

iFluor™ 810 Anti-human CD52 Antibody
HI186Catalog number: 10520000, 10520001
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 IgG2b
Immunogen	CD52 (CAMPATH-1, HE5)
Clone	HI186
Conjugate	iFluor™ 810

Biological Properties

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 810 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 810
Excitation Wavelength	811 nm
Emission Wavelength	822 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 iFluor™ 810 (ex/em = 811/822 nm).