

Recombinant Human IL-26

Interleukin 26 (IL-26) is a cytokine protein and intricate member of the immune response that is overexpressed in herpesvirus samimiri-transformed T cells, and thought to contribute to the transformed phenotype of T cells after herpesvirus infection. Roles in mucosal immunity and may be a pro-inflammatory. Inhibited by heparin, and may be implicated in inflammatory bowel disease.

Materials Provided

Catalog Number:	92-1274
Quantity Supplied:	1 vial(s), 25 µg per vial

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:	17700 g/mol
Source:	Recombinant
Purity:	≥ 95%
Endotoxin Level:	< 0.1EU/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:	0.25 mL
Reconstitution Solvent:	NaH ₂ PO ₄ , NaCl and Glycerol

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