

Recombinant Human G-CSF

Granulocyte-colony stimulating factor (G-CSF) is a glycoprotein growth factor, cytokine, and hormone that stimulates bone marrow to produce granulocytes and stem cells and release them into the bloodstream. Also functions in the differentiation, proliferation, and survival of neutrophils via the JAK/STAT, Ras/MAPK, and PI3K/Akt signal transduction pathways.

Materials Provided

Catalog Number:	92-1143
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:	18700 g/mol
Source:	Recombinant
Purity:	≥ 98% by HPLC
Endotoxin Level:	< 0.1 ng/µg (<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:	0.100 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