

Recombinant Human IFN-g

IFN- γ (interferon-gamma) is produced by CD4 and CD8 T lymphocytes and activated NK cells, and it stimulates a number of lymphoid cell functions, including the anti-microbial and anti-tumor responses of macrophages, NK cells, and neutrophils. IFN- γ receptors are present in most immune cells, which respond to IFN- γ signaling by increasing the surface expression of class I MHC proteins. This promotes the presentation of antigen to T-helper (CD4+) cells. IFN- γ signaling in antigen-presenting cells, and antigen-recognizing B and T lymphocytes, regulates the antigen-specific phases of the immune response.

Materials Provided

Catalog Number:	92-1323
Quantity Supplied:	1 vial(s), 20 μ g 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:	16,800 g/mol
Source:	Recombinant
Purity:	>98%
Endotoxin Level:	N/A
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 μ g/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