

## Anti-VISTA Antibody [730804]

Anti-VISTA Antibody [730804] is a mouse monoclonal antibody that detects human VISTA/B7-H5/PD-1H. VISTA (V-domain Ig suppressor of T cell activation), also known as platelet receptor Gi24, is a member of the Ig superfamily and is a transmembrane molecule expressed in bone, on embryonic stem cells (ESCs), and on tumor cell surfaces. Gi24 interacts with BMP-4, potentiating BMP signaling and the transition from an undifferentiated to a differentiated state; and promotes MT1-MMP expression and activity and serves as a substrate for MT1-MMP.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1311</b>
<b>Quantity Supplied:</b>	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

<b>Molecular Weight:</b>	N/A
<b>Source:</b>	Mouse Monoclonal to Human Vista
<b>Purity:</b>	Protein A or G purified from hybridoma culture supernatant
<b>Endotoxin Level:</b>	N/A
<b>Formulation:</b>	N/A
<b>Storage:</b>	-20°C Please avoid multiple freeze/thaw cycles.

### Reagent Preparation

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

<b>Stock Concentration:</b>	0.5 mg/mL
<b>Reconstitution Volume:</b>	N/A
<b>Reconstitution Solvent:</b>	N/A

***For Research Use Only***

Ordering: +1.510.979.1415 option 4 or e-mail [CustomerServiceDRX@eurofins.com](mailto:CustomerServiceDRX@eurofins.com)  
Technical support: +1.510.979.1415 option 5 or e-mail [DRX\\_SupportUS@eurofinsUS.com](mailto:DRX_SupportUS@eurofinsUS.com)  
General product information: [www.discoverx.com](http://www.discoverx.com)