

## Anti-CD33 Monoclonal Antibody

CD33 is a member of the sialic acid-binding receptor family that is expressed on myeloid cells.  CD33 signaling inhibits macrophage phagocytosis and is associated with various leukemias

### Materials Provided

<b>Catalog Number:</b>	<b>92-1355</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>	Monoclonal antibody
<b>Purity:</b>	purified by affinity chromatography
<b>Endotoxin Level:</b>	NA
<b>Formulation:</b>	Phosphate-buffered solution, pH 7.2, containing 0.09% sodium azide
<b>Storage:</b>	4°C Please avoid multiple freeze/thaw cycles.

### Reagent Preparation

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

<b>Stock Concentration:</b>	500 µg/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)