

Recombinant Human IL-18

Interleukin 18 (IL-18) is a pro-inflammatory cytokine protein that is an intricate member of the immune response and a member of the IL-1 superfamily. Induces cell-mediated immunity from microbial infection, and works in combination with IL-12 to inhibit IgE and IgG1 production from IL-4 and enhance production of B cells from IgG2a. Negatively regulated by IL-18 binding protein (IL-18BP).

Materials Provided

Catalog Number:	92-1292
Quantity Supplied:	1 vial(s), 25 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:	18 kg/mol
Source:	Recombinant
Purity:	≥ 90% by SDS-PAGE
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:	0.25 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