

Recombinant Human IL-6

Interleukin 6 (IL-6) is a pro-inflammatory cytokine protein and intricate member of the immune response. Also acts as an anti-inflammatory myokine. Secreted by T cells, macrophages, osteoblasts, and smooth muscle cells for a variety of functions, including stimulating immune response for bacterial infection, burns, trauma, inducing fever, stimulating osteoclast formation, and inducing inflammation. Also inhibits TNF-alpha and IL-1 and activates IL-1ra and IL-10, leading to anti-inflammatory activity. Agonist of the IL-6 receptor, and implicated in multiple inflammatory diseases, including systemic juvenile rheumatoid arthritis and diabetes mellitus susceptibility.

Materials Provided

Catalog Number:	92-1285
Quantity Supplied:	1 vial(s), 20 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:	20900 g/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:	100 µg/mL
Reconstitution Volume:	0.200 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