

Recombinant Human IL-12 (p70)

Interleukin 12 (IL-12) is a pro-inflammatory cytokine protein and intricate member of the immune response that is produced in response to antigenic stimulation. Responds to bacteria, intracellular parasites, and B lymphocytes activity to induce cytokine interferon-gamma and TNF-alpha production from NK and T cells. Mediator of NK and T cell activity, including cytotoxic enhancement.

Materials Provided

Catalog Number:	92-1286
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:	75 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:	100 µg/mL
Reconstitution Volume:	0.100 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