

Recombinant Human IL-31

Interleukin 31 (IL-31) is a cytokine protein and intricate member of the immune response. Produced by type 2-helper T cells, and is an agonist of STAT3, STAT1, and STAT5 through its activity with the interleukin 31A receptor (IL31RA) and OSMR. May function in skin immunity and inflammation.

Materials Provided

Catalog Number:	92-1261
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:	15800 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