

Anti-BCMA Monoclonal Antibody

BCMA is a member of the TNF receptor superfamily and is expressed on B-cells. High BCMA expression is associated with cancer in plasma cells.

Materials Provided

Catalog Number:	92-1356
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:	Monoclonal antibody
Purity:	purified by affinity chromatography
Endotoxin Level:	N/A
Formulation:	Phosphate-buffered solution, pH 7.2, containing 0.09% sodium azide.
Storage:	4°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	500 µg/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