

Anti-CTLA4 Antibody [14D3]

CTLA4, or CTLA-4 is also known as cytotoxic T-lymphocyte-associated protein 4 or CD152. This monoclonal antibody binds human CTLA4, which functions as a immune checkpoint receptor that downregulates the immune system. CTLA4 is expressed on activated T cells and binds the B7 family members, CD80 (B7-1) and CD86 (B7-2), and the protein restricts the progression of T cells to an activated state by inhibiting IL-2 secretion and cellular proliferation.

Materials Provided

Catalog Number:	92-1319
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:	Mouse Monoclonal to Human CTLA4
Purity:	Functional Grade Purified
Endotoxin Level:	<0.001 ng/µg
Formulation:	N/A
Storage:	-20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	see vial
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