

**iFluor™ 450 Anti-human/monkey CD20
Antibody *2H7***Catalog number: 10202040, 10202041
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, monkey
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse igg2b, κ
Immunogen	CD20 (Bp35, B1)
Clone	2H7
Conjugate	iFluor™ 450

Biological Properties

Appearance	Brown liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 450 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 450
Excitation Wavelength	451 nm
Emission Wavelength	502 nm

Applications

The 2H7 monoclonal antibody binds with human/monkey CD20, a 33 - 37 kD transmembrane glycoprotein commonly expressed on the surface of T cells and B cells. CD20 is a component of essential cellular pathways, in particular, the B cell receptor signaling pathway. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands such as Lyn. CD20 is a moderately popular

antibody target, with over 20000 publications in the last decade. CD20 is vital to immunology research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to iFluor™ 450 (ex/em = 451/502 nm). It is compatible with the 445 nm laser and 510/80 nm bandpass filter (for example, as in the BD FACSAria™ III).