

iFluor™ 532 Anti-human CD15 Antibody
HI98Catalog number: 10150070, 10150071
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 IgM
Immunogen	CD15 (3-FAL, X-hapten, Lewis X, SSEA-1)
Clone	HI98
Conjugate	iFluor™ 532

Biological Properties

Appearance	Dark red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 532 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 532
Excitation Wavelength	537 nm
Emission Wavelength	560 nm

Applications

HI98 is an anti-human monoclonal antibody that forms an immune complex with the CD15 antigen. CD15 (also known as X-hapten) is a transmembrane protein that is found on the surface of cells like stem cells, macrophages and granulocytes. CD15 is associated with a variety of biologically interesting macromolecules/ligands, in particular, Selectins. CD15 is a fairly uncommon antibody target, with a little more than 5800

publications in the last decade. Even still, CD15 is vital to neuroscience, cell biology and stem cells research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to iFluor™ 532 (ex/em = 537/560 nm). It is compatible with the 532 nm laser and 575/25 nm bandpass filter (for example, as in the Luminex Guava easyCyte).