                                                                          |
|-------------|----------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with APC/AF750 under optimal conditions |
| Application | Flow Cytometry (FACS)                                                                                    |

### Spectral Properties

|                       |           |
|-----------------------|-----------|
| Conjugate             | APC/AF750 |
| Excitation Wavelength | 756 nm    |
| Emission Wavelength   | 785 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 APC/XFD750 (ex/em = 756/785 nm). XFD750 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 750 (Alexa Fluor® is the trademark of ThermoFisher).