

## Anti-ErbB2 Monoclonal Antibody

ErbB2 is a member of the EGFR family that has no ligand binding domain on its own but forms heterodimers with other ligand-binding family members. ErbB2 overexpression/gene amplification has been associated with many types of cancers.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1358</b>
<b>Quantity Supplied:</b>	1 vial(s), 500 uL 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>	Monoclonal antibody
<b>Purity:</b>	Protein G purified
<b>Endotoxin Level:</b>	N/A
<b>Formulation:</b>	PBS, pH 7.4, with 0.2% BSA
<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>	0.2 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)