

PE/iFluor™ 750 Anti-human CD200 Antibody
OX-104Catalog number: 120001R0, 120001R1, 120001R2
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, κ
Immunogen	CD200 (OX-2)
Clone	OX-104
Conjugate	PE/iFluor™ 750

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 750 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/iFluor™ 750
Excitation Wavelength	566 nm
Emission Wavelength	778 nm

Applications

OX-104 is an anti-human monoclonal antibody that forms an immune complex with the CD200 antigen. CD200 (sometimes referred to as OX-2 or OX2) is a single-pass type I membrane protein that is expressed on the surface of cells like stem cells, dendritic cells, B cells and endothelial cells. CD200 has been thought to be involved with vital biological processes such as cell-cell adhesion, especially heterotypic cell-cell adhesion. In addition, in many organisms, it suppresses neuron death, is a negative regulator of macrophage activation and is a negative regulator of interleukin-6 secretion. From a research standpoint, it is of biological interest due to its association with important macromolecules/ligands

such as CD200R1. CD200 is a fairly uncommon antibody target, with a little more than 2000 publications in the last decade. Even still, CD200 is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of neuroscience and immunology. This antibody was purified through affinity chromatography and conjugated to PE/iFluor™ 750 (ex/em = 566/778 nm). It is compatible with the 561 nm laser and 780/60 nm bandpass filter (for example, as in the BD FACSARIA™ Fusion).