

Recombinant Human IL-24

Interleukin 24 (IL-24) is a pro-apoptotic cytokine protein of the IL10 family that is an intricate member of the immune response. Is recognized as a tumor-suppressing protein due to its control in cell survival and proliferation through activation of the STAT1 and STAT3 transcription factors. Important roles in wound healing, psoriasis, and cancer tissues.

Materials Provided

Catalog Number:	92-1288
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:	18900 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:	0.200 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