

Recombinant Human IL-9

Interleukin 9 (IL-9) is a cytokine protein and intricate member of the immune response. Produced by CD4+ T helper cells, it stimulates cell proliferation, prevents apoptosis, and regulates hematopoietic cells. Agonist of the interleukin-9 receptor (IL9R) and activates the STAT proteins. Implicated in asthma and the pathogenesis of bronchial hyperresponsiveness.

Materials Provided

Catalog Number:	92-1271
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:	14 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:	0.1 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