

**mFluor™ Red 780 Anti-human CD205
Antibody *HD30***Catalog number: 120500W0, 120500W1
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 kappa
Immunogen	CD205 (DEC205)
Clone	HD30
Conjugate	mFluor™ Red 780

Biological Properties

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Red 780 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Red 780
Excitation Wavelength	629 nm
Emission Wavelength	767 nm

Applications

HD30 is an anti-human monoclonal antibody that targets the CD205 antigen. CD205 (alternatively called DEC205) is a 210 kD glycoprotein that is found on the surface of cells such as B cells. CD205 is associated with a variety of biologically interesting macromolecules/ligands, for example, MIR98. CD205 is a fairly uncommon antibody target, with a little more than 1200 publications in the last decade. Even still, CD205 has a variety

of applications in immunology 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 mFluor™ Red 780 (ex/em = 629/767 nm). It is compatible with the 633 nm laser and 780/60 nm bandpass filter (for example, as in the BD FACSCelesta™).