

Anti-CD200R antibody

The monoclonal antibody OX108 recognizes human CD200R also known as OX2. CD200R is an inhibitory receptor with a similar structure to its ligand, CD200.

Materials Provided

Catalog Number:	92-1342
Quantity Supplied:	1 vial(s), 100 µl per vial

Description

DiscoverX control ligands have been functionally tested and validated for optimal use with all cell line targets and Assay Ready kits.

Product Information

Molecular Weight:	N/A
Source:	Mouse / IgG1, kappa
Purity:	> 90% by SDS-PAGE
Endotoxin Level:	< 0.001 ng/µg antibody, by LAL assay.
Formulation:	storage buffer in PBS, pH 7.2; contains no preservative
Storage:	4°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

To avoid condensation, equilibrate the vial to ambient temperature before opening.

Stock Concentration:	1 mg/mL
Reconstitution Volume:	N/A
Reconstitution Solvent:	N/A

Handling Instructions:

N/A

For Research Use Only

Ordering: +1.510.979.1415 option 4 or e-mail CustomerServiceDRX@eurofins.com
Technical support: +1.510.979.1415 option 5 or e-mail DRX_SupportUS@eurofinsUS.com
General product information: www.discoverx.com