

Recombinant Human CD27 Ligand (TNFSF7)

CD27 Ligand also known as TNFSF7 and CD70 is a 30 kDa type 2 transmembrane glycoprotein that belongs to the tumor necrosis factor (TNF) ligand family. This cytokine is a ligand for TNFRSF27/CD27, expressed as a surface antigen on activated, but not on resting, T and B lymphocytes. It induces proliferation of co-stimulated T cells, enhances the generation of cytotoxic T cells, and contributes to T cell activation.

Materials Provided

Catalog Number:	92-1362
Quantity Supplied:	1 vial(s), 100 µg per vial
Reconstitution Buffer:	1 mL

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:	23,000
Source:	Recombinant
Purity:	>95% (SDS-PAGE)
Endotoxin Level:	<0.10 EU per 1 µg
Formulation:	Lyophilized from a 0.2 µm filtered solution in PBS
Storage:	-20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	200 µg/mL
Reconstitution Volume:	0.500 mL
Reconstitution Solvent:	Reconstitution Buffer

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