

**mFluor™ Violet 450 Anti-human CD22
Antibody *HIB22***Catalog number: 102200Z0, 102200Z1
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	CD22 (BL-CAM, Siglec-2)
Clone	HIB22
Conjugate	mFluor™ Violet 450

Biological Properties

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

Spectral Properties

Conjugate	mFluor™ Violet 450
Excitation Wavelength	406 nm
Emission Wavelength	445 nm

Applications

HIB22 is an anti-human monoclonal antibody that recognizes the CD22 antigen. CD22 (also known as BL-CAM or Lyb8) is a 120 - 130 kD glycoprotein that is located on the surface of cells like B cells and stem cells. CD22 plays a role in essential cellular pathways, namely, the negative regulation of B cell receptor signaling pathway. Moreover, in many organisms, it is a negative regulator of calcium-mediated signaling,

suppresses immunoglobulin secretion and is involved in the negative regulation of B cell receptor signaling pathway. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands like PI3-kinase, PLCg1, p53/56lyn and CD45. CD22 is a fairly uncommon antibody target, with a little more than 4000 publications in the last decade. Even still, CD22 is vital to cell biology 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 mFluor™ Violet 450 (ex/em = 406/445 nm). It is compatible with the 405 nm laser and 450/50 nm bandpass filter (for example, as in the BD FACSARIA™ III).