

iFluor™ 488 Anti-human CD231 Antibody
B2DCatalog number: 12310050, 12310051
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™ 488

Biological Properties

Appearance	Orange-red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 488 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 488
Excitation Wavelength	491 nm
Emission Wavelength	516 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™ 488 (ex/em = 491/516 nm). It is compatible with the 488 nm laser and 530/30 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).