

Recombinant Human IL-23

Interleukin 23 (IL-23) is a cytokine protein and intricate member of the immune response. Composed of IL-23A and the p40 subunit of IL-12 (IL-12B). Agonist of the receptor formed by IL-12RB1 and IL-23R. Responsible for activation of STAT4 and production of IFNG, acting preferentially on CDR(+) T cells. Activates the JAK-STAT pathway, promoting production of pro-inflammatory cytokines. May be implicated in autoimmune inflammatory diseases and tumorigenesis.

Materials Provided

Catalog Number:	92-1266
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:	53500 g/mol
Source:	Recombinant
Purity:	≥ 95%
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:	0.1 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