

Anti-ErbB2 Monoclonal Antibody (trastuzumab) □

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

Catalog Number:	92-1359
Quantity Supplied:	1 vial(s), 200 ug 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:	Protein A affinity purified
Endotoxin Level:	NA
Formulation:	PBS with 0.02% Proclin 300
Storage:	4°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