

Anti-TIM3 Antibody [344801]

TIM-3 (T cell immunoglobulin and mucin domain-3) is an immunosuppressive protein that enhances tolerance and inhibits anti-tumor immunity.

Materials Provided

Catalog Number:	92-1313
Quantity Supplied:	1 vial(s), 25 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:	Rat Monoclonal to Human TIM-3.
Purity:	Protein A or G purified from hybridoma culture supernatant
Endotoxin Level:	N/A
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:	0.5 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