

Anti-mCD80 Monoclonal Antibody

The monoclonal antibody recognizes mouse CD80, a ~55 kDa cell surface glycoprotein. CD80 (also known as B7-1, B7, and BB1) is expressed on monocytes, macrophages, and dendritic cells in mice.

Materials Provided

Catalog Number:	92-1360
Quantity Supplied:	1 vial(s), 100 ug 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:	Monoclonal antibody
Purity:	Greater than 90%, as determined by SDS-PAGE.
Endotoxin Level:	Less than 0.001 ng/μg antibody, as determined by LAL assay
Formulation:	PBS, pH 7.2
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

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