

Anti-PD1 Antibody [NAT105]

Anti-PD1 antibody [NAT105] is a mouse monoclonal antibody that detects PD1 (programmed cell death 1) protein (also known as CD279). The antibody inhibits the inhibitory receptor PD1, potentiates antitumor T-cell response in tumor cells, and prevents PD1 anti-apoptotic activity through inhibition of T-cell anti-tumor activity.

Materials Provided

Catalog Number:	92-1290
Quantity Supplied:	1 vial(s), 0.1 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:	Mouse mAb
Purity:	Immunogen affinity purified
Endotoxin Level:	N/A
Formulation:	PBS, 0.02% azide, 6.97% arginine
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 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