

## Recombinant Human CD40L

CD40 ligand is a type II transmembrane glycoprotein that belongs to the tumor necrosis factor (TNF) superfamily and is predominantly expressed on T cells, B cells and platelets. Binding of CD40 ligand to CD40 on antigen presenting cells, stimulates B cell proliferation, immunoglobulin class switching, antibody secretion, and T cell-dependent humoral responses.

### Materials Provided

<b>Catalog Number:</b>	<b>92-1348</b>
<b>Quantity Supplied:</b>	1 vial(s), 20 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>	21,600
<b>Source:</b>	Recombinant
<b>Purity:</b>	>95%
<b>Endotoxin Level:</b>	<0.01 EU per 1 µg of the protein
<b>Formulation:</b>	Lyophilized from a 0.2 µm filtered solution in PBS and EDTA with BSA as a carrier protein
<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>	21,600 g/mol
<b>Reconstitution Volume:</b>	0.200 mL
<b>Reconstitution Solvent:</b>	Reconstitution Buffer

***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)