

Recombinant Human IL-17 (IL-17A)

Interleukin 17A (IL-17, IL-17A) is a pro-inflammatory cytokine protein produced by activated T cells. Potent mediator in delayed-type inflammatory reactions that increases chemokine production and recruits monocytes and neutrophils to sites of inflammation. Agonist of the interleukin receptors IL-17RA, IL-17RB, and IL-17RC. High levels are associated with several chronic inflammatory diseases, such as rheumatoid arthritis, psoriasis, and multiple sclerosis.

Materials Provided

Catalog Number:	92-1229
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:	31300 g/mol
Source:	Recombinant
Purity:	≥ 98% by HPLC
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:	0.250 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