

Anti-PD-L1 antibody

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

Materials Provided

Catalog Number:	92-1340
Quantity Supplied:	1 vial(s), 100 mg 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 to human PD-L1; Isotype= IgG1k
Purity:	N/A
Endotoxin Level:	<1 EU per 1 µg
Formulation:	8 mM phosphate, 110 mM NaCl, 2.2 mM KCl, pH 7.4, and 20% glycerol
Storage:	-20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	1.69 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