

Recombinant Human soluble RANK ligand

Soluble RANK ligand (sRANKL) is a type II protein and member of the tumor necrosis factor (TNF) superfamily. Controls bone regeneration and remodeling, and affects the immune system. Key factor in osteoclast differentiation and activation, dendritic cell survival, and T cell-dependent immune response.

Materials Provided

Catalog Number:	92-1282
Quantity Supplied:	1 vial(s), 10 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:	20 kg/mol
Source:	Recombinant
Purity:	≥ 98%
Endotoxin Level:	< 1EU/µg
Formulation:	N/A
Storage:	-20°C Please avoid multiple freeze/thaw cycles.

Reagent Preparation

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

Stock Concentration:	10 µM
Reconstitution Volume:	0.050 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