

Anti-mPD-L1 Monoclonal Antibody

Interaction of the IgSF family ligand, mPD-L1, with its receptor, mPD-1, inhibits T-cell proliferation and cytokine production, and down-regulates the immune response.

Materials Provided

Catalog Number:	92-1353
Quantity Supplied:	1 vial(s), 100 µg 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:	>90% pure as determined by SDS-PAGE analysis
Endotoxin Level:	NA
Formulation:	10 mM NaH ₂ PO ₄ , 150 mM NaCl, 0.09% NaN ₃ , pH7.2
Storage:	4°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	0.5 mg/mL
Reconstitution Volume:	NA
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