

PE/Texas Red® Anti-human CD116 Antibody
4H1Catalog number: 111601R0, 111601R1, 111601R2
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 IgG1 kappa
Immunogen	CD116 (GM-CSFR α chain)
Clone	4H1
Conjugate	PE/Texas Red®

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/Texas Red® under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/Texas Red®
Excitation Wavelength	567 nm
Emission Wavelength	615 nm

Applications

The 4H1 monoclonal antibody binds to human CD116, a 70 - 85 kD glycoprotein commonly located on the surface of endothelial precursors and macrophages. CD116 is associated with a variety of biologically interesting macromolecules/ligands, for instance, CD131 and GM-CSF. CD116 is a relatively rare antibody target, with fewer than 100 publications in the last decade. Even still, CD116 is typically used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of immunology. This antibody was purified through affinity chromatography and conjugated to PE/Texas Red® (ex/em = 567/615 nm). It is compatible with the 561 nm laser and 610/20 nm

bandpass filter (for example, as in the BD FACS Aria™ Fusion).