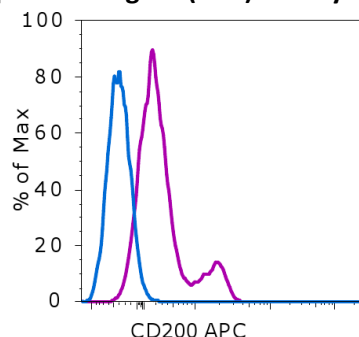


## CD200 APC

**Catalog Number:** 8017-9200

Also known as: OX2

**Analyte Specific Reagent (ASR):** Analytical and performance characteristics are not established.



Fluorescence profiles of normal human peripheral blood lymphocytes unstained (blue histogram) or stained with CD200 APC (purple histogram).

### Product Information



**Contents:** CD200 APC

**Catalog Number:** 8017-9200

**Clone:** OX104

**Concentration:** 5 uL (0.25 ug)/test (a test is defined as the amount that will stain 1 x 10<sup>6</sup> cells in 100 uL)

**Host/Isotype:** Mouse IgG1

**Formulation:** aqueous buffer, 0.09% sodium azide, may contain carrier protein/stabilizer



**Temperature Limitation:** Store at 2-8°C. Do not freeze. Light-sensitive material.



**Batch Code:** Refer to vial



**Analyte Specific Reagent**



**Manufacturer:** eBioscience, Inc., 10255 Science Center Drive, San Diego, CA 92121, USA



**Use By:** Refer to vial

**Caution, contains Azide**

### Description

The monoclonal antibody OX104 recognizes human CD200 also known as OX2. CD200 is a member of the Ig superfamily with 2 Ig domains, a transmembrane and cytoplasmic domain. CD200 is expressed on resting and activated B cells, a subset of resting and activated T cells, keratinocytes, peripheral and central nerve cells, follicular dendritic cells and ovarian cells. The interaction with CD200R results in macrophage activation (IL-6 production), inhibition of mast cell degranulation along with reduced TNF alpha and IL-13 secretion and overall attenuation of the activation status of lymphocytes. A role has also been suggested in maternal tolerance as expression of CD200 is also present on the trophoblast.

Barclay AN, Wright GJ, Brooke G, Brown MH. CD200 and membrane protein interactions in the control of myeloid cells. Trends Immunol. 2002;23(6):285-90.

Wright GJ, Jones M, Puklavec MJ, Brown MH, Barclay AN. The unusual distribution of the neuronal/lymphoid cell surface CD200 (OX2) glycoprotein is conserved in humans. Immunology. 2001;102(2):173-9.

### References

Wright GJ, Cherwinski H, Foster-Cuevas M, Brooke G, Puklavec MJ, Bigler M, Song Y, Jenmalm M, Gorman D, McClanahan T, Liu MR, Brown MH, Sedgwick JD, Phillips JH, Barclay AN. Characterization of the CD200 receptor family in mice and humans and their interactions with CD200. J Immunol. 2003;171(6):3034-46.