

PE/iFluor™ 700 Anti-human CD52 Antibody
HI186Catalog number: 105201X0, 105201X1, 105201X2
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 IgG2b
Immunogen	CD52 (CAMPATH-1, HE5)
Clone	HI186
Conjugate	PE/iFluor™ 700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 700 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/iFluor™ 700
Excitation Wavelength	566 nm
Emission Wavelength	708 nm

Applications

HI186 is an anti-human monoclonal antibody that is specific for the CD52 antigen. CD52 (also known as HE5) is a 7 kD transmembrane glycoprotein that is located on the surface of cells such as epithelial cells, NK cells and T cells. In some organisms, CD52 is an enhancer of cytosolic calcium ion concentration, and is associated with a variety of biologically interesting macromolecules/ligands, in particular, . CD52 is a fairly uncommon antibody target, with a little more than 2600 publications in the last decade. Even still, CD52 is vital to costimulatory molecules research, often serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through

affinity chromatography and conjugated to PE/iFluor™ 700 (ex/em = 566/708 nm). It is compatible with the 561 nm laser and 710/50 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).