

XFD488 Anti-human CD52 Antibody *HI186*

Catalog Number: 10520150, 10520151

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	Lot specific (please consult certificate of analysis for given lot)
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	AF488

Biological Properties

Appearance	Orange liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF488 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined.
	<i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate	AF488
Excitation Wavelength	499 nm
Emission Wavelength	520 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 XFD488 (ex/em = 499/520 nm). XFD488 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 488 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 488 nm laser and 533/30 nm bandpass filter (for example, as in the BD FACSCanto™).