

iFluor™ 488 Anti-human CD20 Antibody
HI20aCatalog number: 10201050, 10201051
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 IgG2a
Immunogen	CD20 (Bp35, B1)
Clone	HI20a
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

HI20a is an anti-human monoclonal antibody that recognizes the CD20 antigen. CD20 (alternatively called MS4A1, Bp35 or Leukocyte surface antigen Leu-16) is a 33 - 37 kD transmembrane protein that is expressed on the surface of cells such as B cells. CD20 plays a role in important cellular pathways, namely, the B cell receptor signaling pathway. From a research standpoint, it is of biological interest due to its association

with key macromolecules/ligands such as Lyn. CD20 is a moderately popular antibody target, with over 20000 publications in the last decade. CD20 has a variety of applications in costimulatory molecules and immunology research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. 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 528/65 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).